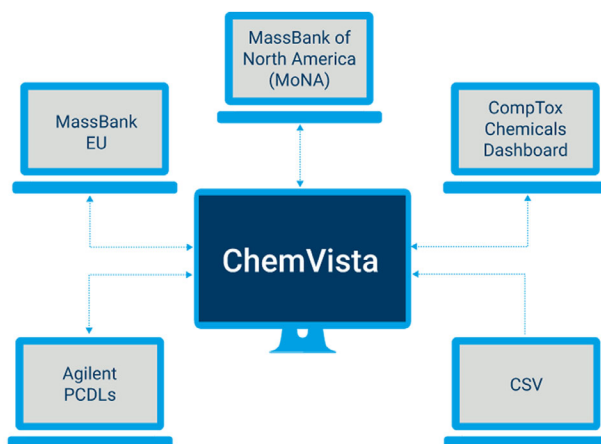


ChemVista Library Manager 1.0

Introduction Workbook



Notices

Document Identification

D0030672
April 2023 Revision A.00

Copyright

© Agilent Technologies, Inc. 2023

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara, CA 95051

Software Revision

This guide is valid for ChemVista Library Manager 1.0 program or higher and compatible ChemVista Library Manager 1.0 programs, until superseded.

Software Manufacturing



Manufactured for Agilent
Technologies
5301 Stevens Creek Blvd
Santa Clara, CA 95051

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

Video Support

This Introduction Workbook has supplemental video support available. Each chapter uses step by step instructions supported by on line videos to view and review the material as needed. Scan the code below or use this link to access the videos (<https://aglt.co/ChemVistaIntro>).



Agilent Community



Agilent Community

To get answers to your questions, join over 10,000 users in the Agilent Community. Review curated support materials organized by platform technology. Ask questions to industry colleagues and collaborators. Get notifications on new videos, documents, tools, and webinars relevant to your work.

<https://community.agilent.com/>

Contents

1	Introduction	7
	About this Training Workbook	8
	How This Training Workbook Works	8
	Notes and Alerts	9
	What This Training Workbook Covers	9
	Requirements	10
2	User Interface and General Navigation	11
	Launching ChemVista Library Manager	12
	User Interface	13
	Pages	15
	Home Page	15
	Search Page	17
	Import Page	18
	Jobs Page	20
	Lists Page	21
	System Page	22
3	Importing Data	23
	Import Files	24
	Introduction	24
	Import data from PCDL	24
	Import data from Mass Bank	26
	Import data from PCDL.CSV	27
4	Lists	31
	Working with Lists	32
	Tag Search	32
	Create a Targeting List from Search Results	33
	Manage Lists from Lists page	35

5	Working with Data	37
	Editing Compounds	38
	Adding Compounds to List	38
	Create a New Substance	39
	Editing Spectra	41
	Add New Spectrum	41
	Verifying Spectra	44
	Update RT	45
6	Exporting Data	47
	Exporting a List	48
7	Appendix	51
	Tips	52
	Using exported PCDLs in MassHunter data analysis	52
	Data Format	52
	3rd Party Resources	53
	MassBank (EU)	53
	MassBank of North America (MoNA)	53
	EPA CompTox Chemicals Dashboard	54
	Common regex searches for Pattern Search	55



1

Introduction

About this Training Workbook	8
How This Training Workbook Works	8
Notes and Alerts	9
What This Training Workbook Covers	9
Requirements	10

About this Training Workbook

This Training Workbook provides instructions on the features of ChemVista Library Manager 1.0.

For additional information on the software and detailed instructions on the features not covered in this workbook, refer to the Online Help.

Use the following exercises to experience how to utilize ChemVista Library Manager to import library data, organize via search and list creation, edit structural and compound metadata, and export to desired formats. Example data is provided with the installation of the software to introduce these steps.

How This Training Workbook Works

This training workbook is your introductory guide for the set-up and execution of ChemVista Library Manager 1.0. This workbook is divided into several chapters, each building upon the last, so we recommend that each chapter is completed in succession. During each chapter, lessons are guided step by step.

Instructions look like this:

Tasks or items needed to complete tasks look like this.

If you are expected to enter any information or if something is important, it is set in italicized type like this:

Type *Blank One* in the field.

If you are expected to press a key on the keyboard or button on the software screen, the key is displayed in bold like this:

press **Enter**.

Cross references appear in blue:

(For example, [Link](#))

Instructions for procedures are presented in a numbered step format:

- 1 Start the software by clicking the icon on the desktop. The main software window appears.

Notes and Alerts

NOTE

The Note text appears here.

CAUTION

The Caution text appears here.

WARNING

The Warning text appears here.

What This Training Workbook Covers



In this learning experience, the goal is to get up and running using the software as quickly as possible. After completing this learning ChemVista Library Manager event, you will have an introductory level of experience in the use of ChemVista Library Manager 1.0.

This learning experience introduces basic concepts in a learning-by-doing, guided manner. Each chapter uses step by step instructions and is supported by on line videos to view and review the material as needed. Scan the code to the right or use this link to access the videos (<https://aglt.co/ChemVistaIntro>). At any time if you have a question or get stuck, to find your local sales and support contact, visit the following page using the link below.

- <https://www.agilent.com/en/contact-us/page>

For technical support, visit the following page:

- <https://www.agilent.com/en/support>

Requirements

To complete the chapters in this training workbook, you need to have installed the software to the most recent release recommended. The software is found on the install media provided in the install kit, along with specially prepared data sets to load onto the software system before the start of this learning exercise.

Please refer to the Online Help or Installation Guide for further information. As always, feel free to contact Agilent Support for additional assistance.



2

User Interface and General Navigation

Launching ChemVista Library Manager 12

User Interface 13

Pages 15

Home Page 15

Search Page 17

Import Page 18

Jobs Page 20

Lists Page 21

System Page 22

Launching ChemVista Library Manager

- 1 After software installation, double-click the ChemVista Library Manager icon or **Click Start > All Programs > Agilent Technologies > ChemVista Library Manager**.



- 2 The software splash screen shows the software initialization.

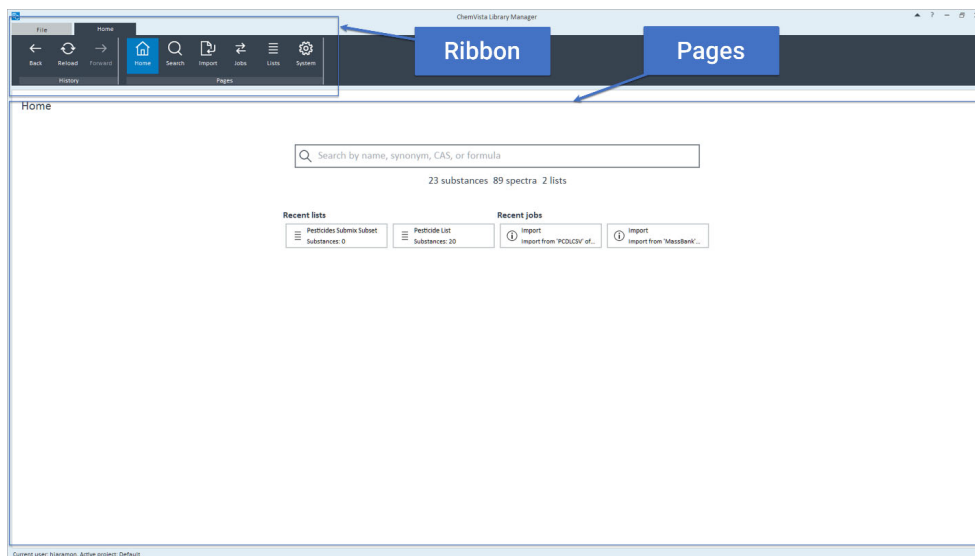


NOTE

The splash screen displays the build number of the software, which is needed when contacting **support**. This information is also available in the **File>About** section of the software.

