

HIGHSCORE AND HIGHSCORE PLUS QUICK START GUIDE





HIGHSCORE AND HIGHSCORE PLUS QUICK START GUIDE

EDITION NOTICE: 4022 339 14083, April 2020

This is the original publication of Edition 3 of this document, to be used with the HighScore and HighScore Plus version 4.9 or higher.



DISCLAIMER

Although diligent care has been used to ensure that the information in this material is accurate, nothing herein can be construed to imply any representation or warranty as to the accuracy, correctness or completeness of this information and we shall not be liable for errors contained herein or for damages in connection with the use of this material. Malvern Panalytical reserves the right to change the content in this material at any time without notice.

COPYRIGHT NOTICE

© 2020 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.

Ma	alvern Panalytical Ltd.	Malvern Panalytical B.V.
Gr	ovewood Road, Malvern,	Lelyweg 1, 7602 EA Almelo,
W	orcestershire, WR14 1XZ,	The Netherlands
Ur	nited Kingdom	
Te	l. +44 1684 892456	Tel. +31 546 534 444
Fa	x. +44 1684 892789	Fax. +31 546 534 598
ofo G) maly arrange tigal com	

info@malvernpanalytical.com www.malvernpanalytical.com



TABLE OF CONTENTS

Chapter 1. Introduction	. 5
1.1 Introduction	. 5
1.2 HighScore Plus	. 5
Chapter 2. Get started	6
2.1 Introduction	. 6
2.2 Start the software	. 6
2.3 Set the desktop to default settings	. 7
2.4 Customize the desktop layout	. 9
2.5 Get PANalytical Example Database prepared	11
Chapter 3. Load and show data	13
3.1 Introduction	13
3.2 Load a scan	13
3.3 Show a scan	14
3.3.1 About display panes	14
3.3.2 Use display panes.	16
3.4 Retrieve a reference pattern.	18
3.5 Show a reference pattern	20
Chapter 4. Use pattern treatments.	23
4.1 Introduction	23
4.2 Find background	23
4.3 Search peaks	25
Chapter 5. Do an automatic profile fitting	28
5.1 Introduction	28
5.2 Prepare for an automatic profile fitting	28
5.3 Start an automatic profile fitting	29
Chapter 6. Do search - match - identify	35
6.1 Introduction	35

6.2 Search and match	35 36
Chapter 7. Change scores.	40
7.1 Introduction.	40
7.2 Change scores.	40
Chapter 8. Use a user batch	42
8.1 Introduction	42
8.2 Use a user batch	42
Chapter 9. Phase identification strategy and troubleshooting	44
9.1 Introduction	44
9.2 Pattern treatment sequence	45
9.2.1 Find background	45
9.2.2 Search peaks with a high significance	46
923 Convert intensity	46
	-0
9.2.4 Strip Ka ₂	46
9.3 Identify.	46 46
 9.2.4 Strip Kα₂ 9.3 Identify 9.4 Troubleshooting 	46 46 47
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus).	46 46 47 49
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction.	46 46 47 49
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction. 10.2 load a measurement. 	 46 46 47 49 49 49
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction. 10.2 load a measurement. 10.3 Search peaks. 	 46 46 47 49 49 50
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction. 10.2 load a measurement. 10.3 Search peaks. 10.4 Search and refine a unit cell. 	 46 46 47 49 49 50 51
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction. 10.2 load a measurement. 10.3 Search peaks. 10.4 Search and refine a unit cell. 10.5 Save results. 	46 46 47 49 49 50 51 53
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction. 10.2 load a measurement. 10.3 Search peaks. 10.4 Search and refine a unit cell. 10.5 Save results. Chapter 11. Do an automatic Rietveld refinement (HighScore Plus).	46 46 47 49 49 49 50 51 53 54
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction. 10.2 load a measurement. 10.3 Search peaks. 10.4 Search and refine a unit cell. 10.5 Save results. Chapter 11. Do an automatic Rietveld refinement (HighScore Plus). 11.1 Introduction.	46 46 47 49 49 50 51 53 54
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction. 10.2 load a measurement. 10.3 Search peaks. 10.4 Search and refine a unit cell. 10.5 Save results. Chapter 11. Do an automatic Rietveld refinement (HighScore Plus). 11.1 Introduction. 11.2 Load data.	46 46 47 49 49 50 51 53 54 54 54
 9.2.4 Strip Kα₂. 9.3 Identify. 9.4 Troubleshooting. Chapter 10. Search and refine a unit cell (HighScore Plus). 10.1 Introduction. 10.2 load a measurement. 10.3 Search peaks. 10.4 Search and refine a unit cell. 10.5 Save results. Chapter 11. Do an automatic Rietveld refinement (HighScore Plus). 11.1 Introduction. 11.2 Load data. 11.3 Do an automatic refinement.	46 46 47 49 49 49 50 51 53 51 53 54 54 54



CHAPTER 1 INTRODUCTION

1.1 Introduction

This Quick Start Guide helps you to get started with HighScore and HighScore Plus quickly.

The examples in this Quick Start Guide show you how to use the software to do simple tasks.

You will use the measurement "Mixture3.xrdml" for many examples in the Quick Start Guide. The sample used for this measurement is an artificial mixture of 3 minerals. The PANalytical Example Database, supplied with HighScore and HighScore Plus, is applicable to this measurement.

For all the tasks in the examples, there can be other possible procedures to complete them, but then you must experiment with HighScore or HighScore Plus and learn the other possibilities yourself.

There can be differences between the example screens in this Quick Start Guide and what you see on your screen. If that occurs, follow the instructions on your screen.

1.2 HighScore Plus

The software packages HighScore and HighScore Plus are almost the same, only that HighScore includes more functionality, such as Rietveld, Pawley, LeBail refinements and crystallographic analysis.



CHAPTER 2 GET STARTED

2.1 Introduction

In this chapter, you will do these tasks:

- Start the software.
- Set the desktop to default settings.
- Customize the desktop layout.
- Select PANalytical Example Database, which is supplied with the software.

2.2 Start the software

To start HighScore, use one of these procedures:

- On the desktop, double-click the icon 🛗
- Click the Windows Start button and then on the list of apps or programs, go to PANalytical HighScore > HighScore.

To start HighScore Plus, use one of these procedures:

- On the desktop, double-click the icon 🎬.
- Click the Windows Start button and then on the list of apps or programs, go to PANalytical HighScore Plus > HighScore Plus.



Figure 2.1 Initial window of HighScore Plus

2.3 Set the desktop to default settings

Before you start with the examples in this Quick Start Guide, we recommend that you set the desktop back to its default settings.

NOTE: After you completed the examples, you can change the settings to your preferences.

- 1. On the menu bar, go to **View > Toolbars > Customize** to open the **Customize** window.
- 2. Go to the **Options** tab.
- 3. Clear the Show full menus after a short delay check box.
- 4. Clear the Menus show recently used commands first check box.

Customize	X
Toolbars Commands Options	
Personalized Menus and Toolbars	
Menus show recently used commands first	
Show full menus after a short delay	
Reset my usage data	
Cutter	
Show Tool <u>T</u> ips on toolbars	
Show shortcut keys in ToolTips	
Menu animations: (None)	

Figure 2.2 Customize window

- 5. Click Close.
- 6. On the menu bar, go to **Customize > Program Settings** to open the **Program Settings** window.
- 7. Click Reset All to Default.

ulation Graphics General Reh	erence Patterns	Automatic Processing	Fitting/Rietveld	d Skinning
5ave	Display -			
🗸 Auto save	Show <u>m</u> ulti line tabs		📝 <u>G</u> ene	rate .BAK files
Auto <u>s</u> ave time [minutes]	📝 Show sta	itus bar <u>h</u> ints	Put in	Recent Files
5 💲	<u>T</u> oolbar drav	v style:	🔘 Las	t opened files
Number of Undo/Redo steps:	Use Skins	-		the second files
10 🗘			O Las	st saved <u>fi</u> les
Brems Cryst HKI D-spacing [Å]			Number	of recent files:
3,136 🌲	Edit Lis	t Item Digits		10 🌲
			Auto bro	wse display time [ms]:
				1000 🖵
Use the same recent folder for	opening and for	inserting <u>d</u> ata		
Folder for t <u>e</u> mplates:				
C:\Users\luyao.zhang\AppData\R;	oaming\Microsoft	:\Windows\Templates\	•	Browse
Full User name:				
			B	eset to Default

Figure 2.3 Reset All to Default

8. When the **Confirm** windows opens, click **Yes**.



Figure 2.4 Confirm to overwrite all personal settings

- 9. In the **Program Settings** window, click **OK**.
- 10. On the menu bar, go to **Customize > Defaults** to open the **Default** window.
- 11. Click Restore Factory Defaults.

refaults				
Instrument Settings Sample	/Container Settings	Global Sett	ings Phase Settings External Stand	dards
📃 Incident beam monoch	romator		<u>G</u> oniometer radius [mm]:	240.00 💲
<u>A</u> node material:	Copper (Cu)	•	Distance focus-div. slit [mm]:	91.00 💲
K-a <u>1 [</u> Å]:	1.540598	\$	Divergence <u>s</u> lit type:	Fixed •
K-a <u>2</u> [Å]:	1.544426	\$	Eixed divergence slit size [°]:	1.00 🗘
<u>K</u> -a [Å]:	1.541874	\$	ADS irradiated length [mm]:	10.00 ‡
к- <u>в</u> [Å]:	1.392250	\$	<u>T</u> ime per step [s]:	1.00 ‡
K-a2 / K-a1 rati <u>o</u> :	0.500000	\$	📝 Incident beam <u>s</u> oller slit present	
K-β filter <u>t</u> hickness [mm]:	0.020	\$	📝 Incident beam mask present	
K-β filter material:	Ni		Inc. beam Soller slit opening [rad]:	0.040 💲
S <u>c</u> an axis:	Gonio	•	Incident beam mask width [mm]:	15.00 💲
Rec. slit si <u>z</u> e [mm]:	0.1000	¢	Incident beam mask position [mm]:	109.00 💲
<u>R</u> estore Factory Defa	ults			OK Cancel

Figure 2.5 Restore Factory Defaults

- 12. Click **OK**.
- 13. On the menu bar, go to **File > New** to open a new empty document.
- 14. On the menu bar, go to **View > Reset all Toolbars**.
- 15. When the Confirm windows opens, click Yes.



