

abcam

# BD Accuri C6 with plate handler

Protocol for multiplex assays

April 2016

This protocol contains instructions for setting up an Accuri C6 with plate handler. If you will be using a different system, consult the appropriate protocol in our flow cytometry protocols for Multiplex assays page, or contact our Technical Support team at [technical@abcam.com](mailto:technical@abcam.com).

For first time use of a system, please follow the cytometer setup/verification. After a system is properly validated, those settings can be saved and used for routine assay readout with the method for assay readout.

**Please read this protocol carefully before cytometer operation. Failure to properly set up a system may result in unusable data, loss of product or system damage. For more detailed instructions on system operation details. Please consult the Accuri C6 operation manual.**

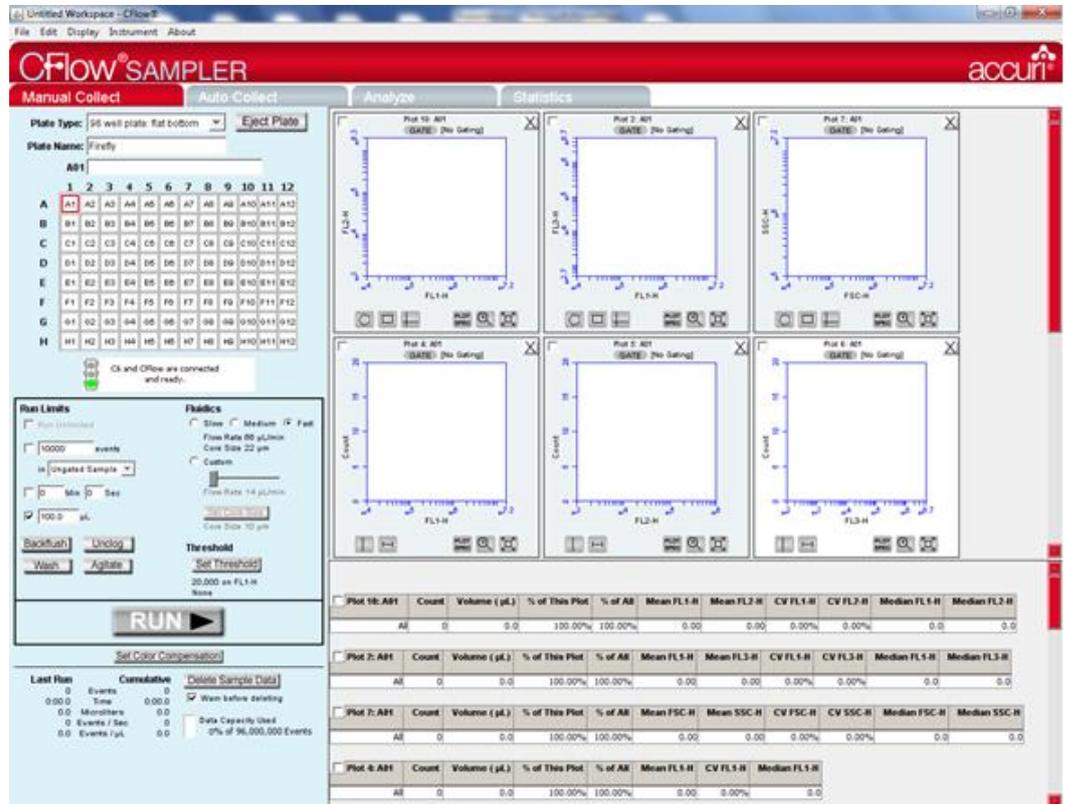
## Requirements

- Multiplex Immunoassay Core Kit OR Multiplex miRNA Assay Core Reagent Kit
- Blade (scalpel or scissors)
- Pipette and tips
- Tungsten cleaning wire (only if system SIP is obstructed)