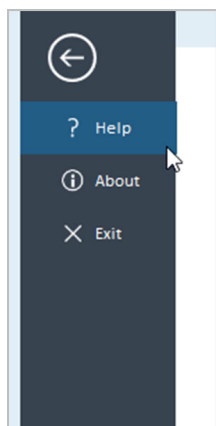
User Interface

Navigation Ribbon



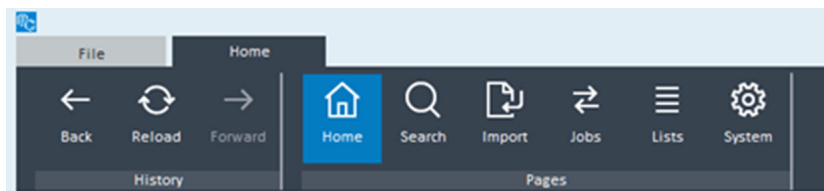
Ribbon Definitions

File Tab - The following commands are available on the File menu.



- Help opens the Online Help file.
- About opens the About Dialog Box
- Exit closes the program.

Home tab - The ribbon has two groups with the following commands and pages:



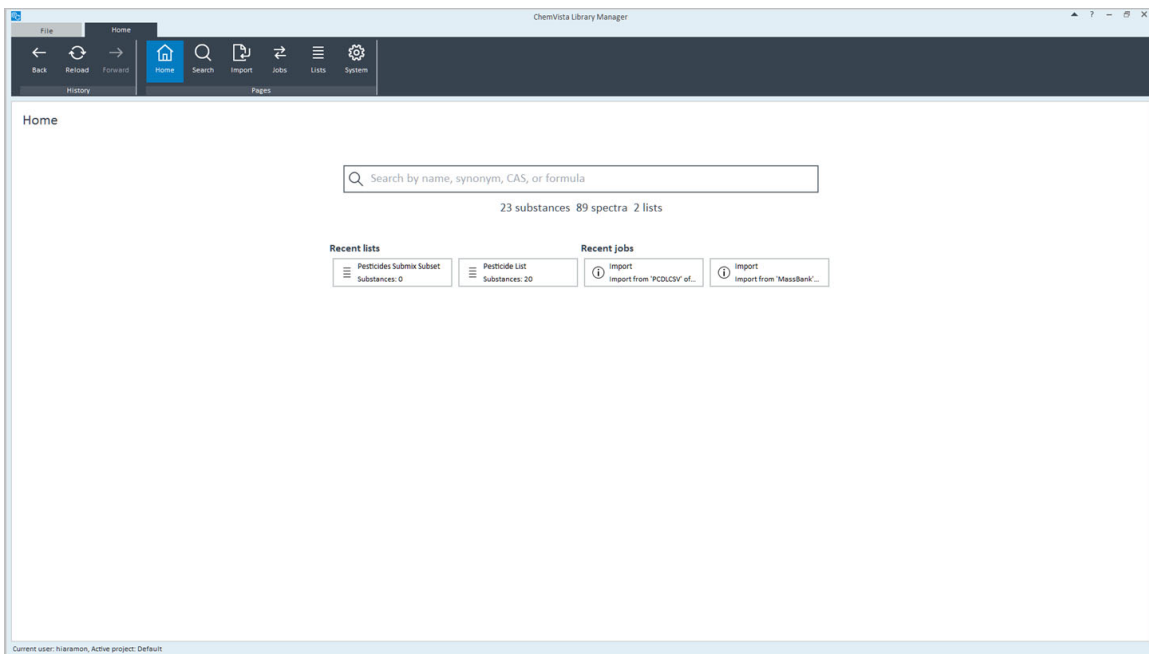
The **Home tab** has two groups with the following commands and pages.

- History
 - **Back** returns to the previous page.
 - **Reload** reloads the current page, updates values, and resets entered text.
 - **Forward** goes to the following page.
- Pages:
 - Home
 - Search
 - Import
 - Jobs
 - Lists
 - System

Pages

The pages and their basic functionality used for these exercises is reviewed in this section. For further definitions or advanced functionality, refer to the Online Help.

Home Page



Home Page Definitions



The Home page is shown when  is clicked on the Ribbon.

Functionality

The Home page shows overall substance and spectral counts and features the ability to search by name, synonym, CAS, or formula. Enter part or all of the name, synonym, CAS, or formula to search for in the field and press Enter on the keyboard to return results.

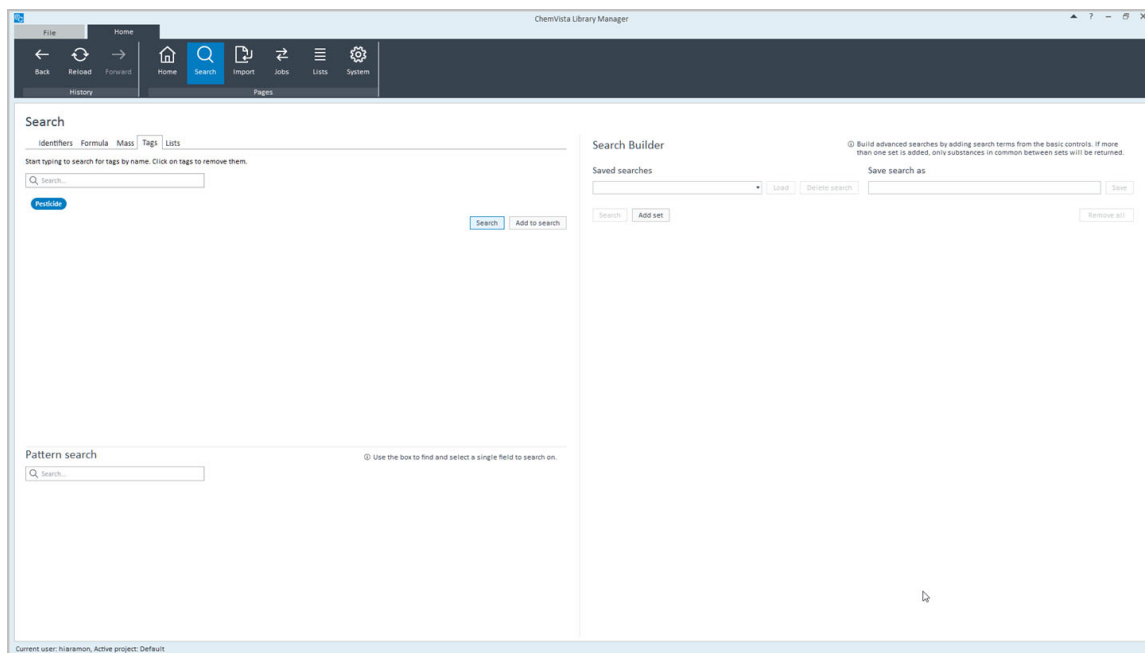
Recent lists

Each box represents a different list, up to 12 recently edited lists are displayed. The box displays the name of the list and how many substances are in the list.