Figure 2.6 Confirm to reset all menu bars and tool bars

- 16. On the menu bar, go to View > Panes Default Setting.
- 17. On the menu bar, go to View and make sure that Lock Pane Positions is selected.

2.4 Customize the desktop layout

- 1. To show a toolbar or a pane on the desktop, on the menu bar, go to **View > Toolbars** and select the pane.
- 2. To hide a toolbar or a pane on the desktop, on the menu bar, go to **View > Toolbars** and clear the selection of the pane.
- 3. To adjust the relative size of a pane, drag the horizontal and vertical splitter bars of the pane.
- 4. To save changes to the desktop layout, do as follows:
 - a. On the menu bar, go to **View > Desktop**.
 - b. If you save the desktop layout under the current name, click **Save Desktop**.
 - c. If you save the desktop layout under a new name, enter the name in the **Desktop Name** field and click **Save Desktop**.

- 5. To put different panes between 2 monitors, do as follows:
 - a. On the menu bar go to **View**.
 - b. Make sure that **Lock Pane Positions** is not selected.
 - c. Put panes between 2 monitors.
 - d. Save it as a desktop layout.
- 6. To set the desktop layout to a pre-programmed desktop layout, for example **Phase-ID**, do as follows:
 - On the **Desktop** toolbar, select **Phase-ID** in the **Desktop Name** field.



Figure 2.7 Select the desktop layout

• Alternatively, on the menu bar, go to View > Desktop and select Phase-ID.



Figure 2.8 Desktop layout "Phase-ID"

You can switch among different desktop layouts and examine the differences in the desktop layouts.

NOTE: You can also set the desktop layout to **<None>** to use no pre-programmed desktop layout. Then the software will save your settings automatically when you close the last document.

For the examples in this chapter, we recommend that you set the desktop layout to **Phase-ID**, because a desktop layout with the **Main Graphics** pane at this size is the most usual layout to start with.

2.5 Get PANalytical Example Database prepared

A very small database PANalytical Example Database is supplied with the software. You will use it for the examples in the Quick Start Guide.

NOTE: Do not use PANalytical Example Database to do phase analysis on unknown samples.

Do not use PANalytical Example Database to do a test of the functionality or examine the phase identification capabilities of the software. Use a large reference database with at least 100,000 patterns instead, for example an ICDD product or the free COD database.

Do not add your own user reference patterns to PANalytical Example Database. Refer to the HighScore Help file for the information about how to make a new, empty reference database.

To get PANalytical Example Database prepared for patterns retrieval, do as follows:

- 1. On the menu bar, go to **Customize > Manage Databases** to open the **Manage Databases** window.
- 2. Examine if **PANalytical Example Database** is in the list.



Figure 2.9 PANalytical Examine Database in the list

- 3. If **PANalytical Example Database** is not in the list, add it to the list:
 - a. Right-click in the Manage Databases window.
 - b. From the pop-up menu, select Add HighScore Database. The Open window opens.
 - c. Open the folder C:\Documents and Settings\user.name\My Documents\PANalytical \HighScore Plus\ExampleDb.
 - d. Select the file "Codes.pdb".

- e. Click Open. PANalytical Manage Database is added to the list.
- 4. In the list, make sure that the **Use** check box is selected for **PANalytical Example Database**.
- 5. In the list, click **Properties** for **PANalytical Example Database**. If the properties are shown, the database can be used.

NOTE: It is not necessary to convert PANalytical Example Database before you use it.

In the **Manage Database** window, the pie chart shows the total number of patterns in the database.

6. Click × or go to **File > Close** on the menu bar to close the document.



CHAPTER 3 LOAD AND SHOW DATA

3.1 Introduction

The most frequent tasks in HighScore and HighScore Plus are to load and show data. In this chapter, you will do some simple examples.

There are more than 1 possible procedures to show data. Different views can also be used together.

3.2 Load a scan

- 1. If the software is not started, start it. Refer to Section 2.2.
- 2. On the menu bar, go to **File > Open**. The Open **window** opens.
- 3. Open the folder C:\Documents and Settings\user.name\My Documents\PANalytical \X'Pert HighScore Plus\Tutorial
- 4. In the **Open** window, select **All files (*.*)** from the drop-down list to show all files in this folder.

Organize 👻 New fold	er	•••• •	0
Favorites Favorites Desktop Downloads ConeDrive Recent Places Libraries Documents Music Pictures Videos	CEO2-NBS.RD CeO2-RietveldFit Ceramic.rd Clinker Clinker Cluster Corundum#1976b Coto.RD Cryst10.RD Cryst50.RD Cryst50+.RD Cryst30+.RD Crystalline + Amorph CSAND.PD3		
 File <u>n</u>	ame: Mixture3 🔹 🖌	ïles (*.*)	•

Figure 3.1 Tutorial folder with all files shown

- 5. Select the file "Mixture3.xrdml".
- 6. Click **Open** to open the file.

A copy of the selected file is loaded into a new document. Because this is the first scan of the document, it automatically becomes the anchor scan.



Figure 3.2 Mixture3.xrdml

3.3 Show a scan

3.3.1 About display panes

There are different panes in the software. Take the desktop **Phase-ID** as an example.

At the left side of the window, the **Main Graphics** pane shows the full anchor scan in **Analyze View**.



Figure 3.3 Main Graphics pane in Analyze view

Below the **Main Graphics** pane is the **Additional Graphics** pane. It shows where the zooming is done on the scan in the **Main Graphics** pane. If there is no zooming, it shows the full range of the scan in opposite colors.



Figure 3.4 Full range in opposite colors

On the right side of the **Main Graphics** pane and **Additional Graphics** pane, there are more panes, for example the **Peak List** pane, the **Refinement Control** pane and the **Scan List** pane. The first time the software is used, the **Pattern List** pane is shown on top. Then the pane that is used the latest is shown on top.

Q Acces	nted Re	f Dattern I	Mone	a Scan Data	- A	Fattern List	х
No.	Visi	Ref. Code	None	Compound t	Va	Chemical Form	Scor
		- 110					
	idatasi	. 110		a			
Cand	idates:	- Mi	(A) c		-	Chaptical Economia	

Figure 3.5 Pattern List pane on top of other panes

3.3.2 Use display panes

- 1. In the **Main Graphics** pane, drag a square on a peak to zoom in. Try to show only part of a peak and examine the changes of the graph in the **Additional Graphics** pane at the same time.
- 2. Double-click in the **Main Graphics** pane to zoom out.
- 3. Zoom in along the y-axis:
 - a. On the **Tool Palette**, click 📠 to switch on the **Zoom Intensity** function.



Figure 3.6 Zoom Intensity button

- b. In the **Main Graphics** pane, drag a square along the y-axis. The graph is zoomed in along y-axis. Try to show only part of the peak and examine the changes of the graph in the **Additional Graphics** pane at the same time.
- 4. Set the scale of the y-axis:
 - a. On the **Tool Palette**, click the small arrow next to \square .
 - b. Select one of these scales:
 - Linear Y-Axis.
 - · Square Root Y-Axis
 - Logarithmic Y-Axis



Figure 3.7 Set y-axis scale

3.4 Retrieve a reference pattern

- 1. On the menu bar, go to **Reference Patterns > Retrieve Pattern by > Restrictions**. The **Restrictions** window opens.
- 2. Go to the **Strings** tab.
- 3. In the Mineral Name field, enter "calcite".
- 4. Click Load.

[onded]			
Subfiles Chemistry	2uality Crystallography 🕨 Strings Mineral/Zeolite Class		
		Exact Match	Load
<u>⊂</u> ompound Name:		··· 🗙 🔳	Save as S <u>u</u> bs
<u>M</u> ineral Name:	calcite	··· 🗙 🔳	
Eormula:		··· 🗙 🔳	
C <u>o</u> lor:		··· 🗙 🔳	
<u>A</u> uthor:		··· 🗙 🔳	
Journal:		··· 🗙 🔳	
	Clea	ł	Close

Figure 3.8 Retrieve a reference pattern by restrictions

- 5. Click **Close** to close the **Restrictions** window.
- 6. Examine these changes in the display panes:
 - The Main Graphics pane shows the reference pattern lines in Analyze View.
 - The Additional Graphics pane shows the reference pattern lines in most views.
 - The **Pattern List** pane shows a summary of the retrieved reference pattern.



Figure 3.9 Reference pattern shown in display panes

 In the Main Graphics pane, go to the Pattern View tab. The reference pattern lines are shown without the scan data. No peaks are shown in Peak List, because the Peak List pane is empty.



Figure 3.10 Reference pattern in Pattern View

3.5 Show a reference pattern

- 1. Go to the **Pattern List** pane.
- 2. Show the details of the reference pattern:
 - Right-click in the Pattern List pane and select Show Pattern from the pop-up menu.



Figure 3.11 Accepted reference patterns pop-up menu

• Alternatively, in the **Pattern List** pane, double-click in the **Accepted Ref. Pattern** grid. The **Reference Pattern** window opens and shows the details of reference pattern.

Reference Pattern: 00-005-0586		x
Name and formula		
Reference code:	00-005-0586	
Mineral name: Compound name: PDF index name:	Calcite, syn Calcium Carbonate Calcium Carbonate	≡
Empirical formula: Chemical formula:	CCaO ₃ CaCO ₃	
<u>Crystallographic parameter</u>	<u>-s</u>	
Crystal system: Space group: Space group number:	Rhombohedral R-3c 167	
a (Å): b (Å): c (Å): Alpha (°): Beta (°): Gamma (°):	4.9890 4.9890 17.0620 90.0000 90.0000 120.0000	
Calculated density (g/cm ³): Measured density (g/cm ³): Volume of cell (10 ⁶ pm ³): Z:	2.71 2.71 367.78 6.00	
RIR:	2.00	
Subfiles and quality		
Subfiles:	Cement and Hydration Product, Common Phase, Educations pattern, Forensic, Inorganic, Mineral, NBS pattern, Pharmaceutical, Superconducting Material	le •
< > Save As ⊆opy Print	Graphics Print All Intensity Scale 🔻 Angle Scale	• •

Figure 3.12 Reference pattern details

- 3. Close the Reference Pattern window.
- 4. Open the **Analyze View** again:
 - In the Main Graphics pane, go to the Analyze View tab.
 - Alternatively, on the menu bar, go to View > Main Graphics > Analyze View.
- 5. On the **Display Mode** toolbar, click k to switch on the **Show Reference Patterns** function. Then you can adjust the display of reference pattern lines in the **Main Graphics** pane.