## Cytometer setup/verification (first time use only)

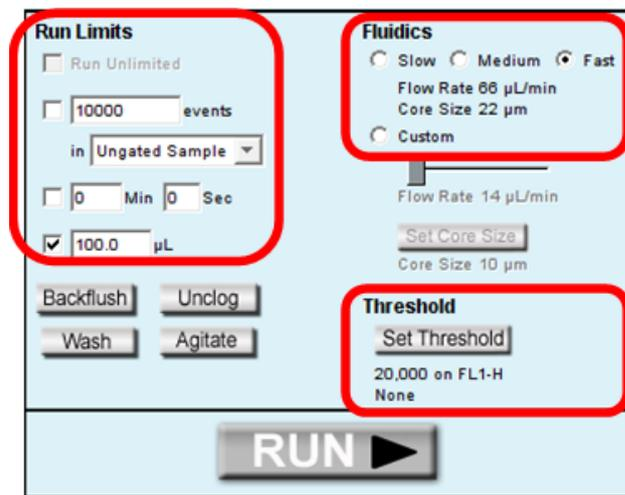
1. If using a Multiplex Immunoassay Core Kit follow the below instructions:
  - Prepare a 1x Capture Particle Master Mix according to the Multiplex Immunoassay Core Kit Booklet.
  - Cut and remove the plate seal from well H12 on the provided filter plate.
  - Add 75µL of 1x Capture Particle Master Mix to the well, and apply vacuum using a vacuum manifold to remove the buffer from the well.
  - Add 100µL of Run Buffer to the well.
2. If using a miRNA Core Reagent Kit follow the below instructions:
  - Cut/remove the plate seal from well H12 on the provided filter plate.
  - Add 35µL of Firefly™ Particles to the well, and apply vacuum to remove all liquid from the well.
  - Add 140µL of Run Buffer to the well
3. Download the Accuri Plate template file from [www.abcam.com/FireflyCytometry](http://www.abcam.com/FireflyCytometry).

4. Start up the Accuri System and open the CFlow Software. Be sure to perform all requisite system startup and cleaning functions for the cytometer.
5. Under the "File Menu", select "Open CFlow File or Template". Select the Accuri template file to load. Once the template has loaded, six plots should be displayed on your screen as illustrated below.



6. Locate the Acquisition Control section located in the middle-left section of the screen. Ensure that the following settings are selected:
  - a. **Threshold** is set to "20,000 on FL1-H".
  - b. **Run Limits** is set to "100 µL".
  - c. **Fluidics** is set to "Fast".

Failure to load the correct settings file will lead to loss of data.



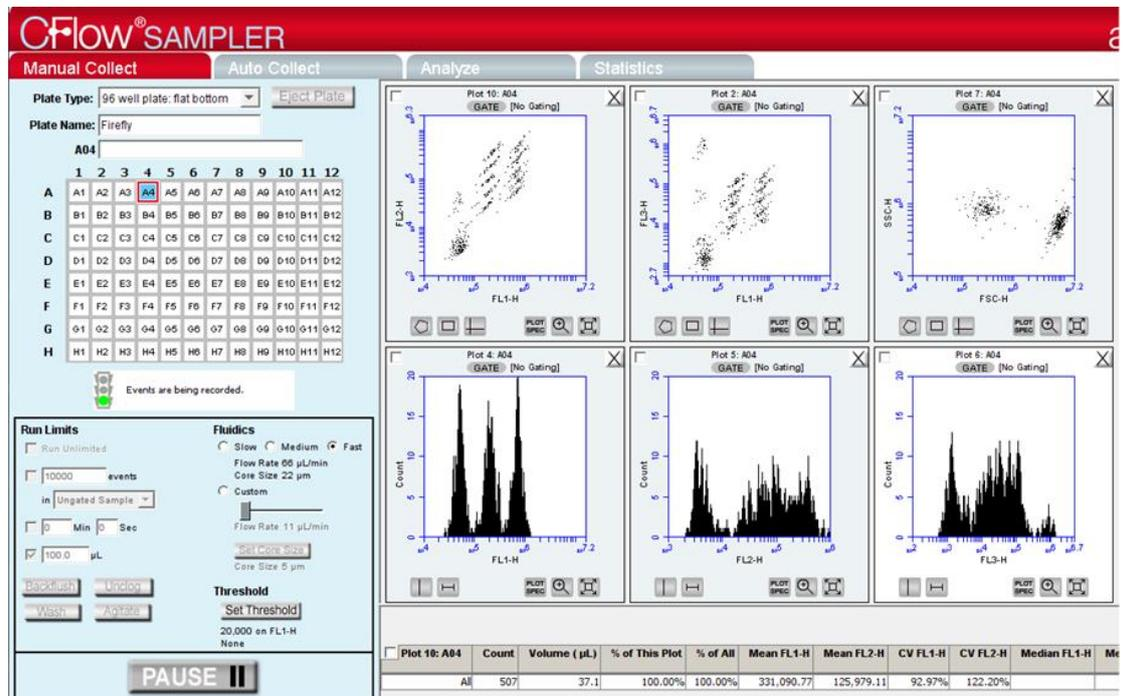
- Click the **"Eject Plate"** button in the top left corner of the control screen to load the plate.

