



User Manual

Sample Attribute Editor

**For research use only.
Not for use in diagnostic procedures.**

Trademarks

©Affymetrix, Inc. All rights reserved. Affymetrix® Axiom®, Command Console®, CytoScan®, DMET™, GeneAtlas®, GeneChip®, GeneChip-compatible™, GeneTitan®, Genotyping Console™, myDesign™, NetAffx®, OncoScan™, Powered by Affymetrix™, Procarta®, and QuantiGene® are trademarks or registered trademarks of Affymetrix, Inc. All other trademarks are the property of their respective owners.

Limited License

Subject to the Affymetrix terms and conditions that govern your use of Affymetrix products, Affymetrix grants you a non-exclusive, non-transferable, non-sublicensable license to use this Affymetrix product only in accordance with the manual and written instructions provided by Affymetrix. You understand and agree that, except as expressly set forth in the Affymetrix terms and conditions, no right or license to any patent or other intellectual property owned or licensable by Affymetrix is conveyed or implied by this Affymetrix product. In particular, no right or license is conveyed or implied to use this Affymetrix product in combination with a product not provided, licensed, or specifically recommended by Affymetrix for such use.

Patents

Cartridge Array Software: **Products may be protected by one or more of the following patents:** 5,733,729; 5,795,716; 5,974,164; 6,066,454; 6,090,555; 6,185,561; 6,188,783; 6,223,127; 6,228,593; 6,229,911; 6,242,180; 6,308,170; 6,361,937; 6,420,108; 6,484,183; 6,505,125; 6,510,391; 6,532,462; 6,546,340; 6,687,692; 6,607,887; 7,062,092 ; 7,451,047; 7,634,363; 7,674,587 and other U.S. or foreign patents.

HT Array Plate/Strip: **Products may be covered by one or more of the following patents:** U.S. Patent Nos. 6,090,555; 6,611,767; 6,687,692; 6,829,376; 7,130,458; 7,451,047; 7,634,363; 7,674,587 and other U.S. or foreign patents.

Copyright

© 2012 Affymetrix Inc. All rights reserved.

Contents

Chapter1	Sample Attribute Editor	1
	Starting the Sample Attribute Editor	1
	Adding or Removing Sample Files From the Sample Attribute Editor	2
	Adding, Editing, or Removing Attributes	2

Sample Attribute Editor

The Sample Attribute Editor enables you to edit or add attributes to an existing sample file (ARR). Topics in this user manual include:

Starting the Sample Attribute Editor

Adding or Removing Sample Files From the Sample Attribute Editor on page 2

Adding, Editing, or Removing Attributes on page 2

Starting the Sample Attribute Editor

1. Select **Affymetrix** → **Sample Attribute Editor** from the Start menu.
2. Select sample files (ARR) in the dialog box that appears and click **Open**.

The Sample Attribute Editor opens and contains the selected sample files (Figure 1).