<u>F</u> ile	<u>E</u> dit	⊻iew	Treat <u>m</u>	lent	Reference <u>P</u> atterns	<u>A</u> nalysis	<u>R</u> eports	<u>T</u> ools
Pos	. [°28]	: 26.643			d-spacing [Å]: 3.3	431	Counts	:
	\mathcal{M}	$\mathrm{dir}\star$	- 4	٩٨	1 da • 🜆 😂 🖊	赵杰	🗛 👻 📊	
	ا 😭	4ixture3		v Show	/ Reference Patterns]		

Figure 3.13 Show Reference Patterns button

- **NOTE:** The high intensity areas of the scan marked in gray are those 'features' of the scan explained by the loaded reference pattern. The first 20 reference patterns are automatically matched and scored against the anchor pattern.
- 6. Save the document:
 - a. On the menu bar, go to **File > Save Document**.
 - b. Use these settings:

File name	Mixture3
Save as type	HighScore Plus (*.HPF)

c. Click Save.

7. Click × or go to **File > Close** on the menu bar to close the document.



CHAPTER 4 USE PATTERN TREATMENTS

4.1 Introduction

Pattern treatment is used to prepare data for phase analysis and crystallographic analysis and sometimes for a structure refinement. With pattern treatment, you can make changes to the data and get additional derived data.

The 2 most important pattern treatments are to find background and peaks.

For phase analysis with the measured net profile data, it is very important to find background correctly. If you include peak data in the search-match-identification process or if you do profile fitting or indexing, peak search is necessary.

4.2 Find background

- 1. If the software is not started, start it. Refer to Section 2.2.
- 2. Open the document "Mixture3.xrdml". Refer to Section 3.2.

NOTE: Be careful that you do not open the document "Mixture3.hpf".

3. On the menu bar, go to **Treatment > Determine Background** to find the background. The **Determine Background** window opens. The background is immediately found and shows as a bright green line in the **Main Graphics** pane.



Figure 4.1 Anchor scan with the background

NOTE: Usually the software starts to process data only after you click a button, but after you go to **Treatment > Determine Background**, the software finds the background immediately.

4. Read the title bar of the **Determine Background** window. The title bar shows the parameter set that you use.



Figure 4.2 Use the Identify1 parameter set

- 5. Make sure that you use the **Identify1** parameter set.
- 6. If not, change the parameter set to **Identify1**:
 - a. Click More to expand the window.
 - b. In the Select Parameter Set field, select Identify1 from the drop-down list.
- 7. Change the bending factor and examine how the background changes:
 - a. On the **Automatic** tab, write down the value in the **Bending factor** field. You will set **Bending factor** back to this value later.
 - b. Move the slider below the **Bending factor** field to change its value and examine how the background changes in the **Main Graphics** pane at the same time.
 - c. If necessary, zoom in to examine small changes.
- 8. Change the granularity and examine how the background changes:
 - a. On the **Automatic** tab, write down the value in the **Granularity** field. You will set **Bending factor** back to this value later.
 - **NOTE:** This parameter changes the number of intervals which are used to find the background. The default value "20" can be used for most scans.
 - b. Move the slider below the **Granularity** field to change its value and examine how the background changes in the **Main Graphics** pane at the same time.
 - c. If necessary, zoom in to examine small changes.
- 9. Read the title bar of the **Determine Background** window again. At this time, the title is changed to **Determine Background [Untitled]**. This shows the parameter set is changed and not saved under a specific name.
- 10. Set the **Bending factor** and **Granularity** fields back to the initial values that you wrote down in steps 7 and 8.
- 11. Click **Accept** to accept the background. The accepted background shows in the **Main Graphics** pane.
 - **NOTE:** The color of the background is set on the **Graphic Colors** tab in the **Document Settings** window. Right-click in the **Main Graphics** pane and select **Document Settings** from the pop-up menu to open the **Document Settings** window.
- 12. Examine the anchor scan data:
 - a. Go to the Anchor Scan Data pane.

b. Examine the data in the **Iback (cts)** column. This is the background data that you just made.

Lis	ts Par	ne						Þ
	Qu	uantification	Rel	finement Co	introl	Object	: Inspector	
	Pattern List		ern List Scan List Peak List			Anchor Scan Data		
	No.	Pos. [*2Th.]	lobs [cts]	Icalc [cts]	Iback [cts]	CT [s]	ESD I) sp 🔺
Þ	1	19.9414	906.9789		906.9789	19.6850	4	4.44
	2	19.9584	901.5664		904.0214	19.6850	4	4.44!
	3	19.9754	941.7132		901.0640	19.6850	4	4.44
	4	19.9924	893.4242		898.1065	19.6850	4	4.43
	5	20.0094	906.2788		895.1490	19.6850	4	4.43:
	6	20.0264	910.6924		892.1916	19.6850	4	1.43
	7	20.0434	905.7835		889.2341	19.6850	4	1.421
	8	20.0604	901.7400		886.2767	19.6850	4	4.42:
	9	20.0774	890.2701		883.3192	19.6850	4	4.41:
	10	20.0944	898.5697		880.3618	19.6850	4	.41!
	11	20.1114	879.0623		877.4043	19.6850	4	1.41
	12	20.1284	865.1569		874.4469	19.6850	4	4.40
	13	20.1454	870.3448		871.4894	19.6850	4	4.40
	14	20.1624	921.6106		868.5320	19.6850	4	4.40
	15	20.1794	900.3833		865.5745	19.6850	4	1.391
	16	20.1964	893.8984		862.6171	19.6850	4	1.39:
	17	20.2134	863.9377		859.6596	19.6850	4	1.38
	18	20.2304	876.5692		856.7022	19.6850	4	1.38!
	10	00.0474	000 4070		000 7445	40.0050		00

Figure 4.3 Anchor Scan Data tab

4.3 Search peaks

- 1. On the menu bar, go to **Treatment > Search Peaks** to open the **Search Peaks** window.
- 2. Read the title bar of the **Search Peaks** window. The title bar shows the parameter set that you use.

Search Peaks	[Identify]		×
Mi <u>n</u> imum signific	ance:	2.00	Search Peaks
Minim <u>u</u> m tip wid	th [°2Th.]:	0.01	Accept
Maximum tip <u>w</u> io	th [°2Th.]:	1.00	
Pea <u>k</u> base width	n [°2Th.]:	2.00	
Met <u>h</u> od:	Minimum 2nd de	erivative 🔽	Close
Tr <u>i</u> al:		~	<u>M</u> ore >>

Figure 4.4 Use the Identify parameter set

- 3. Make sure that you use the **Identify** parameter set.
- 4. If not, change the parameter set to **Identify**:
 - a. Click **More** to expand the window.
 - b. In the **Select Parameter Set** field, select **Identify** from the drop-down list.
- 5. Click **Search Peaks**. The peaks are found and are shown in the **Main Graphics** pane:



Figure 4.5 Anchor scan with peaks and background data

- Solid lines: Kα₁and Kα_{mixed} peaks.
- Dashed lines: Kα₂ peaks.
- V marks above: peaks that are not explained by a reference pattern
- Blue line: the calculated profile

NOTE: The default color of the calculated profile is blue. The color can be changed on the Graphic Colors tab in the Document Settings window. Right-click in the Main Graphics pane and select Document Settings from the pop-up menu to open the Document Settings window.

This is only a preview of the result. If you close the window, no peaks are written into the diffraction document. You can do a check of the quality of the peak search by the peak lines and the profile before you accept any result. You can zoom in for more details.

- 6. In the Search Peaks window, click Accept to add the peaks that are found to the document.
- 7. Change the display of peaks:
 - On the **Display Mode** toolbar, click the small arrow next to <u>u</u> to select the display of peaks.



Figure 4.6 Display of peaks

- Alternatively, click 🔤 to change the display of peaks. Each time you click, the display of peaks is changed.
- 8. Go to the **Peak List** pane to examine the details of each peak. Peaks derived from the $K\alpha_2$ wavelength are shown with gray background.

L	ists F	ane									×
		Quantification	Rel	⁻ inement Ca	ntrol		0	bjec	t Inspector		
	F	Pattern List	Scan List	Pea	ık List		And	hor	Scan Data		
	No	Pos. [*2Th.]	FWHM [°2Th.]	Area calc.	Assignmen	t h	k	1	Multiplicity	Fo	
	► T	1 23.07813	0.066912	0.0000							
ľ		2 24.50766	0.117096	0.0000							
ľ		3 28.27435	0.102	0.0000							
ľ		4 28.36265	0.0408	0.0000							
ľ		5 29.40404	0.0612	0.0000						\square	
ľ		5 29.49417	0.0408	0.0000							
ľ		7 31.42904	0.0816	0.0000							
ľ		33.60781	0.0816	0.0000							
ľ		35.97685	0.0612	0.0000							
ľ	1	36.20969	0.0816	0.0000						\square	
ľ	1	1 39.41504	0.0612	0.0000							
ľ	1	2 39.75338	0.0612	0.0000							
ľ	1	3 41.49192	0.0612	0.0000							
ľ	1	4 41.60407	0.0612	0.0000							
ľ	1	5 43.1655	0.0816	0.0000						\square	
ľ	1	6 43.28667	0.0612	0.0000							
ľ	1	7 44.21186	0.0816	0.0000							
ľ	1	47.00898	0.0816	0.0000							
ľ	1	9 47.14141	0.0612	0.0000							
ľ	2	47.50092	0.0816	0.0000							
ľ	2	1 47.63182	0.0612	0.0000							
ľ	2	2 48.50757	0.0816	0.0000							
ľ	2	3 48.64017	0.0612	0.0000							
ľ	2	4 50.22773	0.0816	0.0000							
ľ	2	50.36435	0.0612	0.0000							
	2	54.85233	0.0816	0.0000							
	2	7 55.00677	0.0816	0.0000							
	2	3 55.76505	0.0816	0.0000							
	2	9 55.91818	0.0612	0.0000							
1						_	_	_		_	

Figure 4.7 Peak List pane

- 9. Save the document:
 - a. On the menu bar, go to **File > Save As**.
 - b. Use these settings:

File name	Mixture3
Save as type	HighScore Plus (*.HPF)

- c. Click Save.
- d. If the Confirm Save As window opens, click Yes.



Figure 4.8 Confirm Save As window

The document is saved with the background data and the peak data.



CHAPTER 5 DO AN AUTOMATIC PROFILE FITTING

5.1 Introduction

In this chapter, you will fit a profile with an automatic fit parameter set. This is called non-phase peak fitting or single peaks profile fitting.