**Touching the plate handler before it has stopped moving may cause the arm to require recalibration.**

- With the arm still in the ejected position, place a Kimwipe under the Sample Injection Port (SIP) tube and click on the **"Backflush"** button. Observe flow out of the SIP tube. The flow from the instrument should briefly be a steady and continuous stream for part of the procedure versus a closely spaced set of distinct droplets. The flow will only be a stream for around one second of the backflush if the instrument is working properly.
  - If the flow is a continuous drip without a period of steady flow insert a tungsten cleaning wire (VWR 72310-068) approximately 2 inches into the SIP to clean the flow cell. Repeat the backflush to see if the flow has improved to a steady stream.
  - If performance does not improve after cleaning the flow cell, ensure peristaltic pump tubing has been changed as per manufacturer recommendations.

**Proceeding without a properly flowing instrument will yield poorly resolved codes with potential data loss.**

- Load the plate in the handling arm, aligning the **"A1"** well of your plate with the indicated position on the plate platform. Click the **"Load Plate"** button to retract the arm into the instrument.
- Select the well in the plate grid and click the **"Run"** button directly below the Acquisition Control panel. When prompted, save the .c6 file to a hard drive location and the data acquisition will begin. Data acquisition should appear similar to the image below.



11. After the well has been acquired, you must export the data as an FCS file to be able to analyze in the Firefly™ Analysis Workbench. Select the well, then go to “**File Menu**” and select “**Export All Files as FCS**”. Locate the FCS file and transfer it to a USB drive to import it to your personal computer.
12. Download the Firefly™ Analysis Workbench onto your personal computer from <http://www.abcam.com/kits/firefly-analysis-workbench-software-for-multiplex-mirna-assays>. Reference your Core Kit instructions for detailed instructions on how to use the Analysis Workbench software.
13. Open the Firefly™ Analysis Workbench and load the FCS file saved in Step 11. When prompted, load the PLX file for your particle mix. Once the data loads, check the color of the well, a Green well indicates a successful setup.

**If the quality indicator is Red, perform a thorough system clean, and repeat this protocol. If the score remains low, contact our Technical Support team at [technical@abcam.com](mailto:technical@abcam.com).**

# Multi-well Auto Collect sample acquisition

**Important – only run this protocol after you have performed “Cytometer Setup/Verification” protocol on your BD Accuri C6 cytometer using particles. This protocol also assumes a completed assay in a 96-well plate, with particles suspended in Run buffer.**

1. Place your plate on a shaker for one minute to resuspend the particles. After shaking, particles are stable in suspension for 30 minutes before scanning.
2. Start up the Accuri System and open the CFlow software. Be sure to perform all requisite system startup and cleaning functions.
3. Under the “**File Menu**”, select “**Open CFlow File or Template**”. Select the Accuri template file to load. Once the template has loaded, six plots should be displayed on your screen as illustrated in Step 5 of the Cytometer Setup/Verification protocol.