Recent jobs

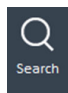
Each box shows jobs that have been started in the last 24 hours.

Search Page



Search Page Definition

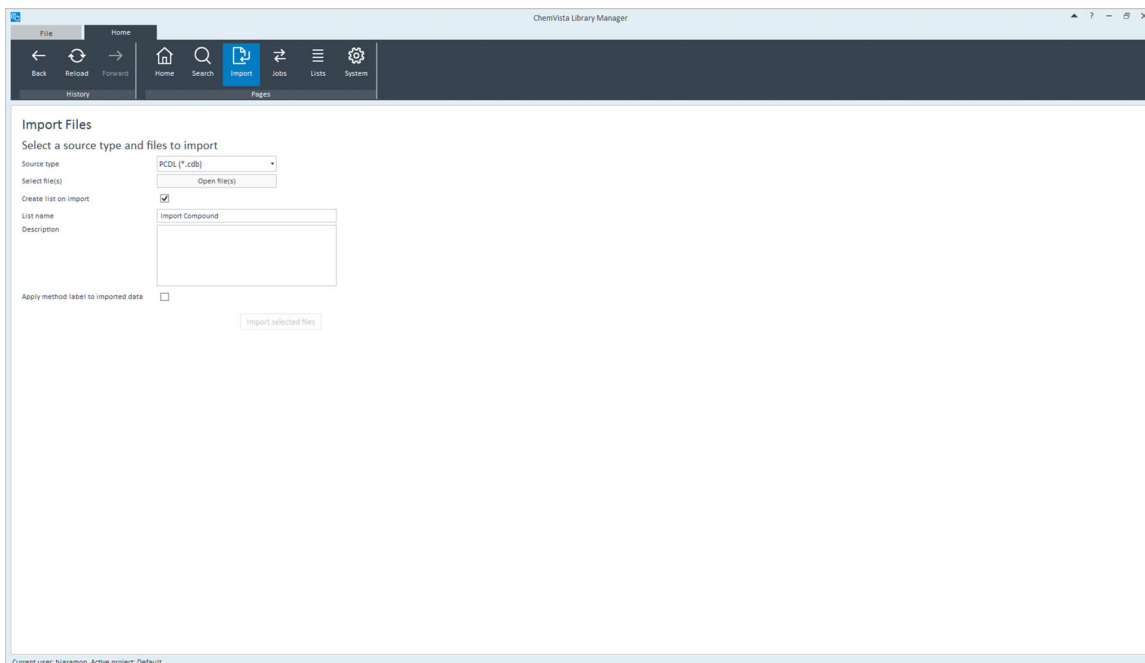


Click  on the Ribbon to show the Search page.

Functionality

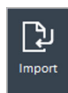
This page is used to search within ChemVista by Identifiers, Formulas, Mass, Tags, or Lists and to view the results of a search.

Import Page



Import Page Definition



Click  on the Ribbon to show the Import page.

Functionality

This page is used to import data into ChemVista from multiple formats.

Source Types

The type of file to be imported. The options include:

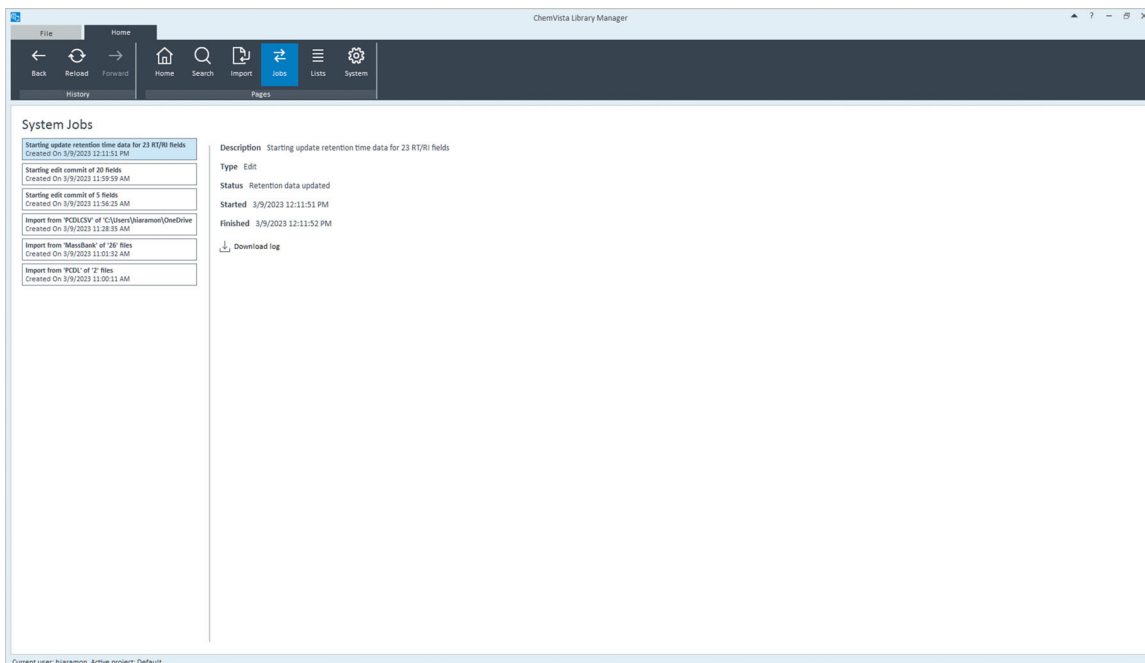
- PCDL (*.cdb)
- PCDL CSV (*.csv)
- MassBank (*.txt)

- SDF (*.sdf)

NOTE

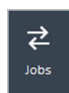
Refer to the Appendix for more information on file types, sources, and examples.

Jobs Page



Jobs Page Definition



Click  on the Ribbon to show the Jobs page.

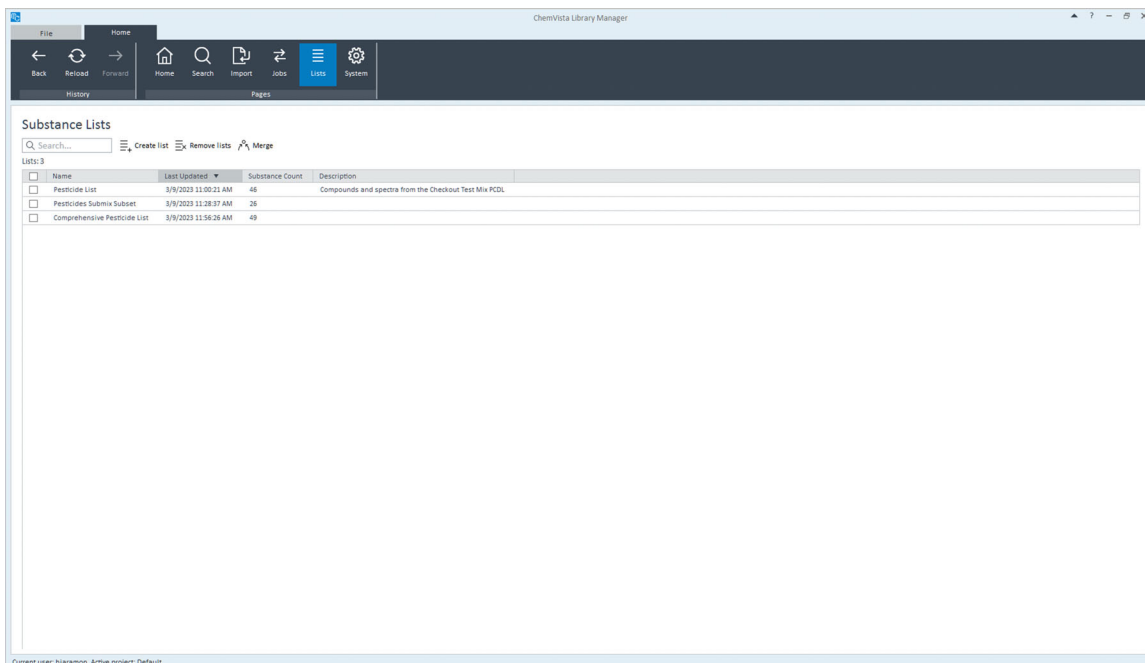
Functionality

This page is used to display information on any data operation (for example, import, add, edit, restore) when created as a job.

