

August 2012

BD FACSCanto™ II Flow Cytometer and BD FACSCanto™ Clinical Software FAQs

What are the flow rates for the BD FACSCanto II flow cytometer?

Low is 10 µL/min
Medium is 60 µL/min
High is 120 µL/min

BD FACSCanto Clinical Software

What is new in BD FACSCanto clinical software v2.4?

BD FACSCanto™ clinical software v2.4 now has 6-color TBNK without BD Trucount™ tubes as a panel, and you can specify a name for the results files.

CAUTION: Do not attempt to read FCS files created with version 2.4 into previous versions of BD FACSCanto clinical software. Previous versions will show incorrect results.

Is it possible to skip samples and run them later?

Skipping samples on a worklist can be done, but you cannot run them later in the same worklist. You will have to make a new worklist and run the samples from it.

Can process control sample results be exported in a single file?

Process controls can be run, and the results exported to a file. The Sample Name has to be control, and the Sample ID should be the lot ID for the control. Running the same lot again appends the new data to the same file. The file location is: C:\Program Files\BD FACSCanto Software\DataFiles. The file name is sample ID.csv.

BD FACS™ 7-color setup beads

What setup settings are printed out?

The setup results printout shows only the Lyse/Wash settings.

If I use optimized settings earlier, are they still applied?

If you have optimized your settings, the optimization is applied to future 7-color setup bead results within the worklist. Each optimization overwrites the previous .opt file.

BD™ Stem Cell Enumeration Kit

What version of the BD FACSCanto clinical software does the BD™ Stem Cell Enumeration kit require?
The BD™ Stem Cell Enumeration kit requires BD FACSCanto clinical software v2.4.



August 2012

BD FACSCanto™ II Flow Cytometer and BD FACSCanto™ Clinical Software FAQs

What samples can be run using BD FACSCanto clinical software?

BD Biosciences provides eight panels. You cannot edit or delete these panels. The number of sample tubes and the reagents associated with the tubes for each panel are shown in the columns to the right in the following table.

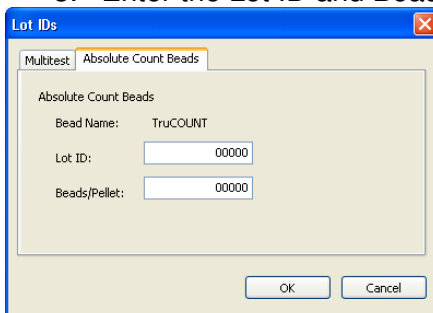
Panel	Tubes	Reagents
4 Color TBNK	1	CD3 FITC/CD8 PE/CD45 PerCP/CD4 APC
	2	CD3 FITC/CD8 PE/CD45 PerCP/CD4 APC
4 Color TBNK + TruC	1	CD3 FITC/CD8 PE/CD45 PerCP/CD4 APC + BD Trucount tube
	2	CD3 FITC/CD16+56 PE/CD45 PerCP/CD19 APC + BD Trucount tube
3/16+56/45/19	1	CD3 FITC/CD16+56 PE/CD45 PerCP/CD19 APC
3/16+56/45/19 + TruC	1	CD3 FITC/CD16+56 PE/CD45 PerCP/CD19 APC + BD Trucount tube
3/8/45/4	1	CD3 FITC/CD8 PE/CD45 PerCP/CD4 APC
3/8/45/4 + TruC	1	CD3 FITC/CD8 PE/CD45 PerCP/CD4 APC + BD Trucount tube
6 Color TBNK + TruC	1	CD3 FITC/CD16+56 PE/CD45 PerCP-Cy5.5/CD4 PE- Cy7/CD19 APC/CD8 APC-Cy7 + BD Trucount tube
6 Color TBNK	1	CD3 FITC/CD16+56 PE/CD45 PerCP-Cy5.5/CD4 PE- Cy7/CD19 APC/CD8 APC-Cy7

What are the default locations for the files?

File Type	Default location
FCS (.fcs)	D:\BDFACSCanto FCSFiles\yyyy\Month\dd
Levey-Jennings (.csv)	C:\Program Files\BD FACSCanto Software\QC
Acquisition Worklist (.wkl)	C:\Program Files\BD FACSCanto Software\Worklists
Analysis Worklist (.wka)	C:\Program Files\BD FACSCanto Software\Worklists
Setup Report (.pdf)	C:\Program Files\BD FACSCanto Software\SetupReports
Optimized Setup Result (.opt)	C:\Program Files\Common Files\BD\Setup Results
Result (.csv)	C:\Program Files\BD FACSCanto Software\DataFiles\yyyy\Month\dd
Process Control Results (.csv)	C:\Program Files\BD FACSCanto Software\DataFiles

How do I enter an Absolute Count Bead Lot ID?

1. Select **Tools > Lot IDs**.
2. In the **Lot IDs** dialog, click the **Absolute Count Beads** tab.
3. Enter the Lot ID and Beads/Pellet, which can be found on the pouch.



August 2012

BD FACSCanto™ II Flow Cytometer and BD FACSCanto™ Clinical Software FAQs

4. Click **OK**.

To obtain accurate results, it is critical that you enter this information accurately. Double-check your entry.

How do I hide the Reviewed By field?

You can hide the *Reviewed by* field that by default is displayed on Setup Reports. To hide this field, follow these steps.