**Failure to load the correct settings file will lead to loss of data.**

4. Click the “**Eject Plate**” button in the top left corner of the control screen to load the plate.

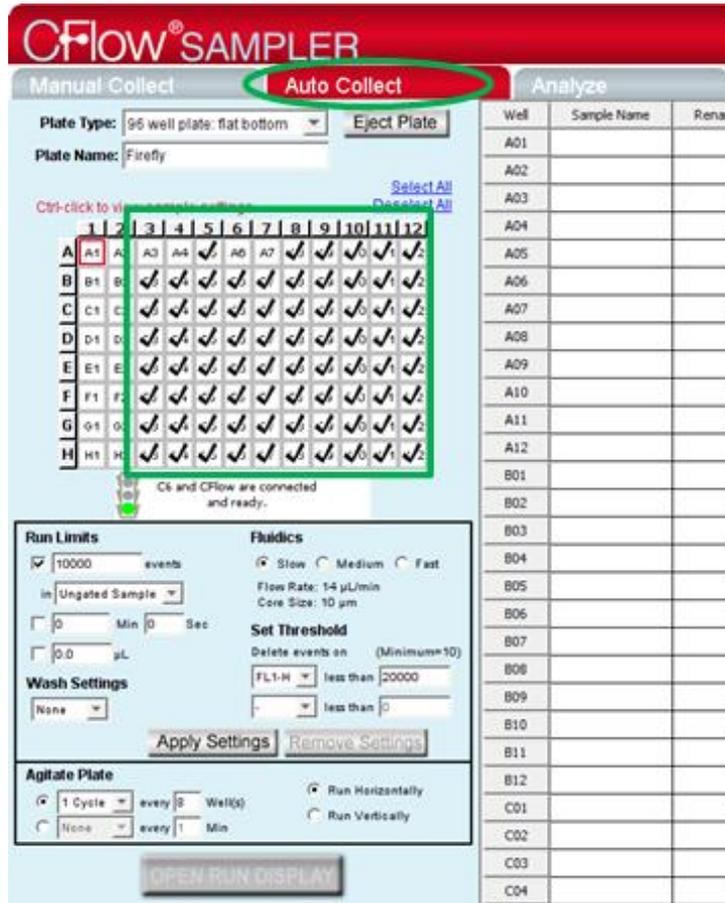
**Touching the plate handler before it has stopped moving may cause the arm to require recalibration.**

5. With the arm still in the ejected position, place a Kimwipe under the Sample Injection Port (SIP) tube and click on the “Backflush” button. Observe flow out of the SIP tube. The flow from the instrument should briefly be a steady and continuous stream for part of the procedure versus a closely spaced set of distinct droplets. The flow will only be a stream for around one second of the backflush if the instrument is working properly.
  - a) If the flow is a continuous drip without a period of steady flow insert a tungsten cleaning wire (VWR 72310-068) approximately 2 inches into the SIP to clean the flow cell. Repeat the backflush to see if the flow has improved to a steady stream.
  - b) If performance does not improve after cleaning the flow cell, ensure peristaltic pump tubing has been changed as per manufacturer recommendations.

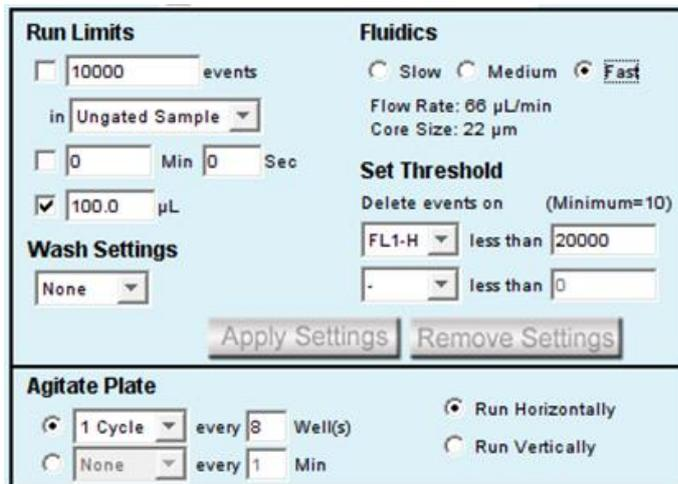
**Proceeding without a properly flowing instrument will yield poorly resolved codes with potential data loss.**

6. Load the plate in the handling arm, aligning the “**A1**” well of your plate with the indicated position on the plate platform. Click the “**Load Plate**” button to retract the arm into the instrument.