To do a non-phase peak fitting is to put a calculated profile around each peak and make each peak follow the original measured scan data as close as possible. This mathematical description of the profile shows peak parameters much better than a simple peak search. It is used for peak deconvolution, unit cell refinements, line profile analysis or structure solution.

For phase identification, it is usually not necessary to do a non-phase peak fitting.

5.2 Prepare for an automatic profile fitting

- 1. If the software is not started, start it. Refer to Section 2.2.
- On the menu bar, go to File > Open to open the document "Mixture3.hpf" that you saved. Refer to Section 4.3. "Mixture3.hpf" contains background data, peak data and profile data, but the profile data is calculated from the peak data and not fitted and not refined.
- 3. In the **Additional Graphics** pane, show the differences between measured scan data and profile data:
 - a. Right-click in the **Additional Graphics** pane.
 - b. From the pop-up menu, go to Show Graphics.
 - c. Select Difference Plot.
- 4. In the Main Graphics pane, show the calculated sum profile in Analyze View:
 - On the **Display Mode** toolbar, click A to switch on the **Show Calculated Profile in Analyze View** function.

<u>V</u> iew Treat <u>m</u> ent	Reference <u>P</u> atterns	<u>A</u> nalysis	<u>R</u> eports	<u>T</u> ools	<u>C</u> ustomize
]: 27.164	d-spacing [Å]: 3.20	302	Counts	8	÷
ilii • 🛌 🗛 A	1 🎄 • 🌆 🗢 🕂	AL 14	🖬 🕶 📊		<u> </u>
Mixture3 🖀 Mixture	es 🎽 🎦 🤌 💈	how Calcu	ilated Prof	ile in Ar	ialyze View

Figure 5.1 Show Calculated Profile in Analyze View button

 Alternatively, on the menu bar, go to View > Display Mode and select Show Calculated Profile.

The the calculated sum profile shows in the **Analyze View** in the **Main Graphics** pane.

5. On the **Desktop** toolbar, set the desktop layout to **Profile Fitting**.

								15	
ntIdeAll	i 🖕 i 🏋	Å	<u>^</u>	<u>^ - 213 - </u> 2	r il il i	🖺 🕸 🧿	🖕 🧯 Profile Fi	tting 🔽 🛃 🕽	k -
		∎						Desktop N	lame 🕨 🕨
			41	68.6642	1.36580	1488,83	0.0816	0.600	
			40	65,5939	1,42210	265.11	0.0816	0.600	
			- 39	65.3087	1,43116	456.24	0.0816	0.600	
			38	65.1205	1,43129	919.43	0.1020	0.600	
		-							

Figure 5.2 Select the desktop layout

- 6. Make sure that the background is used and not changed during profile fitting:
 - a. Go to **Refinement Control** pane.
 - b. Select **Background** to open the **Object Inspector** pane.
 - c. In the row of **Method**, click the arrow at the end to open the drop-down list.
 - d. Select Use available background.

Object Inspector	= ×
Selected object: Glo	bal Settings
Background	
Method	Use avail 🕋
Use Extended Ba	ackgrcPolynomial
Flat Background	Basepoints
1/X Background	Chebyshev I
Variables	Amorphous Sinc Function Damped Amorphous Sinc Function
🖻 Agreement India	ces
R expected	0
P profile	0

Figure 5.3 Use available background

- e. Press **Enter** to save the change.
- 7. Set the fitting mode to **Automatic**:
 - On the Fitting toolbar, select Automatic.

<u>W</u> indow <u>H</u> elp						
: 🗋 🤌 💭 🗋	🍐 🖻 🖑	*06	10 • @	•		本法父友的
		: 🐉 🏦 🔯	10 • 🕅	Automatic	N 7	0 Parameter(s) varied
					Selec	t Profile Fit Mode

Figure 5.4 Select the fitting mode

 Alternatively, on the menu bar, go to Analysis > Fitting and set Fitting Mode to Automatic.

5.3 Start an automatic profile fitting

- 1. Make sure that you prepared for the automatic profile fitting correctly. Refer to Section 5.2.
- 2. Make sure that the profile fitting will be done to the full range of the scan:
 - Fully zoom out the scan in the **Main Graphics** pane. Profile fitting is only done to the part of the range that is shown.



Figure 5.5 Ignore actual Zoom Range for fitting button

- 3. Start an automatic profile fitting:
 - On the **Fitting** toolbar, click **Profile fit> Default Profile Fit**.



Figure 5.6 Select < Profit fit> Default Profile Fit

 Alternatively, on the menu bar, go to Analysis > Fitting > Start Fit and select <Profile fit> Default Profile Fit.

The profile fitting starts. The **R-Value** window opens.





4. Examine how the overall quality of the fitting changes during the fitting process.

After the profile fitting is completed, the profile is fitted to the peaks. The background is used and not changed. The background is high and cuts into the peak feet. This is good for phase identification, but is not optimal for profile fitting.



Figure 5.8 Anchor scan and profile after the first profile fitting

- 5. Examine the agreement indices:
 - a. Go to the **Refinement Control** pane.
 - b. Double-click Global Variables to open the Object Inspector pane.
 - c. Open Agreement Indices.
 - d. Examine the R-values and the value of **Goodness of Fit**.

Re	efinemer	nt Control						Ob	ject Inspector	•	X
	Quantif	fication And	nor Sc	an Data	Pattern	List Peak List	s I	Se	lected object: Global Settings		
ſ	Refi	nement Control	3	X Sca	in List	Structure Plot	4		Background		
	Name	_	In	Refine	Value	Deviation	C		Method	Use avail	
3	THG.	Global Variable	3	The first	T BTOTE	Denation	-		Agreement Indices		
							-		R expected	4.75012	2
									R profile	6.38713	3
									Weighted R profile	9.31467	7
									D-statistics	0.62428	3
									Weighted D-statistics	0	5
									Goodness of Fit	1.96093	3
									Mixture MAC [cm^2/g]	0.00	3

Figure 5.9 Agreement Indices after the first profile fitting

- 6. Change the settings which are used to find the background:
 - a. On the menu bar, go to **Treatment > Determine Background**.
 - b. In the **Determine Background** window, change these settings:

Bending factor	0
Granularity	30
Use smoothed input data	Not selected

Automatic	Manual	By Search <u>P</u> eaks	Subtract
After Sonn	neveld & I	lisser	Save to list
B <u>e</u> nding fa	actor:	0	Net Scan
			Background
<u>G</u> ranularity	y:	30	Accept
	-0		Close

Figure 5.10 New settings

- c. Click **Accept** to save the new settings.
- 7. Set the asymmetry type for peak fitting:
 - a. Go to the **Refinement Control** pane.
 - b. Double-click **Global Variables** to open the **Object Inspector** pane.
 - c. Open Unassigned Peaks Fitting.
 - d. Open Common Peak Fit Settings.
 - e. In the row of **Asymmetry Type**, click the arrow at the end to open the drop-down list.
 - f. Select **Split Width**. This shows the small asymmetry of the peaks at low angles better.

elected object: Global Settings		
Background		\$
Method	Use available background	
Agreement Indices		\$
R expected		5.0388
R profile		12,1180
Weighted R profile		17.1388
D-statistics		0.4115
Weighted D-statistics		1
Goodness of Fit		3,4013
General Fit Properties		\$
Job Type	X-rays	
Weighting Scheme	Against Iobs	
Solver Time Limit [sec]		300
Solver Iteration Limit		100
Solver Nu		
Solver Tolerance		0.00
Max. No of Fit Cycles		2
Peak Base Width for Fit		2
Automatic Cell Constraints		
Automatic Anisotropic Displacement		
Keep ADP's positive definite		
Automatic Atom XYZ Constraints	V	
Specimen Displacement [mm]		
Zero Shift [º20]		-0.0092
Wavelength [Å]		1,540
K-o2 / K-o1 Intensity Ratio		0.
K-β / K-o1 Intensity Ratio		9
Polarisation Correction Coefficient		
Use Brindley Microabsorption Correct		
Calculate Errors		
Unassigned Peaks Fitting		*
Common Peak Fit Settings		\$
Profile Function	Pseudo Voigt	
Use Caglioti Function		
Use Shape Function		
Asymmetry Type	Split Width	
Caglioti Width	No Asymmetry Function	
Peak Shape	Split Width Split Shape Split Width and Shape	
Asymmetry	Finger, Cox, Jephcoat	

Figure 5.11 Select Spit Width

- g. Press **Enter** to save the change.
- 8. Do step 3 to start an automatic profile fitting again with **<Profile fit> Default Profile Fit**. The profile fitting becomes better this time.



Figure 5.12 Anchor scan and profile with better fitting

- 9. Examine the agreement indices:
 - a. Go to the **Refinement Control** pane.
 - b. Double-click Global Variables to open the Object Inspector pane.
 - c. Open Agreement Indices.
 - d. Examine the values of Weighted R profile and Goodness of Fit:
 - The value of **Weighted R profile** must be approximately 6.80.
 - The value of **Goodness of Fit** must be approximately 1.44.

Refinemen	nt Control					0	bject Inspector	= >
Quantif	ication Anch	or Sca	an Data	Pattern Lis	t Peak List	s	elected object: Global Settings	
Refi	nement Control)	< Sca	n List St	ructure Plot		Background	
Name		In	Refine	Value	Deviation	c	Method	Use avail
> B B	Global Variables			March The A	1200001200-001	10	Agreement Indices	
-	Zero Shif		101.	0	0.000000		R expected	4.72034
	Specime			0	0.000000		R profile Weighted D are file	5.02143
	K-α2 / K			0.5	0.000000		D-statistics	0.9923
	К-β / К-α		10%	0	0.000000		Weighted D-statistics	0
+	🗼 Unassign	1	11				Goodness of Fit	1.44041
		1					Mixture MAC [cm^2/a]	0.00

Figure 5.13 Agreement Indices with better fitting

10. Click ➤ or go to **File > Close** on the menu bar to close the document. Do not save the changes.



CHAPTER 6 DO SEARCH - MATCH - IDENTIFY

6.1 Introduction

In this chapter, you will search and match possible candidates and then manually identify the phases of the example document. You will use PANalytical Example Database, which is supplied with the software. This gives you a good example of what to do in real work.

NOTE: Do not use PANalytical Example Database to do a test of the functionality or examine the phase identification capabilities of the software. Use a large reference database with at least 100,000 patterns instead, for example an ICDD product or the free COD database.