NOTE

Any operation that only affects lists (for example creation or editing) does not create a job.

Lists Page



Lists Page Definition

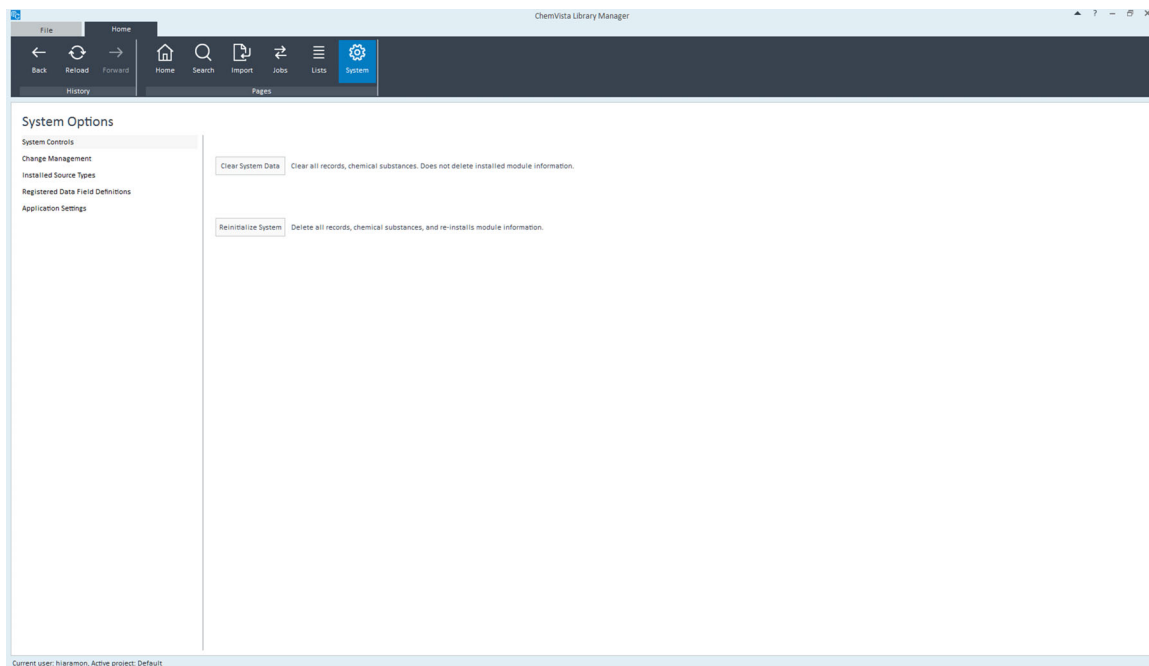


Click  on the Ribbon to show the Lists page.

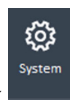
Functionality

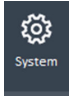
From this page, create, remove, or merge lists, which then display in a table. List names are active links that display the substances in each list when clicked.

System Page



System Page Definition



Click  on the Ribbon to show the System page.

Functionality

From this page clear system data, review registered Data Field definitions, changes, installed source types or review applications settings.

NOTE

These actions should only be done if you understand the effects. Please consult the Online Help for further information.

3 Importing Data

Import Files 24

Introduction 24

Import data from PCDL 24

Import data from Mass Bank 26

Import data from PCDL.CSV 27

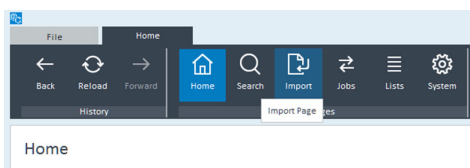
Import Files

Introduction

The following exercises are designed to support the execution of the ChemVista Library Manager features along with video support. The videos are intended to provide visual support for working with the software. We recommend reviewing the video demonstration first, then attempting the processes, using each exercise as a guide. If at any time you have questions or run into an event that is not in alignment with this workbook, please reach out to Agilent Support or your Agilent Consultant.

Import data from PCDL

- 1 Click **Import** on the Ribbon to navigate to the Import Page.



- 2 Select **PCDL(*.cdb)** from the Source type drop-down.

Import Files

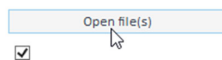
Select a source type and files to import

Source type	PCDL (*.cdb)
Select file(s)	PCDL (*.cdb)
Create list on import	PCDL CSV (*.csv)
List name	MassBank (*.txt)
Description	SDF (*.sdf)

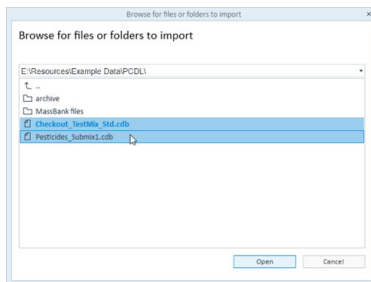
Importing Data

Import data from PCDL

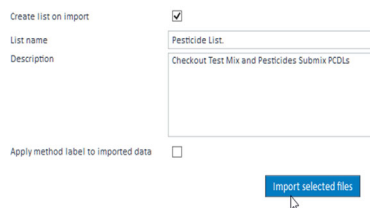
- 3 Click **Open file(s)**.



- 4 Navigate to the folder “\Resources\Example Data\PCDL” and select **Checkout_TestMix_Std.cdb** and **Pesticides_Submix1.cdb**.



- 5 Click **Open**.
- 6 Select **Create List Upon Import**.
- 7 Enter the name *Pesticide List* into the List Name field.
- 8 Enter *Checkout Test Mix and Pesticides Submix PCDLs* into Description.
- 9 Select **Import selected files** to launch the import.



- 10 Click **OK**.

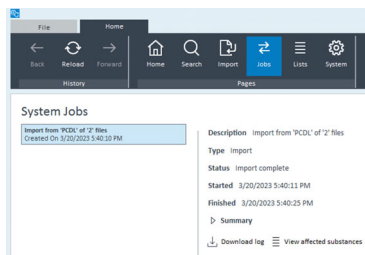
NOTE

The PCDL is not saved within the software, the data is imported from the PCDL and stored in the software's internal format.

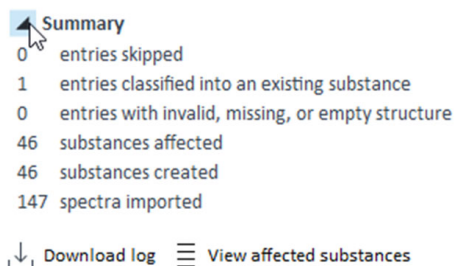
Importing Data

Import data from Mass Bank

- 11 Review the Status of the import from the System Jobs page to confirm it reads Import complete.