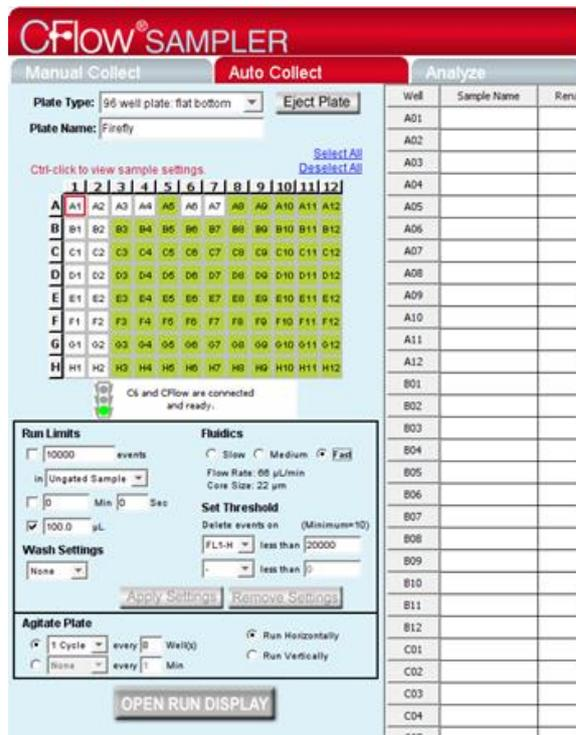
7. Click on the “Auto Collect” tab and select the appropriate assay wells to be acquired; selected wells will be highlighted with a check mark as illustrated below.



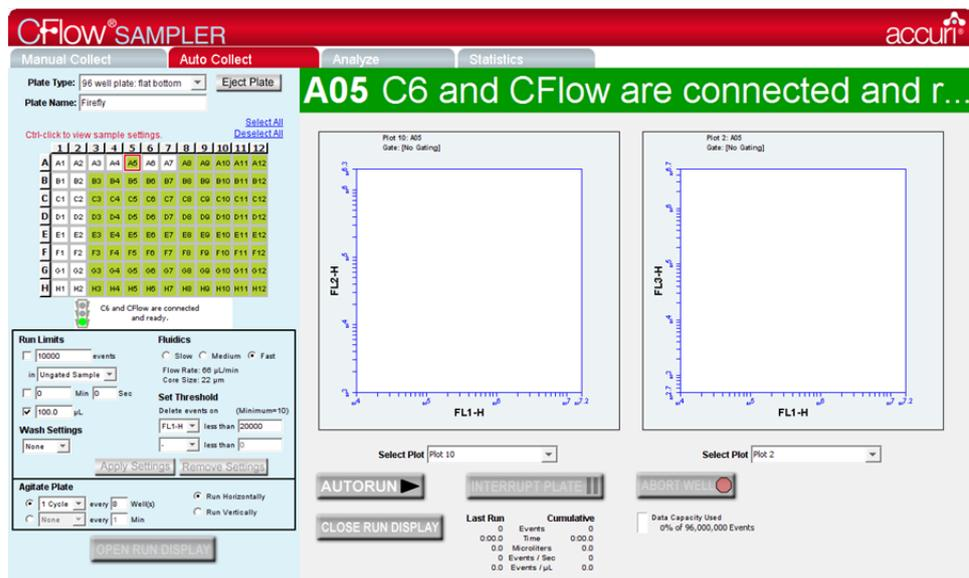
8. Locate the Acquisition Control section located in the middle-left section of the screen. Ensure that the following settings are selected:
- Threshold is set to “20,000 on FL1-H”.
  - Run Limits is set to “100 µl”.
  - Fluidics is set to “Fast”.
  - Agitate Plate is set as “1 Cycle every 8 wells”.



- In the Acquisition Control section, click the **“Apply Settings”** button and then save your \*.c6 experiment file to the hard drive. The check marked wells will turn Green if they have been properly selected for acquisition as illustrated below.