6.2 Search and match

- 1. If the software is not started, start it. Refer to Section 2.2.
- 2. On the menu bar, go to **File > Open** to open the document "Mixture3.hpf" that you saved. Refer to Section 4.3. "Mixture3.hpf" contains background data, peak data and profile data, but the profile data is calculated from the peak data and not fitted and not refined.
- 3. On the **Desktop** toolbar, set the desktop to **Phase-ID**.



Figure 6.1 Set the desktop layout

- On the menu bar, go to Analysis > Search & Match > Execute Search & Match. The Search & Match window opens.
- 5. Read the title bar of the **Search & Match** window. The title bar shows the parameter set that you use.
- 6. Make sure that you use the **Default** parameter set.

Search & Match - [Default]	×
Restrictions Parameters Automatic	
Data source: Peak & Profile Data 💙	Search
Scgring scheme: O Single phase O Multi phase	
 ✓ Auto residue ✓ Demote unmatched strong ✓ Match intensity ✓ Allow pattern shift 	ОК
Known T <u>w</u> o Theta shift [°2Th.]: 0	Cancel More >>

Figure 6.2 Use the Default parameter set

- 7. If not, change the parameter set to **Default**:
 - a. Click **More** to expand the window.
 - b. In the Select Parameter Set field, select Default from the drop-down list.
- 8. In the **Search & Match** window, click **Search**. The scan shows in the **Additional Graphics** pane in the **Compare Mode** view.
- 9. Go to the **Pattern list** pane. The **Candidates** list in the lower part of the pane gives a preview of the candidates.

5	Selected Candidate: 00-038-1479							
	No.	Ref. Code	🐮 S	Compound Name	Chemical Formula	Scale F N		
۲	1	ICOD 00-038-1479	71	Chromium Oxide	Cr2 03	0.157		
	2	ICOD 00-005-0586	64	Calcium Carbonate	Ca C O3	0.375		
	3	ICOD 00-035-0816	60	Calcium Fluoride	Ca F2	0.996		
	4	ICOD 01-077-2041	46	Sodium Erbium Flu	Na Er F4	0.922		
	5	ICOD 00-006-0329	31	Praseodymium Oxide	Pr O1.83	0.806		
	6	ICOD 01-075-0134	25	Uranium Oxide	U O2	0.579		
	7	ICOD 01-073-1667	22	Copper Iron Sulfide	Cu5 Fe S4	0.484		
	8	ICOD 00-027-1402	17	Silicon	Si	0.054		
	9	ICOD 00-033-1161	1	Silicon Oxide	Si O2	0.333		
	10	ICOD 00-046-1045	1	Silicon Oxide	Si O2	0.358		
	11	ICOD 00-006-0694	1	Chromium	Cr	0.008		
4						•		

Figure 6.3 Candidates list

- 10. If necessary, click the column header **Score** to change the sequence of the candidates by score.
- 11. In the Search & Match window, click OK to accept the result.

6.3 Identify

In this section, you will manually accept candidates that have high scores and that match the peaks and features of the measurement. Some views in the **Additional Graphics** pane support a visual comparison of reference pattern lines and the measurement.

- 1. Set the Additional Graphics pane to the Compare Mode view:
 - On the menu bar, go to View > Additional Graphics and select Compare Mode.
 - Alternatively, right-click in the **Additional Graphics** pane, from the pop-up menu, go to **Show Graphics** and select **Compare Mode**.
- 2. Set the display of peaks to Show Peaks Outside:
 - On the **Display Mode** toolbar, click the small arrow next to **und** and select **Show Peaks Outside**.

<u>F</u> ile	<u>E</u> dit	⊻iew	Treat <u>m</u> ent	Reference <u>P</u> atterns	<u>A</u> nalysis	<u>R</u> eports	<u>T</u> ools
Pos	. [°20]:			d-spacing [Å]:		Counts	:
	ЪЛ,	dii	S 14 1	n Au – An 😂 An	A2 A6 1	🕰 👻 📊	
	% a.	4	Show Peaks	Inside and Outside			
	B		Show Peaks	Inside			
S O		1	Show Peaks	Outside		Y	Y
1	Cou		Hide Peaks				

Figure 6.4 Select Show Peaks Outside

- Alternatively, on the menu bar, go to View > Display Mode > Peaks in Main Graphics and select Show Peaks Outside.
- 3. Hide the calculated profile to show the measured scan in a better view:
 - Go to to View > Display Mode and make sure that Show Calculated Profile is not selected.
 - Alternatively, make sure that \overline{M} on the **Display Mode** toolbar is not selected.

<u>V</u> iew Treat <u>m</u> ent	Reference <u>P</u> atterns	<u>A</u> nalysis <u>R</u> eports	<u>T</u> ools <u>C</u> ustomize
]: 27.164	d-spacing [Å]: 3.2	802 Counts	51 -
- Ilii 🗝 🛌 🗛 🗛	1 🌆 • 🌆 🗢 🕂	A 🗛 🕶 🚃	
Mixture3 🛗 Mixture	83 🅦 🎦 🎽	Show Calculated Prof	ile in Analyze View

Figure 6.5 Show Calculated Profile in Analyze View button

- 4. Go to the Pattern List pane.
- 5. Drag the first candidate pattern "00-038-1479, Chromium Oxide" from the Candidates list to the Accepted Ref. Patterns list to accept it. The Accepted Ref. Patterns list is above the Candidates list. When you select this pattern, it is highlighted and its lines are shown for comparison in the Additional Graphics pane.
- 6. Examine these changes in the display panes:
 - Some peaks in the **Main Graphics** pane lost the **V** mark. These peaks are explained by the accepted reference pattern. Peaks marked with **V** marks are still not explained.
 - High intensity areas in the **Main Graphics** pane are shown in gray. These scan features are explained by the accepted reference pattern.

NOTE: If no scan features have gray marks, click M in the **Display Mode** toolbar to switch on the **Show Explained Features** function.



Figure 6.6 Show Explained Features button

• In the **Candidates** list, the candidate "00-005-0586, Calcium Carbonate" moves to the top. The score of this pattern has a small difference than before.

NOTE: If necessary, go to **Edit > Undo** and **Edit > Redo** to make these changes occur again to examine them.

- 7. Drag the top 2 patterns from the **Candidates** list one by one to the **Accepted Ref. Patterns** list to accept them. When you accept the third reference pattern, examine the large changes of the scores of the remaining candidates at the same time.
- 8. Examine if the accepted patterns are minerals and its subfile information:
 - In the Accepted Ref. Patterns list, examine the information in the Subfiles column.
 - **NOTE:** When the information is not fully shown, you can hover the cursor over the cell to see the full information.

.ists Pane							×	
Quantification	Quantification Re			ntrol		Object Inspector		
Pattern List	Scan List Peak List		Anchor Scan Data					
Accepted Ref. Pattern: 00-035-0816								
Compound Name	Chemical	Score	Se	Display Co	lor Dat	abase ID	Subfiles	
Chromium Oxide	Cr2 03	72	13	Blue	C:V	Docume	Alloy, metal or	
Calcium Fluoride	Ca F2	63	60	Lime	C:V	Docume	Nommon Phas	
Calcium Carbonate	Ca C O3	62	27	Gray	/ C:W	Docume	🔃 Common Ph	
						L.		

Figure 6.7 Complete information of the cell

• Alternatively, right-click a reference pattern in the **Accepted Ref. Patterns** list and from the pop-up window, select **Show Pattern** to show the subfile information.

Reference Pattern: 00-038-147	9	x
Name and formula		^
Reference code:	<u></u> ጋዐ-038-1479	≡
Mineral name: Compound name: Common name: PDF index name:	Eskolaite, syn Chromium Oxide chrome green Chromium Oxide	
Empirical formula: Chemical formula:	Cr ₂ O ₃ Cr ₂ O ₃	
<u>Crystallographic param</u>	<u>eters</u>	
Crystal system: Space group: Space group number:	Rhombohedral R-3c 167	
a (Å): b (Å): c (Å): Alpha (°): Beta (°):	4.9588 4.9588 13.5942 90.0000 90.0000	
Gamma (°):	120.0000	•
< > <u>S</u> ave As <u>C</u> opy	Print <u>G</u> raphics Print All	Intensity Scale 🔻 Angle Sc

Figure 6.8 Reference Pattern window

9. In the **Display Mode** toolbar, click M to switch on the **Show Reference Patterns** function. The reference pattern lines show in the **Main Graphics** pane.

<u>F</u> ile	<u>E</u> dit	<u>V</u> iew	Treat <u>r</u>	<u>n</u> ent	Reference <u>P</u> a	tterns	<u>A</u> nalysis	<u>R</u> eports	<u>T</u> ools
Pos	. [°28]	: 26.643			d-spacing [/	Å]: 3.34	431	Counts	:
•	W.	ilit •	~ J	N 14	. Aa - 🗛 🤇	3 M	赵山	🖬 👻 📊	
				15					
	ا 😭	Mixture3	8 🚹	Show	/ Reference Pa	tterns			

Figure 6.9 Show Reference Patterns button

For this time, these peaks or features cannot be explained:

- A small peak around 42 °2theta is not explained, but this can be a real peak or just some noise.
- A second unexplained peak around 79.3 °2theta is a K α_2 peak, which is incorrectly assigned as a K α_1 peak. You can go to **Tools > Spectral Lines** to do a check of it.



Figure 6.10 Unexplained peaks or features

- 10. Save the complete document:
 - a. Go to File > Save As.
 - b. Use these settings:

File name	Mixture3
Save as type	HighScore Plus (*.HPF)

- c. Click **Save**.
- d. If the Confirm Save As window opens, click Yes.



Figure 6.11 Confirm Save As window

The document is saved with background data, peak data, reference patterns and a candidate list.



CHAPTER 7 CHANGE SCORES

7.1 Introduction

In this chapter, you will change the scores of the search and match results in the **Candidates** list.

The scores, shown in the **Candidates** list and the **Accepted Ref. Patterns** list, are related to the parameter sets used for the search and match procedures. These scores can also be changed at any time after the search and match. In real work, if you change scores, other candidates can move to the top of the **Candidates** list and this can help you find more phases.

