

Using Variables to Override Function Settings in a Journal

Abstract

This document describes modifying a journal in the journal editor to enable individual journal functions to use variables or variable formulas instead of the recorded settings. This functionality requires version 6.1 or later of MetaMorph.

Instructions

Document ID
T20030

Product
MetaMorph

Created
17-Jun-2003

Last Reviewed
24-Sep-2004

Note: These procedures must be completed using the Journal Editor; they cannot be recorded.

This document describes the following procedures:

- Selecting a setting to be overwritten
- Assigning a new value to the setting

Note: This document assumes that the Journal drop-in has been loaded using the Meta Imaging Series Administrator. It also assumes that MMVAR has been loaded.

Selecting the Journal Function Containing Settings to be Overwritten

To select a journal function setting to be overwritten, complete the following steps:

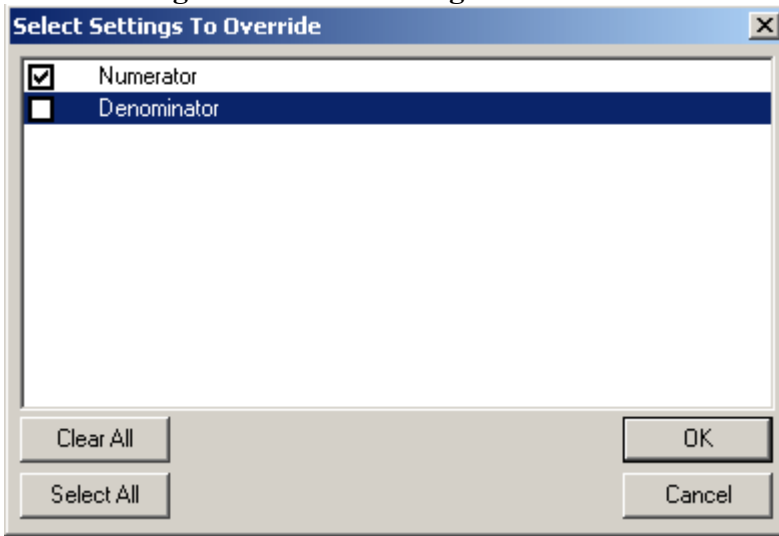
1. Select Edit Journal from the Journal menu. The Select a journal to edit dialog box opens.
2. Select or navigate to the journal you want to edit and click *Open*. The Journal Editor opens with the selected journal.
3. Select the journal function containing the parameters to override, then click *Select Settings to Override*. In this example, the *Descriptions* tab is selected and the *Arithmetic* function is used.

Note: The *Select Settings to Override* option is only available if the selected function can be overwritten.

The Select Settings to Override dialog box opens, as shown in Figure 1:

Keywords: variable journals
Issue Type: Application

Figure 1
Select Settings to Override Dialog Box



4. In the Select Settings to Override dialog box, click the check box next to each setting that you want to override, then click *OK*. The journal function changes in the journal editor to indicate that a parameter has been overridden, as shown below:

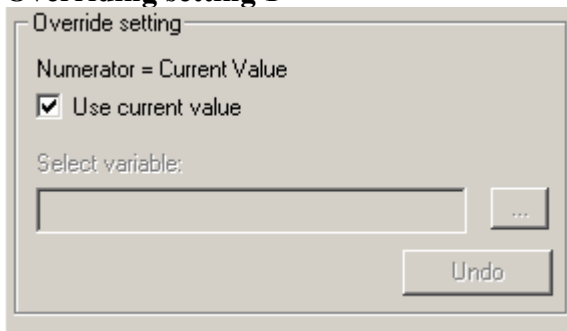
4: New "Add" = [Current At Start] + Overridden
 ↳ Numerator = Current Value

Assigning a New Value to the Function Setting

To override a setting in your journal function, complete the following steps:

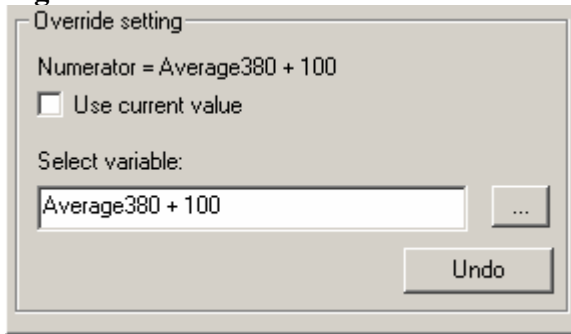
1. Select the line in the journal editor that contains "Current Value." The section below the Functions tab changes to *Override setting*, as shown in Figure 2:

Figure 2
Overriding setting 1




2. Uncheck *Use current value*. The *Select variable* field becomes enabled. Enter a variable or a variable expression to assign to the setting, as shown in Figure 3:

Figure 3



The Journal Editor will update to show the modified version of the journal function, as shown below:

```
1: Overwrite "Ratio" = ("Wavelength 340" * Overridden)/("Wavelength 380" * 1)
   ↳ Numerator = Average380 + 100
```

Note: If your expression requires the use of multiple variables, click the  button to open the Enter Variable Expression dialog box. Use this dialog box to create and assign a custom expression.