- Click the **“Open Run Display”** button located underneath the Acquisition Control. Two plots will appear the right of the plate layout.



- Again, ensure that the Acquisition Controls display the following settings before proceeding.

The screenshot shows the following settings in the CFlow Sampler software:

- Run Limits:** 10000 events, Ungated Sample, 0 Min 0 Sec, 100.0 µL.
- Fluidics:** Slow, Medium, **Fast** (selected), Flow Rate: 66 µL/min, Core Size: 22 µm.
- Set Threshold:** Delete events on (Minimum=10), FL1-H less than 20000, - less than 0.
- Wash Settings:** None.
- Agitate Plate:** 1 Cycle every 8 Well(s), Run Horizontally (selected), Run Vertically.

- Click the “**Auto Run**” button located beneath the first plot. The cytometer will now begin automatic acquisition of the first selected well. Once the well has been acquired, it will be marked with a small blue square and the sampler will move on to the next well.

The screenshot shows the CFlow Sampler software interface with the following elements:

- Header:** CFlow SAMPLER, accuri.
- Navigation:** Manual Collect, **Auto Collect**, Analyze, Statistics.
- Plate Information:** Plate Type: 96 well plate, flat bottom; Plate Name: Firefly.
- Well Plate Grid:** A 96-well plate grid (A1-A12, 1-12) with a 'Preparing to analyze sample' indicator.
- Settings Panel:** Run Limits, Fluidics, Set Threshold, Wash Settings, Agitate Plate (1 Cycle every 8 Well(s), Run Horizontally).
- Plots:** Two plots titled 'A08 Preparing to analyze sample.' Plot 10 (FL2-H vs FL1-H) and Plot 2 (FL3-H vs FL1-H).
- Buttons:** AUTORUN, INTERRUPT PLATE, ABORT WELL, CLOSE RUN DISPLAY, OPEN RUN DISPLAY.
- Statistics:** Last Run and Cumulative statistics for Events, Time, Microliters, Events / Sec, and Events / µL.

- After all samples have been acquired, click the “**Close Run Display**” button.
- Go to “**File Menu**” and select “**Export All Files as FCS**”. Locate the FCS file and transfer it to a USB drive to import it to your personal computer. Load the exported FCS file into the Firefly™ Analysis Workbench for analysis.

File Edit Display Instrument About

Open CFlow File or Template  
 New CFlow File  
 Save  
 Save CFlow File as...  
 Save CFlow Template as...  
 Auto-save Settings...  
 Import FCS File...  
 Export FCS File...  
**Export ALL Samples as FCS...**  
 Export ALL Samples to Third Party...  
 Export Plot Data as CSV...  
 Export Sample Settings...  
 Print Selected Items   
 Quit

**ER**

to Collect Analyze

Load Plate

Select All  
 Deselect All

Well	Sample Name	Reana
A01		
A02		
A03		
A04		
A05		
A06		
A07		
A08		
A09		
A10		
A11		
A12		
B01		
B02		
B03		
B04		
B05		
B06		
B07		
B08		
B09		
B10		
B11		
B12		
C01		
C02		
C03		
C04		

9 | 10 | 11 | 12

A9	A10	A11	A12
B9	B10	B11	B12
C9	C10	C11	C12
D9	D10	D11	D12
E9	E10	E11	E12
F9	F10	F11	F12
G9	G10	G11	G12
H9	H10	H11	H12

C6 and CFlow are connected and ready.

**Run Limits**  
 10000 events  
 in Ungated Sample  
 0 Min 0 Sec  
 100.0 µL

**Fluidics**  
 Slow  Medium  Fast  
 Flow Rate: 66 µL/min  
 Core Size: 22 µm

**Wash Settings**  
 None

**Set Threshold**  
 Delete events on (Minimum=10)  
 FL1-H less than 20000  
 - less than 0

Apply Settings Remove Settings

**Agitate Plate**  
 1 Cycle every 8 Well(s)  
 None every 1 Min  
 Run Horizontally  
 Run Vertically

OPEN RUN DISPLAY