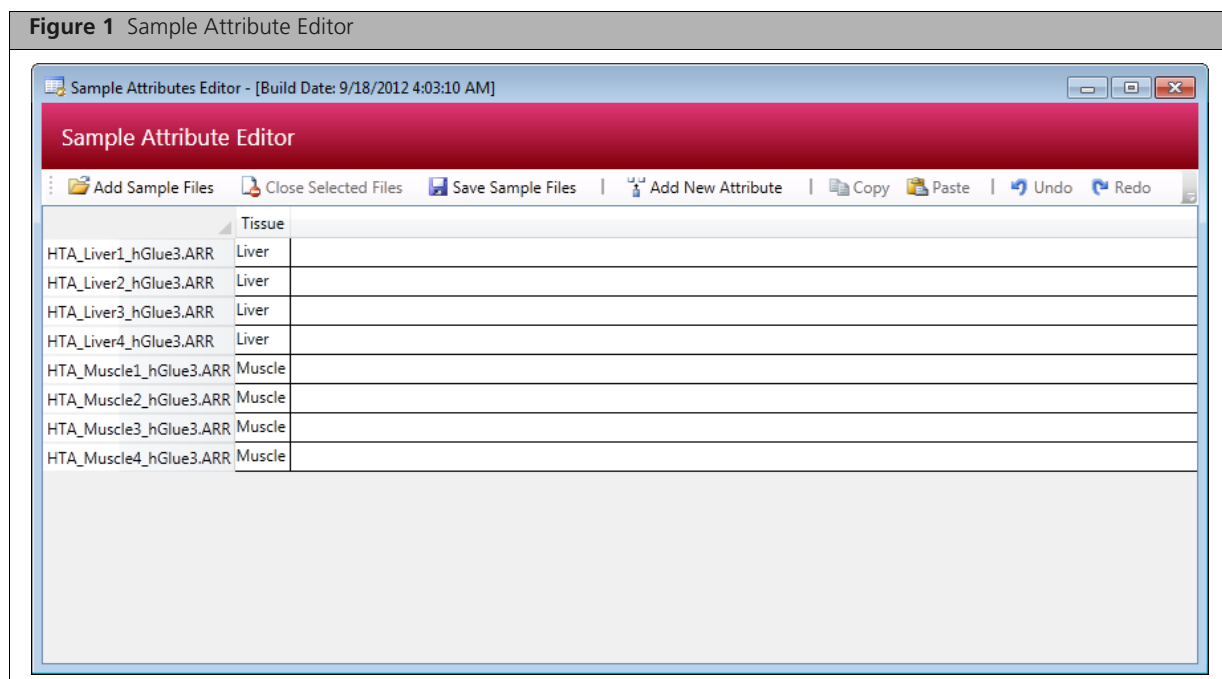


Table 1 Sample Attribute Editor Commands









Item	Description
 Add Sample Files	Opens a dialog box that enables you to select sample files (ARR) to add to the Sample Attribute Editor.
 Close Selected Files	Removes one or more selected ARR files from the Sample Attribute Editor.
 Save Sample Files	After editing sample information, click Save Sample Files to save the changes to the ARR files.

Table 1 Sample Attribute Editor Commands (Continued)

Item	Description
 Add New Attribute	Enables you to create anew column in the Sample Attribute Editor.
 Copy	Copies selected text to the system clipboard. For example: <ul style="list-style-type: none"> ■ Copy a cell and paste it to multiple selected cells ■ Copy multiple cells and paste to multiple cells.
 Paste	Pastes text from the system clipboard.
 Undo	Erases the previous change made in the Sample Attribute Editor (can be performed multiple times).
 Redo	Reverses the undo command (can be performed multiple times).

Adding or Removing Sample Files From the Sample Attribute Editor

To add files:

1. Click **Add Sample Files**.
2. Select a sample file(s) (ARR) from the dialog box that appears.

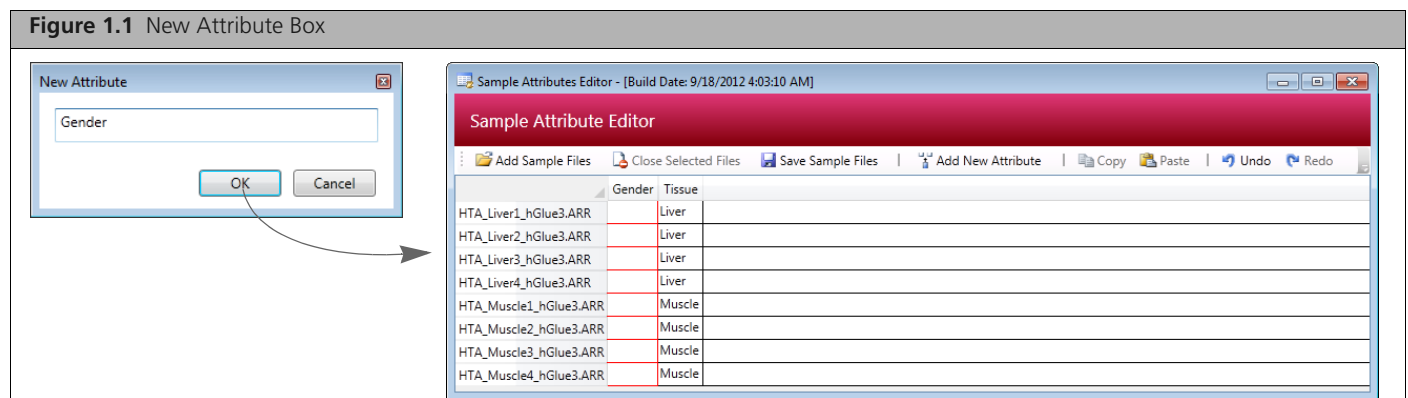
To remove files:

1. Select a file(s) in the Sample Attribute Editor.
2. Click **Close Selected Files**.

Adding, Editing, or Removing Attributes

To add a new attribute:

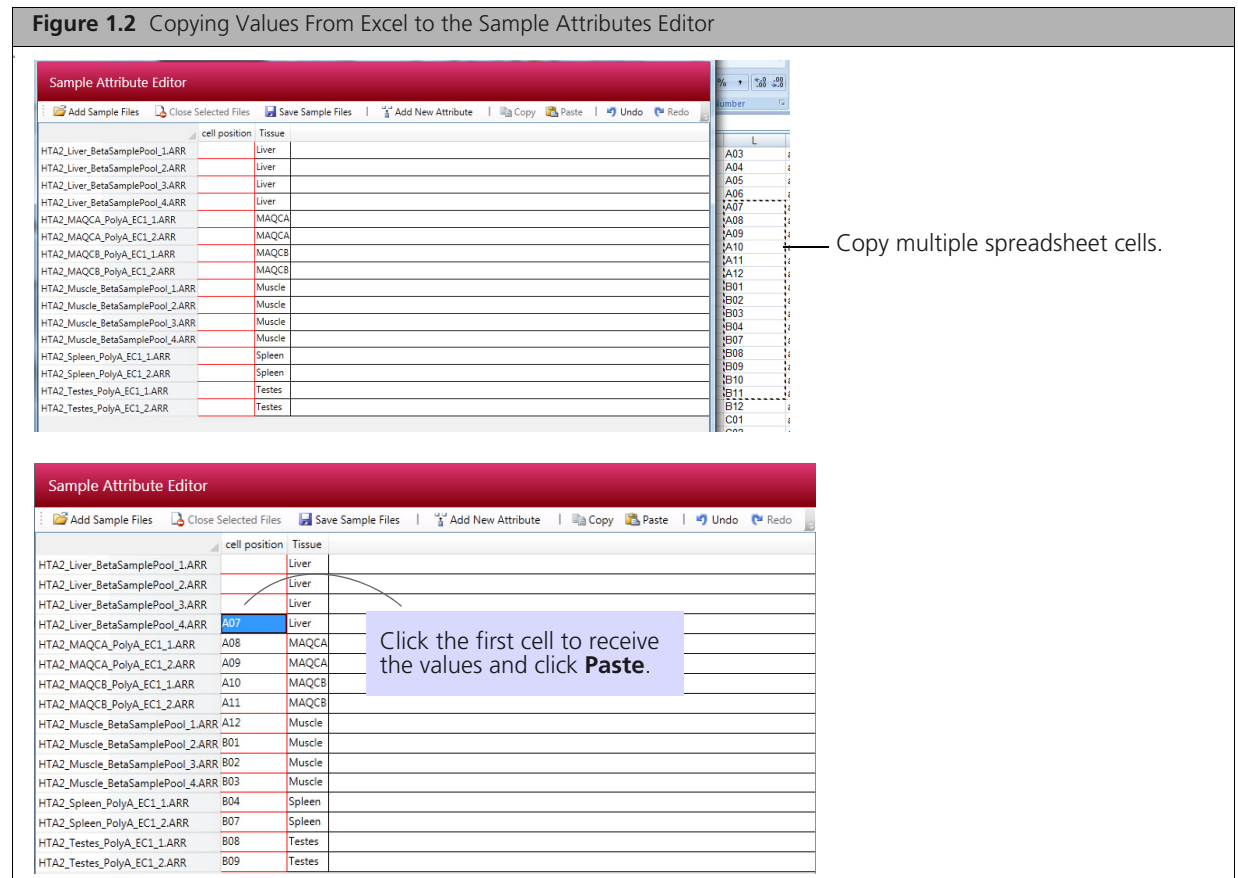
1. Click **Add New Attribute**. Alternatively, right-click an attribute column header and select **New Attribute** on the shortcut menu.
2. Enter a name for the attribute in the box that appears and click **OK**.
The attribute is added to the Sample Attribute Editor.



3. To enter attribute values:
 - A. Select one or more cells.
 - B. Enter text in the last cell in the selection and press **Enter**.
The same value is added to all of the selected cells.
4. Click **Save Sample Files**.

To copy and paste values from a spreadsheet:

1. Copy multiple cells in the spreadsheet.
2. In the Sample Attributes Editor, click the first cell to receive the values and click **Paste** (Figure 1.2).
3. Click **Save Sample Files**.



To edit an attribute value:

1. Double-click the cell.
2. Enter text or right-click the cell to view a shortcut menu of edit commands (Cut, Copy, Paste).

To sort attributes:

1. Click the column header
Attributes are sorted in ascending alphanumeric order.
2. Click the column header again to sort in descending alphanumeric order.

To remove an attribute:

1. Right-click the attribute column header and select **Delete Attribute** on the shortcut menu.
2. Click **Save Sample Files**.