- 12 Expand the Summary to review the number of substances created and spectra imported.



Import data from Mass Bank

- 1 Click **Import** on the ribbon to navigate to the Import Page.

NOTE

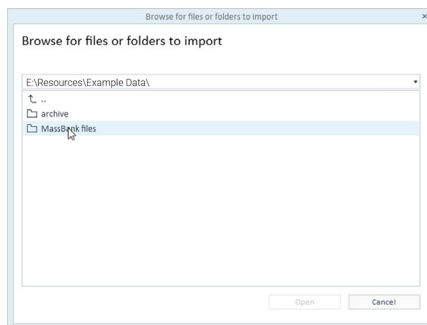
The MassBank files provided with the installation software are only for instruction purposes. It is advised to obtain updated files from MassBank before performing analysis beyond performing these exercises.

- 2 Select **MassBank (*.txt)** from the Source type drop-down.
- 3 Click **Open file(s)**.

Importing Data

Import data from PCDL.CSV

- 4 Navigate to the folder "\\Your Directory\\Example Data\\" and select the **MassBank files** folder.



- 5 Click **Open**.
- 6 Clear the **Create List Upon Import** check box.

Import Files

Select a source type and files to import

Source type	MassBank (*.txt)	
Select file(s)	Open file(s)	26 files selected
Create list on import	<input type="checkbox"/>	
Apply method label to imported data	<input type="checkbox"/>	

- 7 Click **Import Select Files** to launch the import.
- 8 Click **OK**.
- 9 Expand the Summary to review the number of substances created and spectra imported.

NOTE

The MassBank files are not saved within the software, the data is imported from the files and stored in the software's internal format.

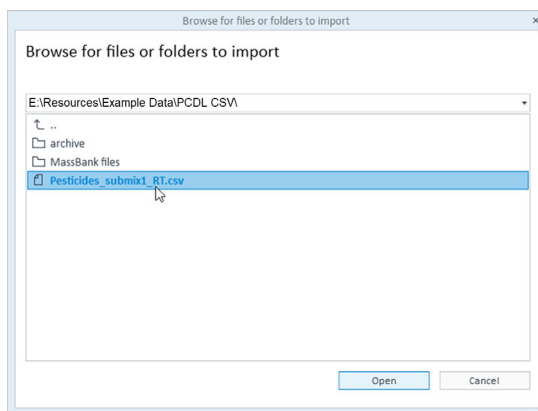
Import data from PCDL.CSV

- 1 Click **Import** in the Ribbon to navigate to the Import Page.
- 2 Select **PCDL CSV (*.csv)** from the Source type drop-down
- 3 Click **Open file(s)**.

Importing Data

Import data from PCDL.CSV

- 4 Navigate to the folder "\\Your Directory\\Example Data\\" and select **Pesticides_submix1_RT.csv**.



- 5 Click **Open**.
- 6 Select the **Create List Upon Import** check box.
- 7 Enter the name *Pesticide Submix Subset* into the List Name field.

List name

- 8 Select the **Apply method label to imported data** check box.

Apply method label to imported data ☒

Provide a label for the imported methods or select from an existing one at right

- 9 Enter *Agilent Eclipse Plus C18, 2.1 mmx150 mm, 01/2021* into the label field.
- 10 Click **Import selected files** to launch the import.

Provide a label for the imported methods or select from an existing one at right

NOTE

Creating a method label is strongly recommended for optimal RT/RI data organization, especially from PCDL and CSV imports. The method label is stored alongside RT/RI data and enhances viewing and downstream use.

- 11 Click **OK**.

Importing Data

Import data from PCDL.CSV

- 12 Expand the **Summary** to review the number of substances created and spectra imported.
- 13 Click **Lists** on the Ribbon to view the Substance List pane.
- 14 View the Substances Lists, noting the two lists created via import display.

Substance Lists

Search...

Create list

Remove lists

Merge

Lists: 2

<input type="checkbox"/>	Name	Last Updated ▼	Substance Count	Description
<input type="checkbox"/>	Pesticide Submix Subset	3/20/2023 5:56:53 PM	26	
<input type="checkbox"/>	Pesticide List	3/20/2023 5:40:24 PM	46	Compounds and spectra from the Checkout Test Mix and Pesticides Submix PCDLs

This page intentionally left blank.

4

Lists

Working with Lists **32**

Tag Search **32**

Create a Targeting List from Search Results **33**

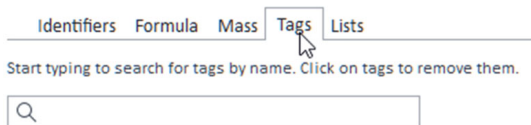
Manage Lists from Lists page **35**

Working with Lists

Tag Search

- 1 Click **Search** on the Ribbon to navigate to the Search Page.
- 2 Click the **Tags** tab.

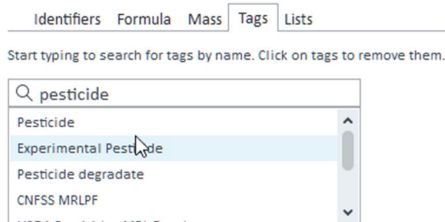
Search



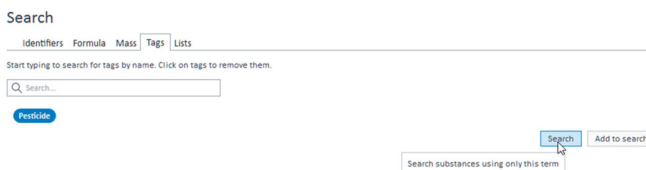
NOTE

Tags come from rich data that's included as part of the Agilent PCDL files. They're parsed from the compound Description and Notes fields and presented as a separate, searchable tag that travels along with the compound information.

- 3 Enter *pesticide* into the search field, then select it from the auto-generated list.



- 4 Click **Search** to load the subset in the Search Results window.



NOTE

Tags are removed by clicking on the text to clear from the search list.

Create a Targeting List from Search Results

- 1 Use the **Select All** check box to select all compounds in the Search results.

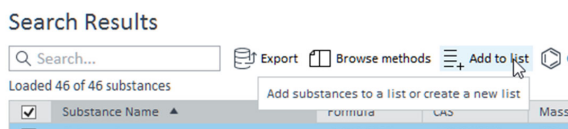
Search Results

Q Search...

