

BD FACSDiva Software Quick Reference Guide for the BD LSR II with HTS Option

This guide contains instructions for using BD FACSDiva™ software version 6.0 and later with BD™ LSR II flow cytometers equipped with the BD™ High Throughput Sampler (HTS) option.

Most of the features for running plate-based experiments on the BD™ HTS option are located in the Plate window. The following figure displays the Setup tab of the Plate window.

The screenshot shows the 'Plate - 96 Well - U bottom' window with the 'Setup' tab selected. The window is divided into several sections:

- Plate Setup Details:** Located at the top left, it includes 'Filter Setup Details' with checkboxes for 'Specimen type', 'Acquisition order', 'Specimen settings', 'First well in group', 'Specimen number', and 'Well settings'.
- Plate Information:** Located at the top right, it includes 'Throughput Mode' (High, Standard) and 'Plate Status' (Loader Status).
- Plate Layout:** The central 8x12 grid shows well types and settings. Callouts indicate specifying well types, creating compensation control wells, and applying cytometer settings.
- Loader Settings:** Located at the bottom right, it includes 'Sample Flow Rate (µL/sec)', 'Sample Volume (µL)', 'Mixing Volume (µL)', 'Mixing Speed (µL/sec)', 'Number of Mixes', and 'Wash Volume (µL)'. Callouts indicate specifying and customizing sample delivery, sample mixing, and between-well washing.
- List of specimens on the plate:** A list on the right side of the window shows 'Setup Controls_001', 'Compensation Controls', and 'Specimen_001'.



Helping all people live healthy lives

Workflow Overview

The following figure shows the steps for daily workflow using BD FACSDiva software.

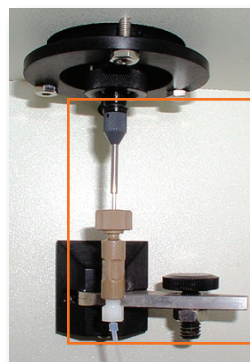


Before starting your daily workflow, ensure that your lab's software administrator has performed all the necessary tasks to set up the software for your use. This guide shows a workflow that uses application settings.

Starting Up the System

- 1 Start up the cytometer, the computer, and the HTS.
- 2 Prepare the fluidics tanks.
- 3 Verify that the optical filters are appropriate for your experiment.
- 4 Place the cytometer in run mode, start BD FACSDiva software, and log in.

The screenshot shows the 'Cytometer - LSR II (1)' status window. The 'Status' section is empty, and the 'Event' section shows 'Homing HTS loader...'. A callout box points to this event with the text 'The HTS initializes.'



After initialization, verify that the sample coupler is properly installed and not leaking.

- 5 Place the cytometer in standby mode.

Checking Cytometer Performance

- 1 Select Cytometer > CST.

Verify the Cytometer Configuration and bead Lot ID.

Clear the checkbox and select the plate type.

If needed, select a new configuration or bead lot ID.

- 2 Place the cytometer in run mode and run the BD™ Cytometer Setup and Tracking beads.
- 3 View the Cytometer Performance Report.
- 4 Close the Cytometer Setup and Tracking window.
- 5 Place the cytometer in standby mode.

Setting Up the Experiment

- 1 Create Browser elements.

Use the Browser toolbar to add elements.

- 2 Right-click Cytometer Settings in the Browser. Select Application Settings > Apply.

Select an application setting.

Name	Owner	Date Created
Application A	User1	07/05/07 12:52:24 PM
Application B	User1	07/05/07 12:53:28 PM

The application settings to be applied do not match the selected cytometer settings.

The following parameters are not in the selected application settings: PerCP-Cy5-5-A, Alexa Fluor 405-A, Alexa Fluor 430-A, DAPI-A, Indo 1 (Blue)-A.

Click Apply to apply PMT Voltage and Threshold values only for matching parameters.
Click Overwrite to replace all parameters and values with those from the selected application settings.

Click Overwrite if necessary.

- 3 Create setup control wells.

Use the Plate toolbar to add wells to the plate layout.

Select the throughput mode.

- 4 Select the first well for the compensation controls, right-click, and select Setup > Create Compensation Controls.

- 5 Create specimen wells.

Rename the specimen.

Verify that the loader settings are appropriate for your sample volume and recorded events.

- 6 Create a global worksheet.

Select Edit > Copy to copy the plots from the normal worksheet.


