



Addendum to BD FACSCalibur IFU

February 2008

Undo Function in BD CellQuest Pro Software

Overview

NOTE This document is an addendum to the *BD FACSCalibur Instructions for Use*. Keep this addendum with the *BD FACSCalibur IFU*.

During recent internal testing of BD CellQuest™ Pro software, we discovered that if the Undo function is utilized to undo a region deletion, the information in the Region Statistics view for the sample is randomly reordered, resulting in a mismatch of results and regions.

This defect is found in BD CellQuest Pro software version 4.02 with the OS 9 operating system and software versions 5.0.1, 5.1, 5.1.1, 5.2, and 5.2.1 with OS X operating systems.

The following pages describe how to avoid using the Undo function after deleting a region, and how to display the data correctly.

Redrawing a Region

As long as you do not use the Undo function to undo a region deletion, the software can be used with full functionality as designed without any mismatch of results.



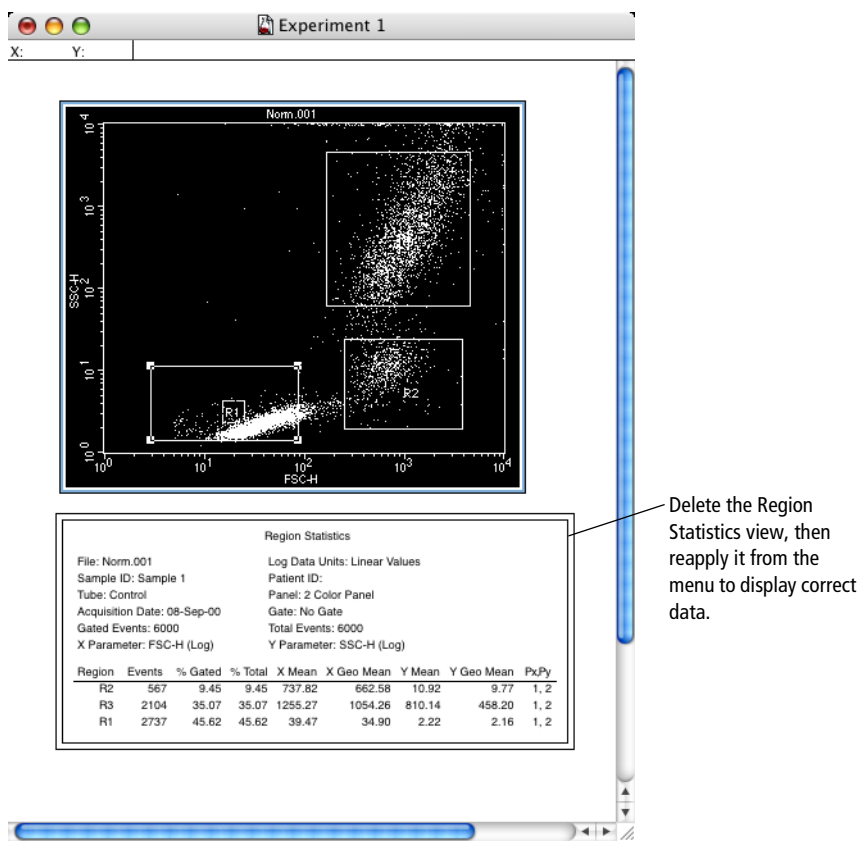
Do not use the Undo function to restore a region after deleting the region. The resulting data that is displayed will be incorrect.

If you deleted a region unintentionally, you need to re-draw the region instead of using the Undo function. Refer to the *BD CellQuest Pro Software Reference Manual* for the details on drawing a region.

Displaying Data Correctly

If you have deleted a region and then used the Undo function to restore the region, the data in the Region Statistics View will be incorrect. Follow this procedure to display the data correctly.

- 1 After deleting a region and using the Undo function to restore the region, select the Region Statistics View and delete it.



Experiment 1

X: Y:

Norm.001

SSC-H

FSC-H

R1

R2

R3

Region Statistics

File: Norm.001
Sample ID: Sample 1
Tube: Control
Acquisition Date: 08-Sep-00
Gated Events: 6000
X Parameter: FSC-H (Log)

Log Data Units: Linear Values
Patient ID:
Panel: 2 Color Panel
Gate: No Gate
Total Events: 6000
Y Parameter: SSC-H (Log)

Region	Events	% Gated	% Total	X Mean	X Geo Mean	Y Mean	Y Geo Mean	Px/Py
R2	567	9.45	9.45	737.82	662.58	10.92	9.77	1, 2
R3	2104	35.07	35.07	1255.27	1054.26	810.14	458.20	1, 2
R1	2737	45.62	45.62	39.47	34.90	2.22	2.16	1, 2

Delete the Region Statistics view, then reapply it from the menu to display correct data.

- 2 Display correct data by selecting Stats > Region Stats from the menu bar. The new Region Stats View now displays the data correctly for the regions in the plot.

Technical Assistance

If you would like more information, please call our Customer Support Center at 877.232.8995. Outside the US, call your local BD Biosciences office or distributor.

© 2008, Becton, Dickinson and Company.

BD, BD Logo and all other trademarks are property of Becton, Dickinson and Company. © 2008 BD

bdbiosciences.com
Part No. 644106 Revision A
February 2008



BD Biosciences

San Jose, CA 95131-1807
USA
Tel (877) 232-8995
Fax (800) 325-9637
facservice@bd.com

Asia Pacific

Tel (65) 6-861-0633
Fax (65) 6-860-1593

Europe

Tel (32) 2 400 98 95
Fax (32) 2 401 70 94
help.biosciences@europe.bd.com

Brazil

Tel (55) 11-5185-9995
Fax (55) 11-5185-9895

Japan

Nippon Becton Dickinson Company, Ltd.
Toll Free 0120-8555-90
Tel 81-24-593-5405
Fax 81-24-593-5761

Canada

Toll Free (888) 259-0187
Tel (905) 542-8028
Fax (888) 229-9918
canada@bd.com

Mexico

Toll Free 01-800-236-2543
Tel (52) 55 5999 8296
Fax (52) 55 5999 8288