Loaded 46 of 46 substances

<input checked="" type="checkbox"/>	Substance Name	Formula	CAS	Mass	InChIKey	Agilent ID	Spectra Count
<input checked="" type="checkbox"/>	Spirooxamine	C18H35NO2	118134-30-8	297.26678	PUYXTUJWRLC	3190	3
<input checked="" type="checkbox"/>	Methamidophos (Metamidophos)	C2H8NO2PS	10265-92-6	141.00134	NNKVPKMPCC	3689	3
<input checked="" type="checkbox"/>	Bentazone	C10H12N2O3S	25057-89-0	240.05686	ZOMSMJXLGFI	132	7
<input checked="" type="checkbox"/>	Proquinazid	C14H17IN2O2	189278-52-4	372.03347	FUVBXVXXML	20022	3
<input checked="" type="checkbox"/>	Dimoxystrobin	C19H22N2O3	149961-52-4	326.16304	WXUZAHNPNV	7025	3
<input checked="" type="checkbox"/>	Atrazine	CBH14CIN5	1912-34-9	215.09377	MXWVTDORC	6328	3
<input checked="" type="checkbox"/>	Acifluorfen	C14H7ClF3NO	50594-66-6	360.99648	NUPNQVOELL	1811	3
<input checked="" type="checkbox"/>	Tralkoxydim	C20H27NO3	87820-88-0	329.19909	DOQPEVARZIQ	3889	3
<input checked="" type="checkbox"/>	Lentacil	C13H18N2O2	2164-08-1	234.13683	ZTMKADLOSYP	1272	9
<input checked="" type="checkbox"/>	Thiabendazole	C10H7N3S	148-79-8	201.05807	WICNZQLZVW	410	3
<input checked="" type="checkbox"/>	2,4,5-TP / Silvex (Fenoprop)	C9H7Cl3O3	93-72-1	267.94608	ZLSWBLPERHF	6668	3
<input checked="" type="checkbox"/>	Spirodiclofen	C21H24Cl2O4	148477-71-8	410.10516	DTDSAWVUFF	7449	3

- 2 Click **Add to list**.



- 3 In the Add to new list section of the dialog box, enter the List name *Comprehensive Pesticide List*.

Create a Targeting List from Search Results

The screenshot shows a dialog box titled "Add to List" with a close button (X) in the top right corner. The dialog is divided into two main sections: "Append to existing list" and "Add to new list".

Append to existing list

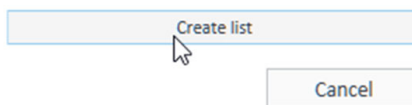
This section contains a label "List:" followed by a dropdown menu currently showing "Pesticide List". Below the dropdown is an "Append" button.

Add to new list

This section contains a label "Name" followed by a text input field containing "Comprehensive Pesticide List.". Below the name field is a label "Description" followed by a large, empty text area with a vertical scrollbar on the right side. At the bottom of this section is a "Create list" button. A mouse cursor is pointing at the "Create list" button.

At the bottom right of the dialog box is a "Cancel" button.

- 4 Click **Create list**.



- 5 Click **OK**.

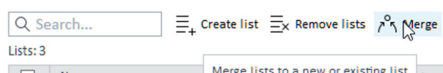
Manage Lists from Lists page

- 1 Click Lists on the Ribbon to navigate to the Lists page.
- 2 Select the Pesticides Submix Subset and the Comprehensive Pesticide List's check box.

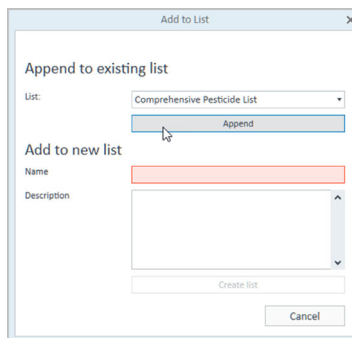
Lists: 3				
<input type="checkbox"/>	Name	Last Updated ▼	Substance Count	Description
<input type="checkbox"/>	Pesticide List	3/9/2023 11:00:21 AM	46	Compounds and spectra from the Checkout Test Mix PCDL
<input checked="" type="checkbox"/>	Pesticides Submix Subset	3/9/2023 11:28:37 AM	26	
<input checked="" type="checkbox"/>	Comprehensive Pesticide List	3/9/2023 11:45:18 AM	46	

- 3 Click **Merge**.

Substance Lists



- 4 In the Add to List dialog box, select the **List** drop down and choose **Comprehensive Pesticide List**.
- 5 Click **Append**.



- 6 Click **OK**.
- 7 The Substance Lists page refreshes and the Comprehensive Pesticidelist displays 47 in Substance count.

Lists: 3

<input type="checkbox"/>	Name	Last Updated ▼	Substance Count	Description
<input type="checkbox"/>	Pesticide Submix Subset	3/20/2023 5:56:53 PM	26	
<input type="checkbox"/>	Pesticide List	3/20/2023 5:40:24 PM	46	Compounds and spectra from the Checkout Test Mix and Pesticides Submix PCDLs
<input type="checkbox"/>	Comprehensive Pesticide List	3/21/2023 12:38:05 PM	47	



5

Working with Data

Editing Compounds 38

Adding Compounds to List 38

Create a New Substance 39

Editing Spectra 41

Add New Spectrum 41

Verifying Spectra 44

Update RT 45

Editing Compounds

Adding Compounds to List

- 1 Click **Search** on the Ribbon to navigate to the Search page.
- 2 Select the **Identifiers** tab.

Search

Identifiers Formula Mass Tags Lists

Select an identifier field: Name

① Separate multiple values with a line break.

- 3 On the Identifiers tab, select the identifier field CAS.

Search

Identifiers Formula Mass Tags Lists

Select an identifier field: CAS

Name
CAS
InChIKey
SMILES

- 4 Enter 2886-65-9 in the search field.

Search

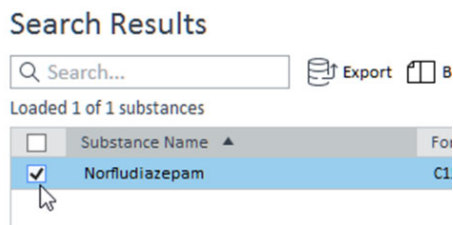
Identifiers Formula Mass Tags Lists

Select an identifier field: CAS

2886-65-9

- 5 Click **Search**.

- 6 Select **Norfluradiazepam**.



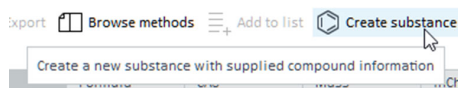
- 7 Click **Add to list**.
 8 Select **Comprehensive Pesticide List**.
 9 Click **Append**.
 10 Click **OK**.

Create a New Substance

- 1 Click **Lists** in the Ribbon to navigate.
 2 Click **Comprehensive Pesticide Screening** list name to view.

<input type="checkbox"/>	Name	Last Updated ▼	Substance Count	Description
<input type="checkbox"/>	Pesticide List	3/9/2023 11:00:21 AM	46	Compounds and spectra from the Checkout Test Mix PCDL
<input type="checkbox"/>	Pesticides Submix Subset	3/9/2023 11:28:37 AM	26	
<input checked="" type="checkbox"/>	Comprehensive Pesticide List	3/9/2023 11:53:09 AM	48	

- 3 Click **Create substance**.



- 4 Enter the following information:
- Name: *Difenoxuron*
 - CAS: 14214-32-5
 - Formula: *C16H18N2O3*
 - SMILES: *COC1C=CC(=CC=1)OC1C=CC(=CC=1)NC(=O)N(C)C*
 - Mass: 286.13174

Dialog box titled "Edit Substance Details" showing fields for substance information:

Field	Value
Name	Difenoxuron
CAS	14214-92-5
Formula	C16H18N2O3
InChI	
InChIKey	
SMILES	COC1C=CC(=CC=C1)OC1C=CC(=CC=C1)N(C)=O[N](C)C
Mass	286.13174
LogP	
Anion	<input type="checkbox"/>
Cation	<input type="checkbox"/>
IUPAC Name	
ChemSpider	
PubChem	
KEGG	
DTXSID	

Buttons: Add custom field, Add to current list, Commit, Cancel

NOTE

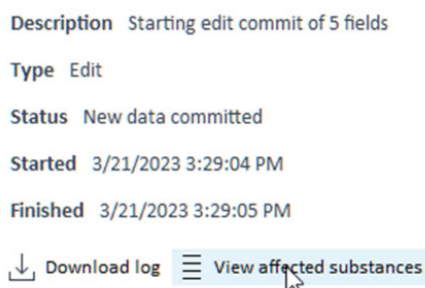
Adding any single structural identifier results in the creation of any missing identifiers, including InChI String, InChIKey, MOL text, and a structural image.

