



CFX Maestro™ SOFTWARE

Converting from CFX Manager™ to CFX Maestro Software: Naming Sample Groups

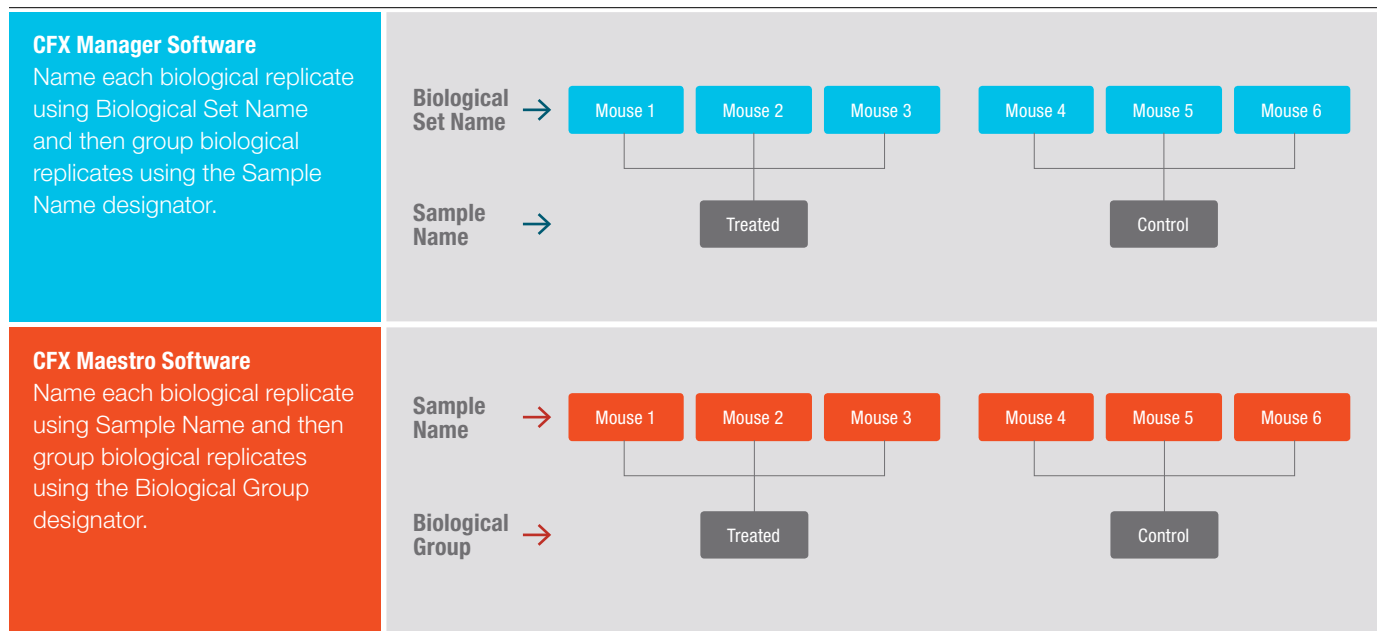
Quick Guide

This guide is only for users who are converting from CFX Manager Software v3.0 and 3.1 to CFX Maestro Software and who were using both Sample Name and Biological Set Name as designators in Plate Setup.

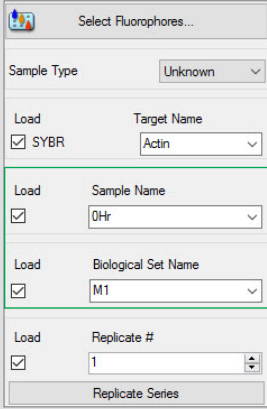
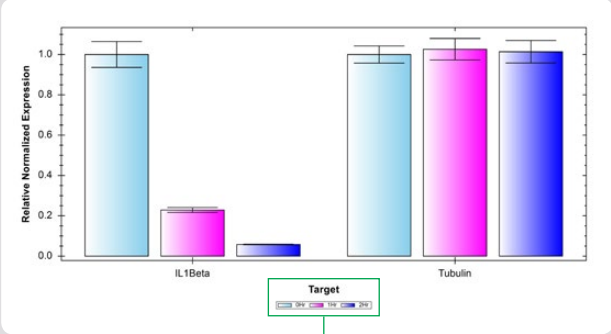
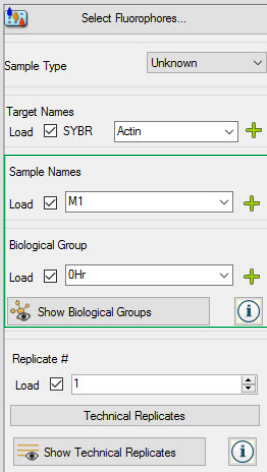
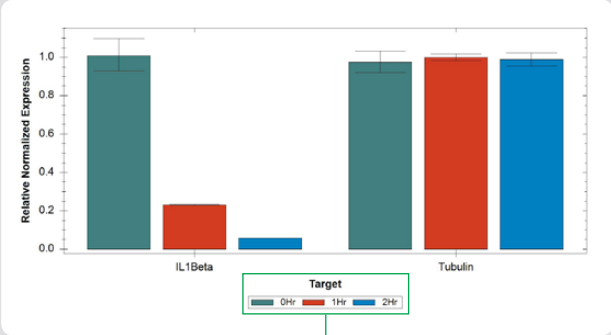
CFX Manager Software v3.0 and 3.1 allow for two hierarchical levels of naming and grouping unknowns: Sample Name and Biological Set Name. Biological Set Name can be used to designate each biological replicate (Mouse 1, Mouse 2, etc.) while Sample Name can be used to designate biological groups (Treated vs. Control).

CFX Maestro Software uses Sample Name to designate a single individual unknown or biological replicate, for example Mouse 1. Biological Group is now used to designate a group of samples (Treated vs. Control).

Comparison of Naming and Grouping Conventions



Comparison of Naming and Grouping Conventions (continued)

	Plate Setup	Data Analysis
CFX Manager Software		
CFX Maestro Software		

SYBR is a trademark of Life Technologies Corporation.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site bio-rad.com **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 177 **Belgium** 32 (0)3 710 53 00 **Brazil** 55 11 3065 7550
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 45 44 52 10 00 **Finland** 358 09 804 22 00
France 33 01 47 95 69 65 **Germany** 49 89 31 884 0 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300 **Israel** 972 03 963 6050
Italy 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 31 (0)318 540 666 **New Zealand** 64 9 415 2280
Norway 47 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 (0) 861 246 723
Spain 34 91 590 5200 **Sweden** 46 08 555 12700 **Switzerland** 41 026 674 55 05 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311
United Arab Emirates 971 4 8187300 **United Kingdom** 44 020 8328 2000