Select Edit > Paste to paste the plots to a global worksheet.

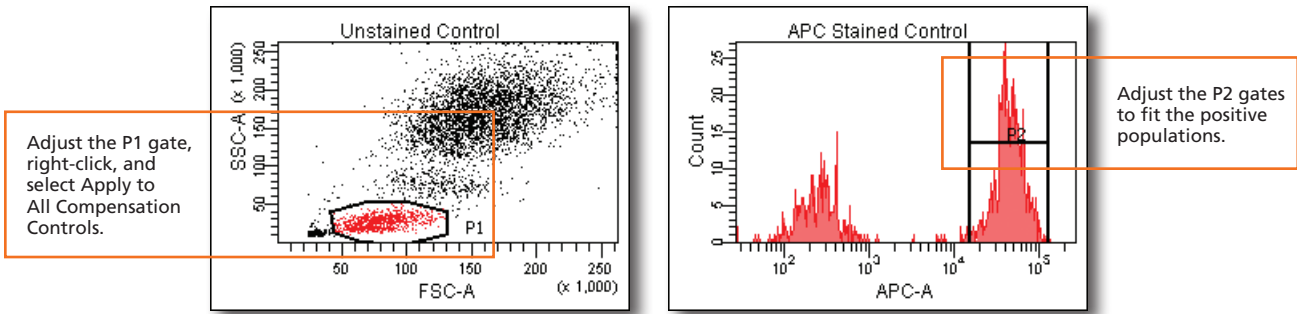
- 7 Install the prepared plate onto the HTS and place the cytometer in run mode.

- 8 Select the Setup Control well and click .

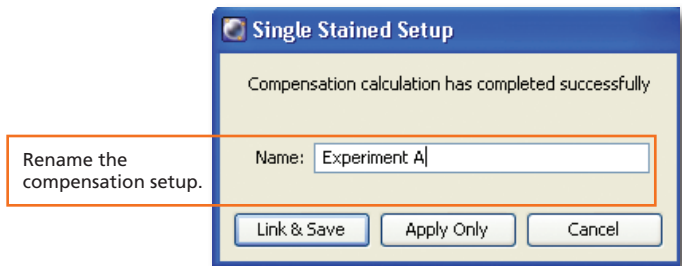
Verify that the FSC, SSC, and threshold settings are appropriate.

Parameter	Voltage	A	H	...
FSC	485	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
SSC	251	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>
FITC	466	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>
PE	479	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>
PE-Cy7	621	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>
APC	579	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>
APC-Cy7	568	<input checked="" type="checkbox"/>	<input checked="" type="checkbox"/>	<input type="checkbox"/>

- 9 Select all the compensation control wells and click  .
- 10 View the recorded data in the normal worksheets and gate the positive populations.

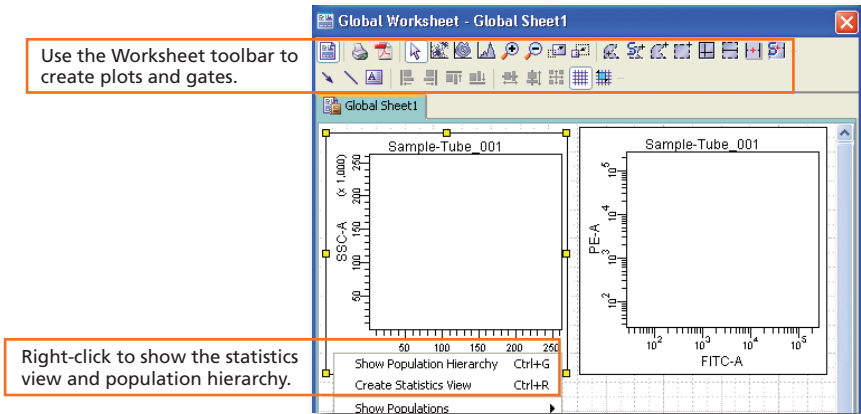



- 11 Select Experiment > Compensation Setup > Calculate Compensation.