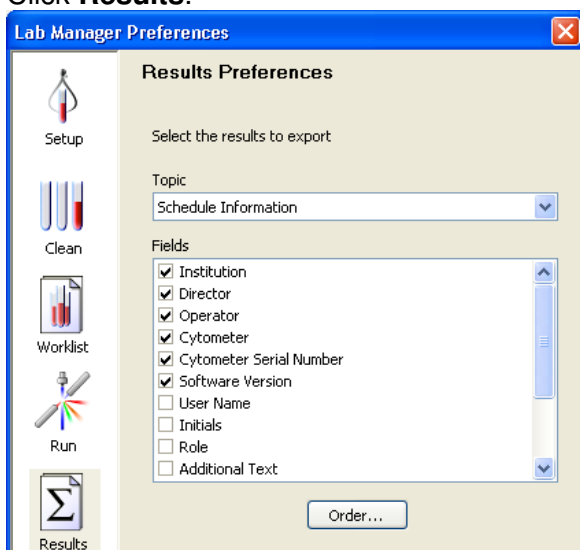
1. Select **Tools > Lab Manager Preferences**.
2. Click **Setup**.
3. Clear the **Review field on Setup Report** checkbox.
4. Click **OK**.

How do I set results preferences?

You can choose which results are exported to the results file that is created on a daily basis. First, specify the results and the order.

Caution: Changing the order affects the CSV output file. If the CSV result file is sent to an LIS, the file might not be read correctly.

1. Select **Tools > Lab Manager Preferences**.
2. Click **Results**.

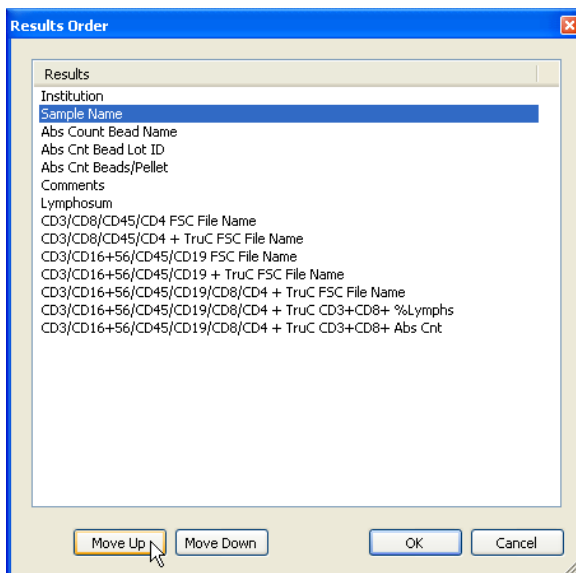


3. Select a topic from the menu.
4. Select or clear fields associated with the topic.
The information for selected fields will be exported.
The information for cleared fields will not be exported.

August 2012

BD FACSCanto™ II Flow Cytometer and BD FACSCanto™ Clinical Software FAQs

5. Select another topic.
6. Select or clear fields associated with the topic.
7. Continue until all needed topics and fields have been included for export.
8. Select the order for exported information.
 - a. Click **Order**.
 - b. Select a line of information.



- c. Click **Move Up** or **Move Down** until the information is in the preferred order.
- d. Repeat as needed.
- e. Click **OK** when you are done.

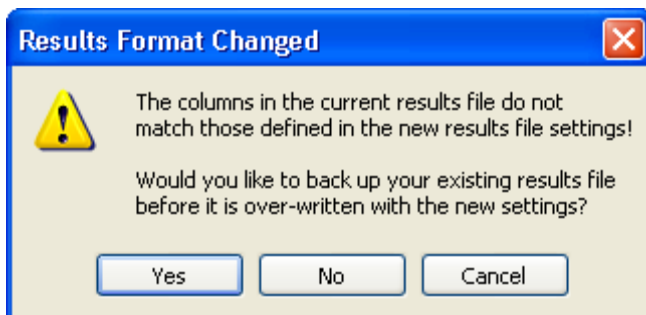
Next, save the results and back up the existing file, if necessary. When you change the type of information included in the results file, headers in the new file are different from the previous file. Because the new format is incompatible with the old one, you need to either back up your previous file and start a new file, or overwrite the previous file with the new one.

1. Click **OK** to save the new results for export.
2. If prompted, decide whether to back up your existing file.



August 2012

BD FACSCanto™ II Flow Cytometer and BD FACSCanto™ Clinical Software FAQs



- Click **Yes** to back up your existing file.
- Click **No** to overwrite your previous file with the new file.
- Click **Cancel** to return to the **Results Preferences** window.

I imported BD FACSCanto clinical software files into BD FACSDiva™ software. Where are the Name/ID and Case Number?

BD FACSCanto clinical software FCS files have to be imported as files. They cannot be imported as an experiment. The following table shows what happens to information from the original FCS files.

BD FACSCanto clinical field	BD FACSDiva field	Comments
Name	Specimen Name	
ID	\$SMNO in the keyword list	
Case Number	Case number	
Panel	Becomes a selectable keyword	The tube is the reagent name.
QC messages	QC codes	Cannot be displayed in BD FACSDiva software

All of the fields in the table can be displayed in the statistics view. In BD FACSDiva software, the population hierarchy displays the Tube name, # events, %Total, % Parent, and the population relationships. You can display the Specimen name, Tube name, and populations at the top of a plot. It is also possible to have custom titles. However, the Specimen name, Tube name, and population cannot be used to title the worksheet automatically.

Cy is a trademark of Amersham Biosciences Corp. Cy dyes are subject to proprietary rights of Amersham Biosciences Corp and Carnegie Mellon University and are made and sold under license from Amersham Biosciences Corp only for research and in vitro diagnostic use. Any other use requires a commercial sublicense from Amersham Biosciences Corp, 800 Centennial Avenue, Piscataway, NJ 08855-1327, USA.

