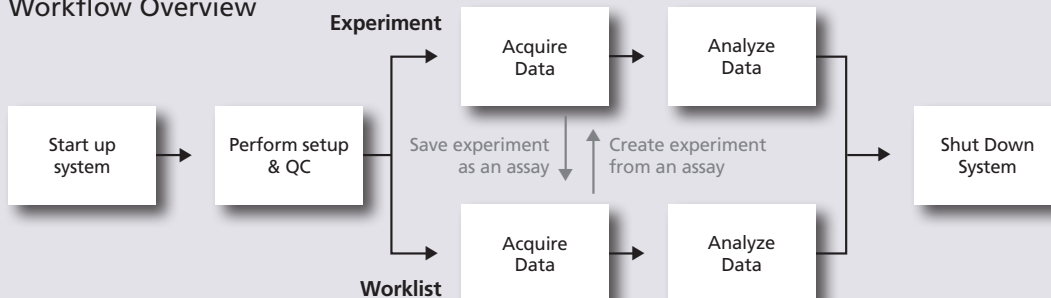


# BD FACSuite Software Quick Reference Guide for the Worklist Workflow

This guide contains instructions for using BD FACSuite™ software with the BD FACSVerser™ flow cytometer using the worklist workflow. Data can be acquired and analyzed in an experiment or in a worklist, as shown in the following diagram.

## Workflow Overview



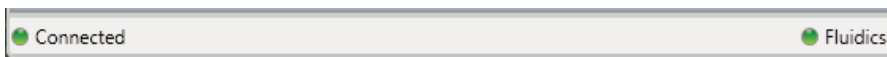
Before starting your daily workflow, ensure that your lab's software administrator has performed all the initial setup tasks as specified in the *BD FACSVerser System Reference*.



Helping all people  
live healthy lives

## Start up the system

- 1 Log in to BD FACSuite software.
- 2 Check the fluid levels.
- 3 Check the connection and fluidics status.



- 4 Verify that the 20-minute laser warmup has been completed.

# Performance QC and Assay Setup

## Performance QC

- 1 Open the **Setup & QC** workspace.

Select **Setup & QC**.

Verify the bead lot ID.

Click **Start**.

The image shows a 'Setup' menu with 'Setup & QC...' selected. To the right, the 'Setup & QC Options' dialog box is open, showing 'Task: Performance QC' and 'CS&T Bead Lot ID: 12161 (RUO, Expires: 5/31/2015)'. The 'Start' button is highlighted.

- 2 Run the BD FACSuite™ CS&T Research beads.
- 3 View the **Setup and QC Report** and **QC Tracking** tabs.

Last Performance QC (Normal Fluidics Mode):  Passed

Last Performance QC (High Sensitivity Fluidics Mode):  Passed

Verify that the Performance QC passed.

The image shows a summary of performance QC results with two 'Passed' status indicators.

## Assay and Tube Settings Setup

- 1 Select the **Setup & QC** tab.

Select the **Assay & Tube Settings Setup** task.

Click **Select**.

The image shows the 'Setup & QC Options' dialog box with 'Task: Assay & Tube Settings Setup' selected. The 'Assays & Tube Settings' section has a 'Select...' button highlighted.

- 2 Select the **Assays** tab.

Select the assays that you want to set up.

Completed	Name	Type	Author	Tube Settings
<input checked="" type="checkbox"/>	4 Color Beads: UD	UserDefined	BDAdministrator	4 Color Beads

The image shows a table with one row of assay data. The 'Completed' checkbox is checked.

- 3 Run the BD FACSuite CS&T Research beads.

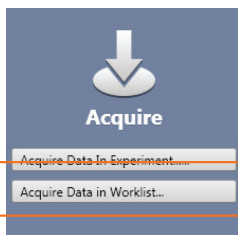
Verify the Lot ID.

Click **Start**.

The image shows the 'Assay & Tube Settings Setup' dialog box with 'Task: Assay & Tube Settings Setup' and 'CS&T Bead Lot ID: 12161 (RUO, Expires: 5/31/2015)'. The 'Start' button is highlighted.