7.2 Change scores

- 1. If the software is not started, start it. Refer to Section 2.2.
- On the menu bar, go to File > Open to open the document "Mixture3.hpf" that you saved. Refer to Section 6.3. "Mixture3.hpf" contains background data, peak data, reference patterns and a candidate list.
- 3. Drag the reference pattern "00-035-0816, Calcium Fluoride" from the **Accepted Ref. Patterns** list to the **Candidates** list.
- 4. Make sure that the **Pattern** toolbar is shown:
 - a. Go to **View > Toolbars**.
 - b. Select Pattern Toolbar.





5. Click ^{the} to change the data source and examine the changes of the score of the Fluorite reference pattern at the same time.



Figure 7.2 Click Data Source button

NOTE: The look of the **Select Data Source** button is related to the selected data source:

- 쓴: Profile
- 🛄: Peak List
- 🏠 : Both
- 6. Click $\frac{1}{2}$ to include or exclude the quality of matching relative intensity in the score and examine the changes of the scores at the same time.

NOTE: Steps 5 and 6 are frequently used to change the scores of a **Candidates** list.

7. Click 🕍 to change the scoring scheme and examine the changes of the scores at the same time.



Figure 7.3 Click Scoring Scheme button

- **NOTE:** The look of the **Select Scoring Scheme** button is related to the selected scoring scheme:
 - 🛓: Single Phase
 - ៉: Multi Phase
- 8. Click to switch on and off the pattern shift and examine the changes of the scores and scale factors at the same time.
 - **NOTE:** Steps 7 and 8 are almost never used to change the scores. Usually, the scoring scheme is set to **Multi Phase** and the pattern shift is switched off.
- 9. Click ▼ or go to **File > Close** on the menu bar to close the document. Do not save the changes.



CHAPTER 8 USE A USER BATCH

8.1 Introduction

All actions that you did about pattern treatments and search, match and identification can be done automatically with pre-programmed pattern restrictions. The default user batches in the software have pre-programmed pattern restrictions. With user batches, you can do a full analysis with just a click of a button.

Your knowledge about the sample is the most powerful tool that you have for phase identification. On the **Restrictions** tab of the **Search & Match** window, you can use some reference patterns for 1 or more conditions which will be used for search and match.

In this example, you will not use any restrictions, but search fully through the small PANalytical Example Database.

8.2 Use a user batch

- 1. If the software is not started, start it. Refer to Section 2.2.
- On the menu bar, go to File > Open to open the document "Mixture3.xrdml". Refer to Section 3.2.

NOTE: Be careful that you do not open the document "Mixture3.hpf".

- 3. Start the pre-programmed batch "IdeALL":
 - On the **Batches** toolbar, click **IdeAll**.

🖭 ClipAllToZoom 🖭 Default 🛄 IdeAll 💷 IdeCom 🖭 IdeMin 🖭 Merge PDF scans 🖭 MinorMinerals 🖭 MultiRiet 🖭 Overlay Scans 🖭 PrintIdeAll 🥃

Figure 8.1 User batch IdeAll

• Alternatively, on the menu bar, go to **Tools** and select **IdeAll**.

Table 8.1 What IdeALL does

Sequence	Step
1	Find the background.
2	Search peaks.
3	Convert divergence slit from automatic to fixed divergence slit intensities. NOTE: This step is not used in this example, because the measurement "Mixture3.xrdml" was done with fixed divergence slit. Batch step not executed. For more information please have a look at the process log. <i>Figure 8.2 A batch step is not used</i>
4	Use the full PANalytical Example Database without restrictions for search and match.
5	Automatically identify candidates that have high scores.

- 4. After the operation of the user batch is completed, examine the peaks or features that cannot be explained in the **Main Graphics** pane.
- 5. Examine if the accepted patterns are minerals and its subfile information:
 - Go to the **Pattern List** pane and in the **Accepted Ref. Patterns** list, examine the information in the **Subfiles** column.

NOTE: When the information is not fully shown, you can hover the cursor over the cell to see the full information.

- Alternatively, right-click a reference pattern in the **Accepted Ref. Patterns** list and from the pop-up window, select **Show Pattern** to show the subfile information.
- 6. Examine the details of the batch:
 - a. Go to File > Properties to open the Properties of Mixture 3 window.
 - b. Go to the **Process Log** tab.

Process log shows that a batch step is not used.



Figure 8.3 This batch step is not used

- 7. Save the document:
 - a. On the menu bar, go to File > Save Document.
 - b. Use these settings:

File name	MixtureBatch
Save as type	HighScore Plus (*.HPF)

c. Click **Save**.

The document is saved with the identified phases.



CHAPTER 9 PHASE IDENTIFICATION STRATEGY AND TROUBLESHOOTING

9.1 Introduction

There is no recipe for phase identification. However, this phase identification strategy can give you some simplified and schematic steps on how to identify an unknown sample in the software.

NOTE: In the identification examples, at the start you alway search the full reference database with no restrictions. This is done to show the capabilities of HighScore in the 'worst scenario', but it is not necessarily the best approach to all phase identification problems.



Figure 9.1 Phase identification strategy

Table 9.1 Description of the strategy steps

Step	Description
Start	Load a measurement.
Pattern Treatments	 Find the background. This helps before you search peaks and it is very important when you use profile data for identification. Search peaks with a high significance. Do not try to use very small peaks. Also use profile data for input. Convert intensities to fixed slit intensities when an automatic theta-compensating divergence slit was used. This step is optional, but it makes measured intensities match the reference data better.
	• Strip $K\alpha_2$. This step is optional. When you do not use the Strip K Alpha2 function($\frac{l}{M}$), this step is done by the search and match algorithm. The process is done in the background and not shown to you. Most users find it is better not to see the process because then they can see the original measured data in the graphics.
Search and Match	Start search and match to get a new Candidates list.
Identification	 Use score and scale factor values to identify and accept candidates. Do a check of the graphics. Examine the lists for more textual information.
Ide Batch Programs	 You can use the user batch programs with names that start with "Ide" to do the steps above automatically.
Change Range/ Peaks Included	• Examine the unexplained peaks or features with the Track Graphics Range function (A) or exclude the explained peaks and start search and match again to get a new Candidates list.
Change Restrictions	 Change restrictions. It is possible that the restrictions that you used were too tight and excluded some phases in the sample, or they were too wide and included many isotypical patterns and made the Candidates list full.
Change scores, group candidates	 Change scoring parameters to put other patterns to the top of the Candidates list, or put the candidates together to make similar patterns into a group. The idea is to identify as many phases as possible from the Candidates list before you start search and match again.
Select standard	Select 1 identified pattern with simple, fixed chemistry as an internal standard.
All Correctly Explained?	 Do a check of all the patterns that were identified automatically. An automatic identification is not always correct.
End	Save your result as a diffraction document in HDF format.

It is always good to include all information about the sample that you have. Use the subfiles made by ICDD or give special restrictions for the samples and analytical problems that occur to you.

9.2 Pattern treatment sequence

9.2.1 Find background

You must find the background before you search peaks. When there is no background data, peak search automatically makes its own background data.

When profile data is used for identification, it is very important to find the background correctly. If you are not sure, usually for phase identification, it is better to have the background too high than too low.

9.2.2 Search peaks with a high significance

Do not try to find all small peaks. When you select profile data for phase identification, all intensity above the background is used as input, which also includes very small peaks that are difficult to find.

9.2.3 Convert intensity

Convert intensity to fixed slit intensity when an automatic theta-compensating divergence slit was used. This is optional. The converted intensity will match the reference data better, which gives better scores when the **Match Intensity** function (^{±±}) is used.

9.2.4 Strip Ka₂

If you do not use the **Strip K Alpha2** function (\swarrow) to strip Ka₂, Ka₂ stripping is done by the search and match algorithm. The process is done in the background and not shown to you. Most users find it is better not to see the process because then they can see the original measured data in the graphics.

9.3 Identify

Some tools are available to get information about the match of a candidate or accepted pattern:

- For information about the explained scan regions, click **Show Explained Feature** button to switch on the **Show Explained Feature** feature.
- For information about matched peak, do a check of the **V** marks in the **Main Graphics** pane and the **Matched** column in the **Peak List** pane.
 - NOTE: If the Matched column is not shown in the Peak List pane, you can right-click in the Peak List pane, select Customize Peak List from the pop-up menu to open the Customization window, and then double-click Matched to add it as a column in the pane.
- For information about matched peaks by reference codes, do a check of the **Matched by** column in the **Peak List** pane.
- For information about the matched reference pattern lines, right-click in the **Pattern List** pane and select **Analyze Pattern Lines** from the pop-up menu.
- For information about the number of matched lines, do a check of the the **Matched Lines** column in the **Accepted Ref. Patterns** list in the **Pattern List** pane.

When you cannot identify all phases from the **Candidates** list of the initial search and match, do as follows:

- 1. Change the scores to find more candidates from the **Candidates** list.
 - Change the data source with the Change Data Source button 4, 11 or 4 on the Pattern toolbar.
 - Match the parameter with the **Match Intensity** button $rac{11}{200}$ on the **Pattern** toolbar.
 - When you look for minor phases, do not select the **Demote Unmatched Strong** button A on the **Pattern** toolbar.

- If you do not have strong solid-solution effects or an incorrect sample height, do not select the **Allow Pattern Shift** button in on the **Pattern** toolbar. For the usual small pattern shifts, this parameter is usually more applicable to patterns that do not fit very well to the measurement.
- If necessary, change the scoring scheme with the **Select Scoring Scheme** button ²⁴ and do not select the **Auto Residue** button ²⁶ on the **Pattern** toolbar.
- 2. Start a search and match again with new, different restrictions or different parameters. It is possible that the restrictions that you used were too tight and excluded some phases in the sample, or they were too wide and included many isotypical patterns and made the **Candidates** list full.
- 3. If necessary, do step 1 again with the new **Candidates** list. This can solve many identification problems.
- 4. Optionally, when you identified 1 or several phases , but cannot identify more components in a mixture, this can be caused by a sample height or an alignment problem. If this occurs, it helps to select an identified pattern with a well defined, stable chemistry and use it as an internal standard:
 - a. Right-click in the Accepted Ref. Pattern list.
 - b. From the pop up menu, select **Correct Scan with Pattern**. The measurement with peaks and background is then shifted to fit optimally to the standard selected pattern.
 - **NOTE:** Be careful when you do this step. When you select a pattern with variable chemistry (= solid solutions possible) as internal standard, it can have a bad effect on the correct identification of more phases.

Make sure that the **Allow Pattern Shift** button is not selected on the **Pattern** toolbar when you select an internal standard.