Recording Specimen Data

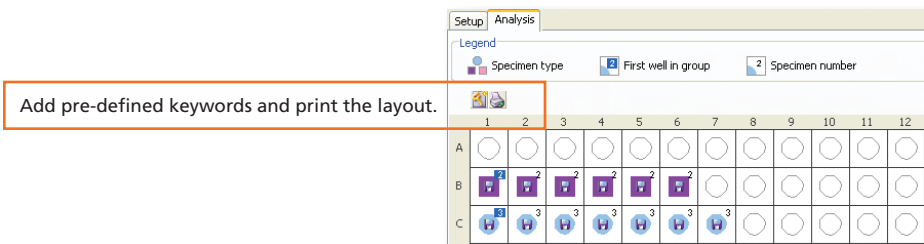
- 1 Create plots, gates, and statistics needed for recording.



- 2 Select the first specimen well and click  .
- 3 When recording is complete, place the cytometer in standby mode.

Analyzing Data

- 1 Under the Analysis tab of the Plate window, select a recorded well.



- 2 Create plots, gates, and statistics needed for analysis on a global worksheet.

Create new global worksheets.

Customize plots using the Plot Inspector.

Create custom text and graphics.

- 3 Perform quality control of the analysis.

Verify that gates are set appropriately for all samples.

Population	#Events	%Parent	%Total
All Events	10,000	###	100.0
Parent	1,867	18.7	18.7
Child A	128	6.9	1.3
Child B	219	11.7	2.2

Use the population hierarchy to verify parent/child relationships.

- 4 Right-click a specimen and select Batch Analysis.

Select to print, save as a PDF, or export the statistics as needed.

Shutting Down the System

- 1 Create a new experiment in the Browser.
- 2 Select HTS > Clean.

Select the Daily Clean template.

- 3 Install the prepared plate and click OK to begin cleaning.
- 4 Select File > Quit.
- 5 Turn off the cytometer and computer.

HTS Loader Settings Overview

HTS loader settings are specified under the Setup tab of the Plate window. Ensure that the loader settings are appropriate for your sample volume, sample concentration, and the specified events to record.

Default Loader Settings

Loader Settings	Loader Settings	Loader Settings	Loader Settings
Sample Flow Rate (µL/sec) 1.0	Sample Flow Rate (µL/sec) 1.0	Sample Flow Rate (µL/sec) 0.5	Sample Flow Rate (µL/sec) 1.0
Sample Volume (µL) 10	Sample Volume (µL) 3	Sample Volume (µL) 200	Sample Volume (µL) 10
Mixing Volume (µL) 100	Mixing Volume (µL) 50	Mixing Volume (µL) 100	Mixing Volume (µL) 100
Mixing Speed (µL/sec) 180	Mixing Speed (µL/sec) 200	Mixing Speed (µL/sec) 180	Mixing Speed (µL/sec) 180
Number of Mixes 2	Number of Mixes 2	Number of Mixes 2	Number of Mixes 2
Wash Volume (µL) 400	Wash Volume (µL) 200	Wash Volume (µL) 400	Wash Volume (µL) 400

Specimen wells using Standard Throughput mode

Specimen wells using High Throughput mode

Setup Control wells

Compensation Control wells

Loader Setting	Description	Important Considerations
Sample Flow Rate	Amount of sample (in µL per second) that is delivered to the flow cell. Select a rate between 0.5 and 3.0 in increments of 0.5 µL per second.	The larger the value entered, the shorter the plate running time, but this increases the sample core, causing more variation of data.
Sample Volume	Amount of sample (in µL) aspirated from the well and delivered to the flow cell. Select a volume between 2 and 200 µL.	For High Throughput mode, the system aspirates a set amount of 22 µL of sample, but records data for a volume between 2 and 10 µL. For Standard Throughput mode, the system aspirates the sample volume amount plus 20 µL. This value does not include the system default volume or the plate-dependent dead volume.
Mixing Volume	Amount of sample (in µL) aspirated and dispensed from the well to resuspend the particles.	To avoid introducing bubbles into the fluidics, this value should be half the total well volume.
Mixing Speed	Rate (in µL per second) that the mixing volume sample is aspirated and dispensed.	The faster the rate, the more likely that cell shearing occurs, especially for delicate cells. A faster rate can introduce bubbles in the sample delivered to the cytometer and compromise the separator bubble.
Number of Mixes	The number of times the mixing volume sample is aspirated and dispensed at the mixing speed. Select a number between 0 and 5 mixes.	The larger the number, the longer the plate running time.
Wash Volume	Amount of sheath fluid (in µL) drawn through the HTS fluidics between wells. Select a volume between 200 and 800 µL.	Enter a higher value to reduce cross contamination between wells. Enter a lower value to decrease the plate running time.