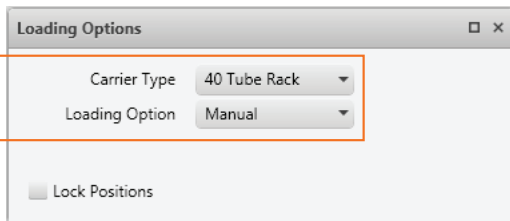
## Acquire and Analyze Data

- 1 Open the **Worklist** workspace.



Select **Acquire Data In Worklist**.

- 2 Modify the loading options, as needed.



Select the Carrier Type and Loading Option.

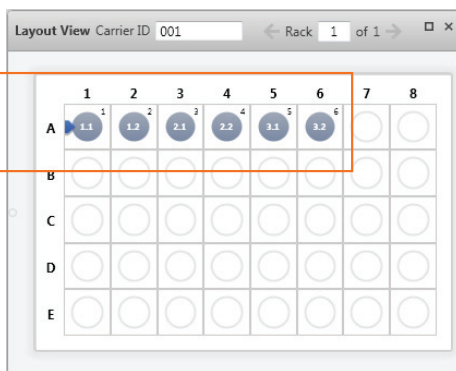
- 3 Create Worklist entries.

	Sample ID	Task	Status	Location	Sample Carrier
▶ 1	Donor 1	6-color Stimulation UD	Ready	A1-A2	001
▶ 2	Donor 2	6-color Stimulation UD	Ready	A3-A4	001
▶ 3		Perform Daily Cleaning	Ready	A5-A6	001

Enter a Sample ID.

Under **Task**, select an assay or fluidics task.

- 4 If using the Loader, place the sample carrier on the Loader tray. Click **Load**.



Place your tubes in the positions designated in the **Layout View**.

- 5 Click **Run All**.



6 Verify the gates, PMT voltages, and threshold values.

To make modifications:

- a. Click **Stop Timer** to pause the acquisition delay timer.

Click **PMTV** to enable the sliders in the worksheet to make adjustments to the PMT voltages.

Verify that the gates and populations are set appropriately.

Verify that the voltage and threshold settings are appropriate.

- b. Click **Resume**.

7 Perform quality control of the analysis report.

To make modifications:

- a. Click **Stop Timer** to pause the acquisition delay timer.

Verify that the gates and populations are set appropriately.

- b. Click **Resume**.

## Shut Down the System

- 1 Open a worklist.

	Sample ID	Task	Status
▶ 1		Perform Daily	Ready
▶ 2		Shutdown	Ready

If needed, add the daily clean and shutdown tasks to the worklist.

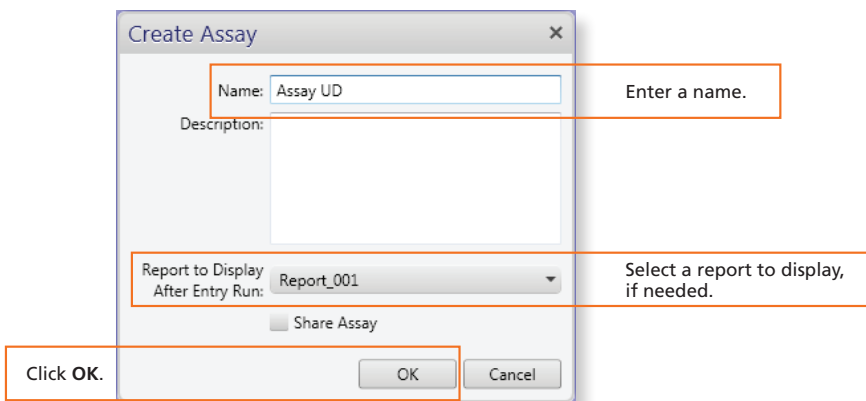
- 2 Run the worklist.
- 3 Log out of the software, if necessary.

# Additional Information

Any experiment can be used to create a user-defined assay and used in a worklist to acquire and analyze data. Additionally, experiments can be created from assays and used to acquire and analyze data in the Experiment workspace or saved as an assay to be used in a worklist.

## Creating an Assay from an Experiment

- 1 Open an existing experiment or create a new one.
- 2 Select **File > Create Assay**.



## Creating an Experiment from an Assay

- 1 Select **File > New Experiment from Assay**.
- 2 Modify the assay as needed.
- 3 Select **File > Create Assay** to save the experiment as an assay and acquire data in a worklist. Alternatively, you can continue with data acquisition and analysis in the **Experiment** workspace.