

# Automated Axiom™ Analysis Software Library File Installation

## Expectations

- User of this procedure has familiarity with Linux command lines and its terminologies.
- The array package is in a zip format.
- Have access to ssh and the system account information to remotely log into the server where Core and Automated Axiom is running.
- Zip package(s) are located at a network location from which the system account on the server has access to the zip files to be installed.

## Installing

1. Obtain the location of library zip package to be installed and make sure it is accessible to the Client's server and system account.
2. Obtain the Client's log in information.
3. "ssh" into the server using system account information.
4. Change directory to: **cd /srv/core/roots/libs/automated-axiom**
5. Depending on what library files need to be installed, follow steps 6 and/or 7.
6. Install the **ax\_threshold** file. Example: **cp ../../install\_v1/libraries/Human.v5.ax\_thresholds ./**
7. Install library package zip files. Repeat step **7c** for each zip package.
  - a. If the library type is a newer revision of one already installed in the directory, the older one must be removed from the directory. If not, analysis issues may occur if the array type is new to the directory.
  - b. Note that array folders sometimes are of the package name which includes a "rX". X denotes the revision number in the directory name, while others may not. Regardless, there may only be one array type in the "automated-axiom" folder.
  - c. Unzip a zip package.
    - 1) Type **unzip [path to zip file] -d ./[zip file name]**
      - i. Where "[path to zip file]" is, enter the path where the zip package is accessible to the system.

- ii. **-d** is a flag to create a folder with name of array library package.
  - iii. Create and name new library destination folders using the same name has the library package zip file minus ".zip". For example, the folder name must be **Axiom\_PMDA.r7** not **Axiom\_PMDA.r7.zip**
8. From the Client's browser (that accesses the Automated Axiom application), do the following:
    - a. Navigate to the **Analysis Setup** page. If already on this page, refresh it.
    - b. Click the **Array Types** drop-down, to verify the installed library packages are listed. **Note:** Library package names may differ from the name displayed in the drop-down.
    - c. If the expected array package(s) are listed in the drop-down, then the installation was a success. If not, verify the missing package folder is in the library directory and that the files reside inside the folder.
  9. Close the ssh terminal, then type **exit**.



Affymetrix, Inc.  
3450 Central Expressway  
Santa Clara, CA 95051

The information in this guide is subject to change without notice.

**DISCLAIMER**

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Revision history:** MAN1000085

| Revision | Date       | Description      |
|----------|------------|------------------|
| A.0      | April 2024 | Initial release. |

**Important Software Licensing Information**

Your installation and/or use of this Automated Axiom Analysis software is subject to the terms and conditions contained in the End User License Agreement (EULA) which is incorporated within the Automated Axiom Analysis software, and you will be bound by the EULA terms and conditions if you install and/or use the software.

**Legal entity**

Affymetrix, Inc. | Santa Clara, CA 95051 USA | Toll Free in USA 1 800 955 6288

**TRADEMARKS**

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

©2024 Thermo Fisher Scientific Inc. All rights reserved.