- 5. If you still cannot identify all phases from the **Candidates** list, do as follows:
 - a. Switch on the **Track Graphics Range** function with \Lambda on the **Pattern** toolbar.
 - b. Zoom in on the low-angle region which contains peaks and features that are not explained and examine the **Candidates** list again. For this time, only the peaks and features in the shown, zoomed-in region are used as input.
 - c. If there are no new matches found in the top of the **Candidates** list, start a new search and match only with the zoomed-in range as input.
 - d. Alternatively, exclude all matched peaks and do a search and match again with only the remaining peak data as input.
 - e. If necessary, do steps 1-3 again with the new Candidates list.

9.4 Troubleshooting

If you still cannot identify all phases from the **Candidates** list or get very bad result, you can do troubleshooting:

- 1. Examine the background. A background level that was found too low has an effect on each reference pattern which match your measurement.
- 2. Examine the document wavelength, which is usually derived from the anchor scan. If the document wavelength is not in the anchor data, it is supplied by the default instrument settings and cannot match the actual scan data.

- 3. Examine which databases are used for search and match. Make sure there is a reference pattern in these databases applicable to your measurement.
 - **NOTE:** All patterns in the pattern list are treated as 'known' when you start search and match. They are kept in the pattern list and the first 20 are scored. When the **Auto Residue** function (A) is used, these first 20 patterns also have an effect on the scores of all candidates.



CHAPTER 10 SEARCH AND REFINE A UNIT CELL (HIGHSCORE PLUS)

10.1 Introduction

In this chapter, you will do these tasks:

- a. Load a measurement, set the correct wavelength, and convert the document format from RFL to RD, a format in which the data about the used wavelength can be saved.
- b. Search diffraction peaks. The first 8 peaks (reflections) are used to start a unit cell search, which is also called indexing, with the TREOR and DICVOL indexing routines.
- c. Refine a good cell candidate with all diffraction peaks as input.
- d. Save the results.

10.2 load a measurement

- 1. If HighScore Plus is not started, start it. Refer to Section 2.2.
- 2. On the **Desktop** toolbar, set the desktop layout to **Structures**.



Figure 10.1 Set desktop layout

- 3. On the menu bar, go to File > Open to open the document "TaSSE.rfl" in this folder C:\Users \user.name\Documents\PANalytical\X'Pert HighScore Plus\Tutorial.
- 4. Set the the wavelength for the data set to Copper $\ensuremath{\text{K}}\ensuremath{\alpha}_1$ only:
 - a. Go to the **Scan List** pane.
 - b. Double-click in the list to open the **Object Inspector** pane.
 - c. Open Instrument Settings.
 - d. Select Incident Beam Monochromator.

Lists Pane				×	0	bject Inspector	×
Quantifica	ation	Ref	inement Control		5	elected object: Scan(s)	
Structure Plot	Fourie	r Map	Distances and Angles	;		Scan Display	۲
Pattern List	Scan List	PeakList	Anchor Scan Dat	a	r	General Scan Info	*
No	Visibl	e	Name	St	r	Scan Statistics	۲
	1	~ ~	TaSSe		h	Peak Statistics	*
		_			h	Instrument Settings	*
					h	Incident Beam Monochromat	
						Spinner used	
						Mode Linear Detector	None
						Length Linear Detector [°2Th	2
						Anode Material	Copper (C
						Tube Current [mA]	0
						Tube Tension [kV]	0
						Divergence Slit Type	Fixed
						Fixed Div. Slit Size [°]	1

Figure 10.2 Select Incident Beam Monochromator

5. Go to **File > Save As** to save the document as a PHILIPS binary scan in the RD format. The wavelength with the scan data is saved in this format.

By this time, the important instrument information is corrected and the scan is converted from RFL file format to RD file format.

10.3 Search peaks

- 1. On the menu bar, go to **Treatment > Search Peaks**.
- 2. Click More to expand the window.
- 3. In the **Select Parameter Set** field, select **Default** from the drop-down list. The title bar of the **Search Peaks** window shows the name of the parameter set that is used.

Search Peaks	[Default]			×
Mi <u>n</u> imum signific	ance:	1.	.00	Search Peaks
Minim <u>u</u> m tip wid	th [°2Th.]:	0.	.01	Accept
Maximum tip <u>w</u> id	dth [°2Th.]:	1.	.00	
Pea <u>k</u> base widtł	n [°2Th.]:	2	.00	
Met <u>h</u> od:	Minimum 2nd d	lerivative	~	Close
Trijal:			~	Less <<
Select Paramete	er Set			
Default		× 6		(🖻 🔌 🗶

Figure 10.3 Use the Default parameter set

- 4. In the Search Peaks window, click Search Peaks to search peaks.
- 5. Click **Accept** to accept the results. The **Main Graphics** pane shows the peaks and a theoretical profile.
- 6. Zoom in around 60 to 65 °2theta. You can see that a peak is incorrectly found at about 61 °2theta.



Figure 10.4 Peak incorrectly found

- 7. Hover the cursor over the peak to see detailed information.
- 8. On the menu bar, go to View and select Peak List Pane.
- 9. Find the only peak that is not correctly at its place: No. 17. This peak is selected when you hover the cursor over it.

- 10. Delete the peak:
 - Press **Delete**.
 - Alternatively, in the **Peak List** pane, right-click in the row of the peak and from the pop-up menu, select **Delete Peak**.

10.4 Search and refine a unit cell

- 1. On the menu bar, go to **Analysis > Crystallography > Search Unit Cell**. The **Search Unit Cell** window opens.
- 2. Go to **General**.
- 3. Set Indexing Method to Treor.

Sea	rch Unit Cell - [Default]				×
E	Peak Parameters		Execute Cell Search	Selected Cell Can	didate
	Minimum Intensity [%]	0.1		» [8]	
	Use the x first peaks	10	Show Detailed Output		
E	Cell Parameters			U[A]	
	FOM better than	12	Show Existing Results	C [A]	
	Maximum Beta [°]	130		Alpha [°]	
	Maximum Axis [Å]	25		Beta [°]	
	Maximum Volume [Å^3]	1000		Gamma [°]	
-1	Crustal System	1000		Volume [Ä^3]	
	Tech Manual Comer Avia	D		Indexing Method	
	Test Monoci, Super Axis	•		FOM	ОК
	Include Monoclinic			Not Indexed	
	Include Triclinic				Close
	General			, ,	
	Indexing Method	Treor			<u>M</u> ore >>

Figure 10.5 Set Indexing Method

- 4. Click **Execute Cell Search**. The **Cell Candidates** window opens. It shows some cell candidates.
- 5. Click Cancel to close the Cell Candidates window.
- 6. In the Search Unit Cell window, set Indexing Method to Dicvol.
- 7. Click **Execute Cell Search**. The **Cell Candidates** window opens. All cell candidates are almost the same. Therefore, there is a very possible unit cell.

(el	l Can	didates										X
		No.	a [Å]	Ь [Å]	c [Å]	Alpha [°]	Beta [°]	Gamma [°]	Cell Volume [Å3]	Indexing Method	Not in	FOM	
	١	1	3.3305	3.3305	21.9013	90.0000	90.0000	90.0000	242.9413	Treor	0	52.0000	
		2	3.3310	3.3310	21.8883	90.0000	90.0000	90.0000	242.8657	Treor	0	44.0000	
		3	3.3310	3.3310	21.8883	90.0000	90.0000	90.0000	242.8657	Treor	0	44.0000	
		4	3.3309	3.3309	21.8875	90.0000	90.0000	90.0000	242.8384	Dicvol	0	39.1000	
		5	21.8559	3.3270	3.3357	90.0000	90.0000	90.0000	242.5581	Dicvol	0	36.7000	
		⊆ор	у [Print	<u>D</u> ele	te selected	Row	Clear <u>T</u> able	<u>R</u> efine Cell		к (Cancel	

Figure 10.6 Cell Candidates window

- 8. Select one of the cell candidates.
- 9. Click Refine Cell. The Refine Unit Cell window opens.

🛣 Refine Unit Cell - [Defau	lt]						
Cell Refinement Calculated an	d Observed Peaks 🛛 Space Group Te	est					
Reflection Conditions		Unit Cell					
Crystal System	Tetragonal	a[Å]	3.3305(8)				
Bravais Type	Primitive (P)	ь [Å]	3.3305(8)				
Space Group		c [Å]	21.90(1)				
Instrument Settings		Alpha [°]	90				
Goniometer Radius [mm]	240.00	Beta [°]	90				
		Gamma [°]	90				
1		Volume [Å^3]	242.94				
		Refinement Results					
Edit <u>P</u> arar	neter Set	2Theta Zero Shift [°]					
Always switch to Calc. and Ob	- Deaks Tab on Definement	Specimen Displacement [mm]					
Miways switch to calc, and ob:	s, reaks tab on Kennement	No. Unindexed Lines					
		No. Indexed Lines					
		Total No. Calculated Lines					
		Chi Square					
		Snyder's FOM					
		Colorb Cont	idata call				
		Select Cand					
Start Refinement	Stop <u>R</u> efinement		OK Cancel				
			.:				

Figure 10.7 Refine Unit Cell window

- 10. Read the title bar of the **Refine Unit Cell** window. The title bar shows the parameter set that you use.
- 11. Make sure that you use the **Default** parameter set.
- 12. Click Start Refinement.
- 13. Go to the **Calculated and Observed Peaks** tab. It shows the calculated and indexed peaks and their deviation from the theoretical values.



Figure 10.8 Cell refine results

- 14. Go to the **Unindexed Peaks** tab. It must be empty for this time.
- 15. Click **OK** to accept the results and close the window.
- 16. Examine the refined unit cell parameters:
 - a. On the menu bar, go to **View** and select **Refinement Control Pane**.
 - b. Open the Search Unit Cell Result 1 phase.
 - c. Open Unit Cell. The refined unit cell parameters is shown.
- 17. Examine the derived data:
 - a. Double-click Search Unit Cell Result to go to the Object Inspector pane.
 - b. Examine the derived data, for example the cell volume and the estimated cell volume error. It also shows the error of the unit cell axis or angle selected in the refinement control.
- 18. When a phase color is accidentally red, change the display color:
 - a. Double-click the phase to open the **Object Inspector** pane.
 - b. Open Phase Display.
 - c. In the row of **Display Color**, click the arrow at the end to open the drop-down list.
 - d. Select a different color.