5 Click **Commit**. The Systems Jobs pane displays.

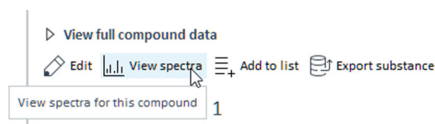
Editing Spectra

Add New Spectrum

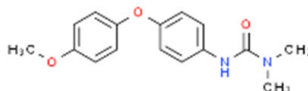
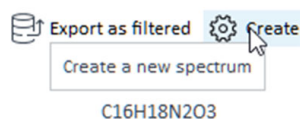
- 1 From Systems Jobs page, click View Affected Substances.



- 2 Click View spectra.



- 3 Click Create.



4 Add the following information into **m/z** and **Intensity**:

m/z	Intensity
46.065126	19.316
72.04439	32.018
123.044056	18.668
124.051881	1.139
287.139019	100

5 Edit the Spectral Details with the following information:

- Separation technique: **LC**
- Mass Analyzer: **QTOF**
- Ionization Technique: **ESI**
- Polarity: **POSITIVE**
- Precursor Type: **[M+H]⁺**
- CE: *10*
- Precursor ion: *287.1390189076*
- MS Level: 2

Working with Data

Add New Spectrum

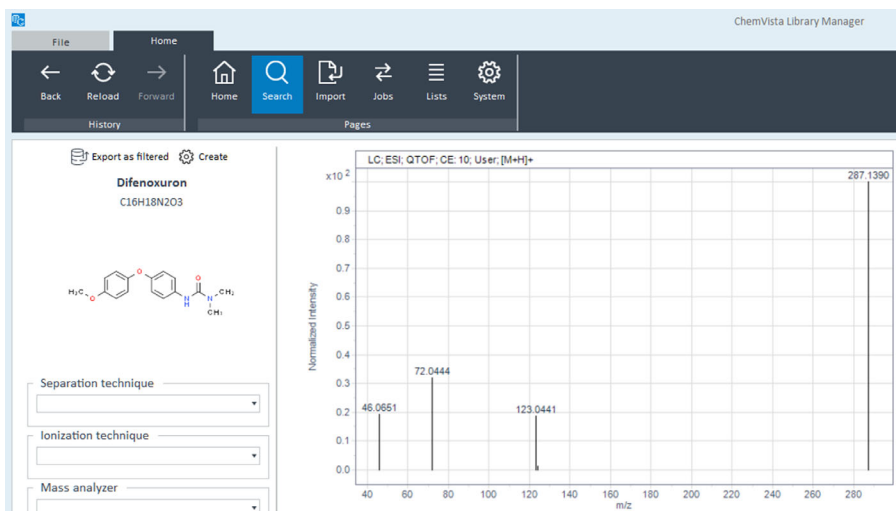
Edit Spectral Details

m/z	Intensity
46.065126	19.316
72.04439	32.018
123.044056	18.668
124.051881	1.139
287.139019	100

Separation technique: LC
 Mass analyzer: QTOF
 Ionization technique: ESI
 Polarity: POSITIVE
 Precursor type: [M+H]⁺
 CE: 10
 Precursor ion: 287.1390189076
 MS Level: 2
 Fragmentor:

Commit Cancel

- Click **Commit**.
- From Systems Jobs, click **View Affected Substances**.
- Click **View Spectra** to verify that the data was added.
- Review the data as needed.



Verifying Spectra

- 1 Click **Lists** on the Ribbon to navigate to the Lists page.
- 2 Click **Comprehensive Pesticide List** to view the list of compounds.
- 3 Sort the list alphabetically by clicking the **Substance Name** column header.

<input type="checkbox"/>	Substance Name ▲	For
<input type="checkbox"/>	2,4,5-T / 2,4,5-Trichlorophenoxyacetic acid	C8
<input type="checkbox"/>	2,4,5-TP / Silvex (Fenoprop)	C9
<input type="checkbox"/>	Acephate	C4
<input type="checkbox"/>	Acifluorfen	C1
<input type="checkbox"/>	Aminocarb	C1
<input type="checkbox"/>	Atrazine	C8
<input type="checkbox"/>	Azoxonazole	C1

- 4 Highlight **Bentazone**.

<input type="checkbox"/>	Bentazone	C10H12N2O3S	25057-89-0	240.05686	ZOMSMJLGF1	132	7
--------------------------	-----------	-------------	------------	-----------	------------	-----	---

- 5 Click **View Spectra**.
- 6 Click **Polarity**.

Polarity

NEGATIVE
POSITIVE

- 7 Select **Negative**.

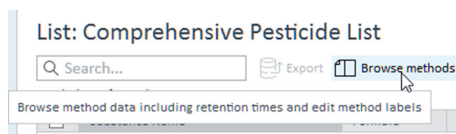
Polarity

NEGATIVE

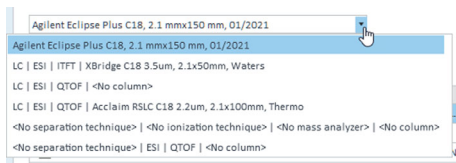
- 8 View Negative mode spectra and confirm 3 spectra are present.

Update RT

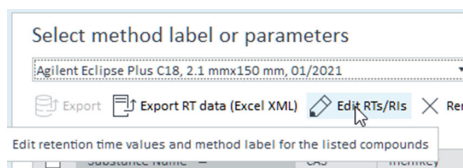
- 1 From the ribbon, click **Back** to return to the prior page.
- 2 Click **Browse Methods**.



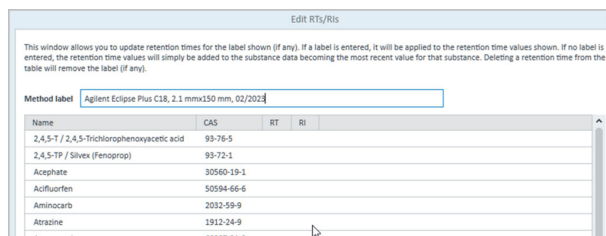
- 3 Click **Select method label or parameters** to view the methods available. Click anywhere to close.



- 4 Click **Edit RT /RIs**.



- 5 Type in *Agilent Eclipse Plus C18, 2.1 mmx150 mm, 02/2023* in the Method Label field.



- 6 Click the **RT** column header to sort the list.

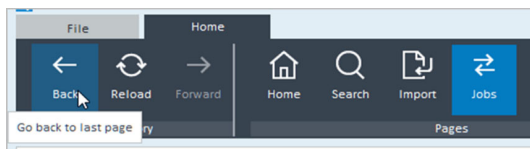
Name	CAS	RT	RI
Fluopicolid	239110-15-7	5.115	

- 7 Add additional Retention times in the **RT** column as needed.

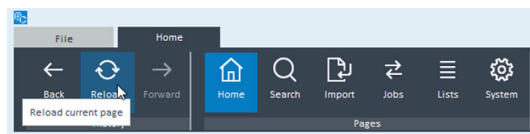
Acephate	30560-19-1	3.189
Thiabendazole	148-79-8	3.209
Lenacil	2164-08-1	

- 8 Click **Commit**.

