

Using Variables to Override Function Settings in a Journal

Abstract

This document describes how to use the journal editor to modify a journal 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 the Discovery-1 System software.

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.

Instructions

Selecting the Journal Function Containing Settings to be Overwritten

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


1. From the Journal editor, open the journal to be modified (Journal menu > Edit Journal).
2. Select the journal function for which you wish to override the parameters. If there are settings that can be overwritten, Select Settings to Override  will be active.
3. Click Select Settings To Override; the Select Settings to Override dialog opens.

Figure 1

Select Settings to Override

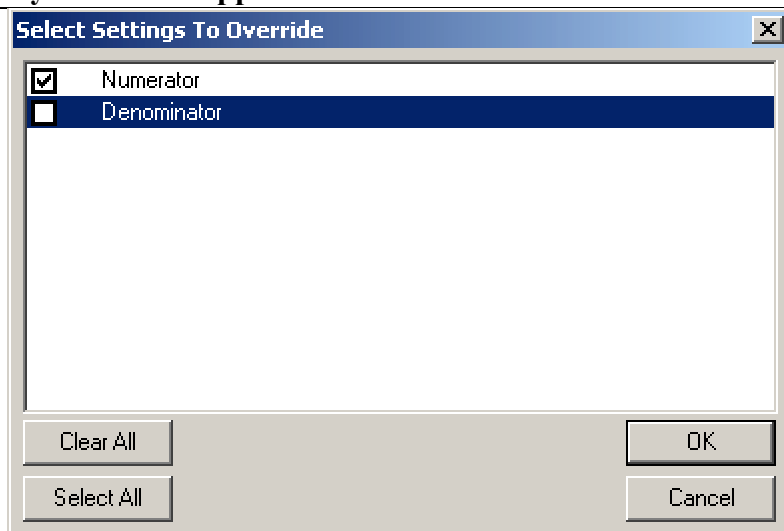
Document ID
D50008

Product
Discovery-1 version 6.1

Created
17-Jun-2003

Last Reviewed
29-Jul-2003

Keywords: variable journals v6.1 v6
Issue Type: analysisjournal



Assigning a New Value to the Function Setting

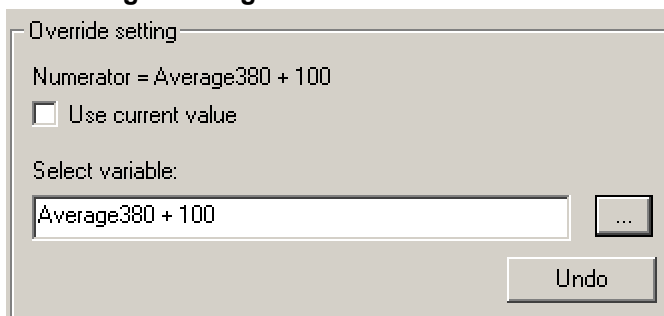
To override a setting in your journal function, complete the following steps:

1. In the Select Settings to Override 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.

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

2. Select the line in the journal editor that has "Current Value." The right bottom corner of the journal editor changes to Override setting.

Figure 2
Overriding a setting





3. Uncheck Use current value. The dialog will allow you to select a variable or a variable expression to assign to the setting. 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. The journal editor will show the assignment that has been made.

Figure 3
Final appearance of the journal function in the editor

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