Lists Pane					X	Object Inspector		×
Structure I	Plot Fourie	r Map 👘 D	istan	ces and	Angles	Selected object: Phase		
Pattern Lis	: Scan List	Anchor Scan Data			Phase Display		_ ی	
Quantification Refin			ement Control			Display Color		-
Name	Ness			Refine	n L Value	Show Phase	Olive	
	ahal Uaviahlar		mio	TTEILITE	value	General Phase Inf	Navy	
	oveb Upit Cell D	ocult 1				Title	Purple	
	ale factor	esult I			0.0001(Use Phase	Teal	
	ale racior oforrod Orioptati	on	0		1.00001	Fitting Mode	Silver	
	overall	on	0		0.00000	Mean Particle Dia	Red	
E F	tinction				0.00000	Standard Weight	Lime	
Ela Ela	at Plate Absorpti	on Correction			0.00000	Pseudo Formula Ma	ISS	-1
Pc	rositu	on concetion			0.00000	Scale Factor		0.0001
B	auabness				0.00000	Overall Displacemer	nt Param	0
- 60	Unit Cell				0.00000	Extinction		0
	Atomic coordir	nates			0 00001	Roughness		0
- A	Profile Variable	35			5.00001	Porosity		0
		-				Flat Plate Absorptio	on Corre	0

Figure 10.9 Display color

e. In the **Main Graphics** pane, examine the peaks. All peaks are assigned to the phase and therefore have the phase color. By this time, there are no non-phase peaks in this example.

10.5 Save results

- 1. Go to File > Save Document.
- 2. Use these settings:

File name	TaSSe
Save as type	HighScore Plus (*.HPF)

- 3. Click Save.
- 4. If the Confirm Save As window opens, click Yes.

The document is saved with scan data, background data peaks, phase data, list of cell candidates and history of the analysis steps.



CHAPTER 11 DO AN AUTOMATIC RIETVELD REFINEMENT (HIGHSCORE PLUS)

11.1 Introduction

This chapter shows you the Rietveld refinement of an artificial mixture of 2 minerals.

You will do an example for quantitative phase analysis. The example has pre-defined refinement steps. An automatic fitting parameter set for phase fits is used for this Rietveld refinement.

11.2 Load data

- 1. If HighScore Plus is not started, start it. Refer to Section 2.2.
- 2. Go to File > Open to open the document "25-75.rd" in this folder C:\Users\user.name \Documents\PANalytical\X'Pert HighScore Plus\Tutorial.
- Go to File > Insert to insert "Example.cry" in this folder C:\Users\user.name\Documents \PANalytical\X'Pert HighScore Plus\Structures. The CIR / CRYSTIN Import Structures opens.
 - **NOTE:** The "Structures" folder contains some crystal structures supplied with HighScore Plus. "Example.cry" contains the missing crystal structure data.
- 4. Select Fluorite and Corundum.

No.	Use	Name	Formula	Space group	Comment	
1		Eskolaite	Cr2 03	R3-cH	Comment not found	
2	V	Fluorite	Ca F2	Fm3-m	Comment not found	
3		Calcite	Ca (C 03)	R3-cH	Comment not found	
4	7	Corundum	Al2 03	R3-cH	Comment not found	
5		Tantalum Sulfide/Se	Ta5(S,Se)2	I4/mmm	Comment not found	

Figure 11.1 Fluorite and Corundum selected

5. Click **OK**. Fluorite and Corundum are inserted into the diffraction document.

11.3 Do an automatic refinement

1. On the **Desktop** toolbar, set the desktop layout to **Rietveld Analysis**.



Figure 11.2 Set the desktop layout

- 2. On the menu bar, go to Analysis > Fitting.
- 3. Set Fitting Mode to Automatic.
- 4. On the menu bar, go to **Analysis > Fitting > Start Fit** and select **<Phase fit> Default Rietveld** to start the refinement.



Figure 11.3 < Phase fit> Default Rietveld selected

The R-Values window opens.



Figure 11.4 Rietveld refinement

After the automatic refinement is completed, the peaks and the calculated profile show in the **Main Graphics** pane.



Figure 11.5 Peaks and the calculated profile after the refinement

- 5. Set the **Additional Graphics** pane to show the difference plot:
 - a. Right click in Additional Graphics pane.
 - b. From the pop-up menu, select **Show Graphics > Difference Plot**.
- 6. Hide the dynamic difference scale to show the small differences better:
 - a. Right-click in the **Additional Graphics** pane.

- b. On the pop-up menu, make sure that **Dynamic Difference Scale** is not selected.
- 7. In the **Main Graphics** pane, zoom in and examine different parts of the profile and the difference plot.
- 8. Examine the agreement indices:
 - a. Go to the **Refinement Control** pane.
 - b. Double-click **Global Variables** to open the **Object Inspector** pane.
 - c. Open Agreement Indices.the values of Goodness of Fit and Weighted R profile.

Quantification	11	Anchor 5	can Data	Pattern	List	Peak L	ist	Selected object: Global Settings
Refinement Co	ntrol		x	Scan List	Stru	ucture Plot	1	🗄 Background
Name	In	Refine	Value	Deviation	Code	Constra	Maxim	Agreement Indices
🕀 🔂 Global Variables							-	R expected 4, 1996
+ (i) Fluorite		1						R profile 16.1367
G Corundum					-			Weighted R profile 20.6787
		Line						D-statistics 0.2006 Weighted D-statistics
								Goodness of Fit 4.9239

Figure 11.6 Agreement Indices

- 9. Examine the phase amounts:
 - Go to **Quantification Pane**. The pie chart shows that the result is 78.3 % Corundum and 21.7 % Fluorite, which is close to the correct weight percentages, 75 % and 25 %.



Figure 11.7 Quantification pane

• Alternatively, go to the **Main Graphics** pane. The phase amounts are usually shown with the phase legend.

11.4 Do a better refinement

- 1. On the menu bar, go to **Analysis > Fitting**.
- 2. Set **Fitting Mode** to **Manual**. In this mode, you can switch on and off the items that can be refined and automatic fitting parameters are no longer used.
- 3. Go to the **Refinement Control** pane.
- 4. Right-click to open the pop-up menu.
- 5. Go to Refine All.
- 6. Select **V's**.

Refine All	•	Scale Factors
Fix All	•	Cells
Show Refine	ed Values / Constraints	W's
Add New Ph	ase (Structure	V's
Add New Ab		U's
Duplicate At	om	Pref. Orientation Parameters
Load Phase	from ICSD by Collection Code	Asymmetry Parameters
		Peak Shape Parameter 1's
Delete Phase	e / Atom	Peak Shape Parameter 2's
Delete All Ph	hases	Peak Shape Parameter 3's
Initialize Glo	bal Variables	
Initialize Pha	ase	
Initialize All F	Phases	
Add All Phas	es to User Reference Database	
Сору То	•	
Paste Phase	es from Dataset	
Duplicate Ph	hase	
Standardize	Phase	-
🚺 Refine Unit (Cell	
Reduce Unit	Cell	
Transform C	iell / Structure	
Create Peak	s from Unit Cell	
Renormalize	Pseudo Formula Masses	
Take as Size	-Strain Standard	
Recalculate	Scales from standard weight percentages	
Calculate Ba	ckground Coefficients from Background	
Expand Roo	t Nodes	
Collapse Roo	ot Nodes	
Autosize Col	lumns	
🔣 Customize R	efinement Control	
눱 Copy List		
칂 Print List		
📕 Save List As	····	

Figure 11.8 V's selected

- 7. Right-click to open the pop-up menu again.
- 8. Go to Refine All.
- 9. Select **U's**.
- 10. Start the refinement again:
 - Go to Analysis > Fitting >Start Fit.
 - Alternatively, on the **Fitting** toolbar, click 膨.

Help									
🗑 🖹 🔌 🖻 🖉	X	0	1	ю	• 6	- 14		Ŧ	
- 🛃 🖕	: *	- 51 -		JD.	\otimes	Man	ual	10223	Ŧ
				5	5				
					Execu	ite Fi	tting		

Figure 11.9 Execute Fitting button

- 11. Right-click in the **Refinement Control** pane to open the pop-up menu.
- 12. Go to Refine All.
- 13. Select Peak Shape Parameter 1's.
- 14. Do step 10 to start the refinement again.
- 15. Change the FWHM to 30:
 - a. Go to the **Refinement Control** pane.

- b. Double-click **Global Variables** to open the **Object Inspector** pane.
- c. Open General Fit Properties.
- d. Change Peak Base Width for Fit from "20" to "30".

Refinement Control			•		Object Inspector	•
Quantification And	hor Scan Data	Pattern List	Peak List		Selected object: Global Settings	
Refinement Control	× So	an List Stru	icture Plot		Background	*
Name	In Refine	Value D	Deviation	c	Agreement Indices	*
🕨 🕀 🔁 Global Variable	s				General Fit Properties	\$
+ 🖻 Fluorite					Job Type	X-rays
±- 🛱 Corundum					Weighting Scheme	Against I
					Solver Time Limit [sec]	300
					Solver Iteration Limit	100
					Solver Nu	2
					Solver Tolerance	0.001
					Max. No of Fit Cycles	20
					Peak Base Width for Fit	30
					Automatic Cell Constraints	V
					Automatic Anisotropic Displacement	v
					Keep ADP's positive definite	v
					Automatic Atom XYZ Constraints	v
					Specimen Displacement [mm]	-0.0746
					Zero Shift [º28]	0

Figure 11.10 Change Peak Base Width for Fit

- e. Press **Enter** to save the change.
- 16. Do step 10 to start the refinement again.
- 17. Examine the agreement indices:
 - a. Go to the **Refinement Control** pane.
 - b. Double-click Global Variables to open the Object Inspector pane.
 - c. Open Agreement Indices.

Ξ,	Agreement Indices	
	R expected	4, 19752
	R profile	8.76649
	Weighted R profile	11.33908
	D-statistics	0.31356
	Weighted D-statistics	0
	Goodness of Fit	2.70138
	Mixture MAC [cm^2/g]	44.98

Figure 11.11 Agreement Indices

- d. Examine the values of Goodness of Fit and Weighted R profile:
 - The value of **Goodness of Fit** is 2.70.
 - The value of Weighted R profile is < 11.34.

You have a better calculated profile with smaller differences to the measurement.

18. Go to Quantification Pane to examine the phase amounts. The phase amounts have small changes: 23.1 % for Fluorite and 76.9 % for Corundum, and they are closer to the given values. The phase quantification is within 2 % of the true values, which is a good result for a phase analysis without standard.

NOTE: A quantitative phase analysis is not always correct with a good fit and low R-values.