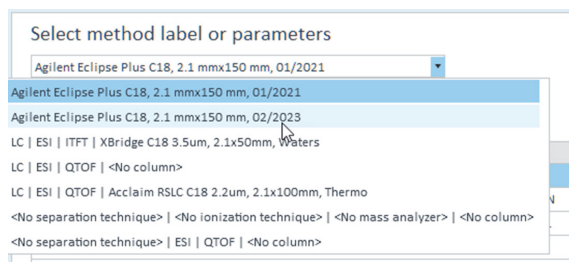
- 9 To confirm the data is loaded in the software, click **Back** on the Ribbon.



- 10 Click **Reload** on the Ribbon of the Methods page.



- 11 Select method label or parameters to confirm the method was applied to the list.





6 Exporting Data

Exporting a List **48**

- 1 Select the **Select All** check box.

2 Click **Export** to navigate to the Export Page and display options.

3 Confirm that PCDL (*.cbd) is selected for Export As.

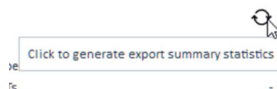
4 Click **Ionization technique** and select **ESI**.

5 Click **Polarity** and select **Positive**.

6 Click **Method label** and select **Agilent Eclipse Plus, C18, 2.1 mmx150 mm, 02/2023**.

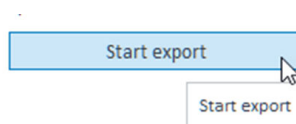
ChemVista Library Manager 1.0 Introduction Workbook

- 7 Click **Summary Load Statistics** to review summary statistics of the data included in the export.

**CAUTION**

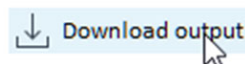
Loading of Summary Statistics may take several minutes if the exported data set is large.

- 8 Click **Start Export**.



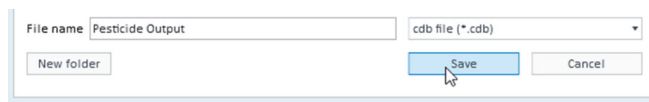
- 9 Click **OK**.

- 10 Click **Download output**.



- 11 In the Save file dialog box, browse to the desired location to save the file.

- 12 Enter a name for the output and click **Save**.



This page intentionally left blank.

Tips **52**

Using exported PCDLs in MassHunter data analysis **52**

Data Format **52**

3rd Party Resources **53**

MassBank (EU) **53**

MassBank of North America (MoNA) **53**

EPA CompTox Chemicals Dashboard **54**

Common regex searches for Pattern Search **55**

Tips

Using exported PCDLs in MassHunter data analysis

Exporting PCDLs from Agilent ChemVista enables screening and identification workflows in downstream MassHunter data analysis software. For guidance on using PCDLs in MassHunter Quantitative Analysis, refer to the MassHunter Quant LC/Q-TOF Screener (M6005-10006).

Spectra and RTs can be extracted in MH Qual and sent to a desired PCDL. Once the desired spectra and RTs are in the PCDL, import the PCDL into ChemVista. The import process will use the structural identifiers and meta data to merge data together and eliminate duplicate compounds and will store the new spectra and RTs accordingly. On import, create a method label to keep RTs organized.

Data Format

Format	Support File Type	Example Data/Access
Agilent PCDL	.cdb	E:\Resources\Example Data
CSV	.csv	E:\Resources\Example Data E:\Resources\Templates

3rd Party Resources

Format	Support File Type	Example Data/Access
MassBank EU	.txt	E:\Resources\Example Data https://massbank.eu/MassBank/ https://github.com/MassBank/MassBank-data
MassBank of North America (MoNA)	.sdf	https://massbank.us/ https://mona.fiehnlab.ucdavis.edu/downloads
EPA CompTox Chemicals Dashboard	.sdf	https://comptox.epa.gov/dashboard/ https://comptox.epa.gov/dashboard/chemical-lists

MassBank (EU)

Spectra from MassBank is viewed on the MassBank web application (<https://massbank.eu/MassBank/>) and downloaded via the MassBank data repository hosted on GitHub (<https://github.com/MassBank/MassBank-data>) as .txt files. The data repository is organized in sub-folders according to the contributor or contributing group. Each .txt file contains a single spectral record and there may be many .txt files mapping to the same chemical substance or compound. Downloaded .txt files are imported directly into ChemVista.

MassBank of North America (MoNA)

Spectra from MoNA is viewed and downloaded from the MoNA web application. Spectra are downloaded from MoNA (<https://mona.fiehnlab.ucdavis.edu/downloads>) in three different formats, but

the format supported by ChemVista is the SDF. The SDF download is organized by spectral record, but all entries in the download are contained in a single downloaded file. Downloaded SDFs are imported directly into ChemVista.

Many subset downloads are available from the MoNA downloads page, the following are recommended downloads to supplement Agilent MassHunter data analysis applications:

- All LC-MS/MS QTOF download
- All LC-MS/MS Agilent QTOF download
- All LC-MS/MS Orbitrap download

EPA CompTox Chemicals Dashboard

The EPA CompTox Chemicals Dashboard is a chemical data resource. Data is downloaded from multiple places in four different formats, but the format supported by ChemVista is the SDF (in either v2000 or v3000 MOL format). Data can be downloaded from pre-curated lists of interest (<https://comptox.epa.gov/dashboard/chemical-lists>) as well as from batch searches (<https://comptox.epa.gov/dashboard/batch-search>) conducted in the Dashboard web application. No spectra are available for download from the Dashboard.

Suggested export filters for MassHunter data analysis

When using third party data in MassHunter workflows, the possibility exists that certain spectra are not compatible with MassHunter data analysis algorithms. For maximum compatibility, the following export filters are recommended:

- MS level = 1 or 2
- Polarity = NEGATIVE and/or POSITIVE
- For optimal library matching, restricting the instrument type to 'QTOF' is recommended.

NOTE

MS Level 2 is for LC-MS/MS and GC-MS/MS; MS Level 1 is for GC-MS.

NOTE

Forcing a polarity to be selected removes spectra that have errant polarity values from third party sources that would cause issues in data analysis.

NOTE

Collision Energy (CE) values that are not supported in the PCDL format (e.g., 'Ramp 21.1-31.6 eV') are written out by ChemVista with the value '9999' in the CE field in the exported PCDL.

NOTE

For maximum compatibility with the PCDL format, the presence of a formula and mass are required for chemical substances to be included in the export. This export option is automatically enforced for PCDL exports. For consistency and compatibility outside of MassHunter workflows, it is recommended to select the same filter when exporting to SDF or MassBank.

Common regex searches for Pattern Search

The Pattern Search option on the Search page in ChemVista supports regular expression (regex) inputs for the selected data field. Some common regex queries that might be helpful are provided below:

Regular Expression	Search Conducted
(?i)	When used prior to a search string, makes the search case insensitive ("(?i)pesticide")
.	Matches anything
\d	Matches any digit 0-9
[a-d]	Matches anything containing any of the letters a, b, c, or d
[0-9]	Matches anything containing any of the values 0, 1, 2, 3, 4, 5, 6, 7, 8, or 9
^([0-9] [1-9][0-9] [1-9][0-9][0-9])\$	Matches any value between 0 and 999

www.agilent.com

© Agilent Technologies, Inc. 2023
D0030672
Revision A.00

April 2023

