

Analyzing Data Collected on the BD Accuri™ C6 Using FlowJo™

Koen Verbrugghe, PhD
Marketing Applications Specialist
BD Biosciences



The BD Accuri C6 Flow Cytometer System

- An affordable, full-featured, easy-to-use flow cytometer
- Two lasers and six detectors



Detection: Wide Dynamic Range



Advantages of Pre-optimized Voltage

- Greatly reduces the risk of lost data due to improper setup
 - Saves time and sample
- No specialist training or dedicated operator required
- Predictable, reproducible analysis relative to the sample type and application
- Attenuation filters (for bright signals) give controlled signal reduction
- Predictable fluorescence spillover
- Focus on the science of measuring fluorescence, not the art of setting voltages

Intuitive Software

Sample Grid

Cytometer Status

Fluidics Controls

Run Criteria

Real-Time Updates

Collect
Analyze
Statistics

A4 HPB CD3-F CD4, CD45, CD8

	1	2	3	4	5	6	7	8	9	10	11	12
A	A1	A2	A3	A4	A5	A6	A7	A8	A9	A10	A11	A12
B	B1	B2	B3	B4	B5	B6	B7	B8	B9	B10	B11	B12
C	C1	C2	C3	C4	C5	C6	C7	C8	C9	C10	C11	C12
D	D1	D2	D3	D4	D5	D6	D7	D8	D9	D10	D11	D12
E	E1	E2	E3	E4	E5	E6	E7	E8	E9	E10	E11	E12
F	F1	F2	F3	F4	F5	F6	F7	F8	F9	F10	F11	F12
G	G1	G2	G3	G4	G5	G6	G7	G8	G9	G10	G11	G12
H	H1	H2	H3	H4	H5	H6	H7	H8	H9	H10	H11	H12

C6 is connected and ready.

Run Limits

Run Unlimited

125000 events

in **Ungated Sample**

0 Min 0 Sec

0.0 μ L

Backflush **Unlog**

Fluidics

Slow Medium Fast

Flow Rate 14 μ L/min

Core Size 10 μ m

Custom

Flow Rate 11 μ L/min

Set Core Size

Core Size 5 μ m

Threshold

Set Threshold

None

80,000 on FSC-H

ADD to A4

Set Color Compensation

Last Run **Cumulative** **Delete Sample Data**

0 Events 229,303 Warn before deleting

0.00 Time 0.00

0.0 Microliters 0.0

0 Events / Sec 0

0.0 Events / μ L 0.0

Data Capacity Used
8% of 10,000,000 Events

Plot 1: Sample HPB CD3-F CD4, CD45, CD8

GATE: [No Gating]

Plot 2: Sample HPB CD3-F CD4, CD45, CD8

GATE: (R1 in (P1 in all))

Plot 3: Sample HPB CD3-F CD4, CD45, CD8

GATE: (P1 in all)

Plot 4: Sample HPB CD3-F CD4, CD45, CD8

GATE: (P1 in all)

Plot 5: Sample HPB CD3-F CD4, CD45, CD8

GATE: (P1 in all)

Plot 6: Sample HPB CD3-F CD4, CD45, CD8

GATE: (R1 in (P1 in all))

Plot	Count	Volume (μ L)	% of This Plot	% of All	Mean CD3 FITC-A	Mean CD4 PE-A	CV CD3 FITC-A	CV CD4 PE-A
Plot 2: Sample HPB CD3-F CD4, CD45, CD8								
Gated on (R1 in (P1 in all))								
This Plot	99,183	0.0	100.0%	43.3%	23,875.3	1,762.1	62.0%	
Q3-UL	147	0.0	0.1%	0.1%	960.4	1,607.0	53.3%	
Q3-UR	59,194	0.0	59.7%	25.8%	32,131.1	2,933.7	28.5%	
Q3-LL	21,455	0.0	21.6%	9.4%	547.0	21.5	47.8%	
Q3-LR	18,387	0.0	18.5%	8.0%	24,701.0	22.5	29.2%	
Plot 3: Sample HPB CD3-F CD4, CD45, CD8								
Gated on (P1 in all)								
This Plot	102,719	0.0	100.0%	44.8%	23,071.3	38,996.7	65.7%	
R1	99,183	0.0	96.6%	43.3%	23,875.3	40,270.2	62.0%	
Q1-UL	21,694	0.0	21.1%	9.5%	540.5	32,690.2	45.5%	
Q1-UR	77,625	0.0	75.6%	33.9%	30,356.9	42,417.0	30.7%	
Q1-LL	3,380	0.0	3.3%	1.5%	447.0	1,141.0	76.5%	
Q1-LR	20	0.0	0.0%	0.0%	6,449.7	2,092.6	97.0%	

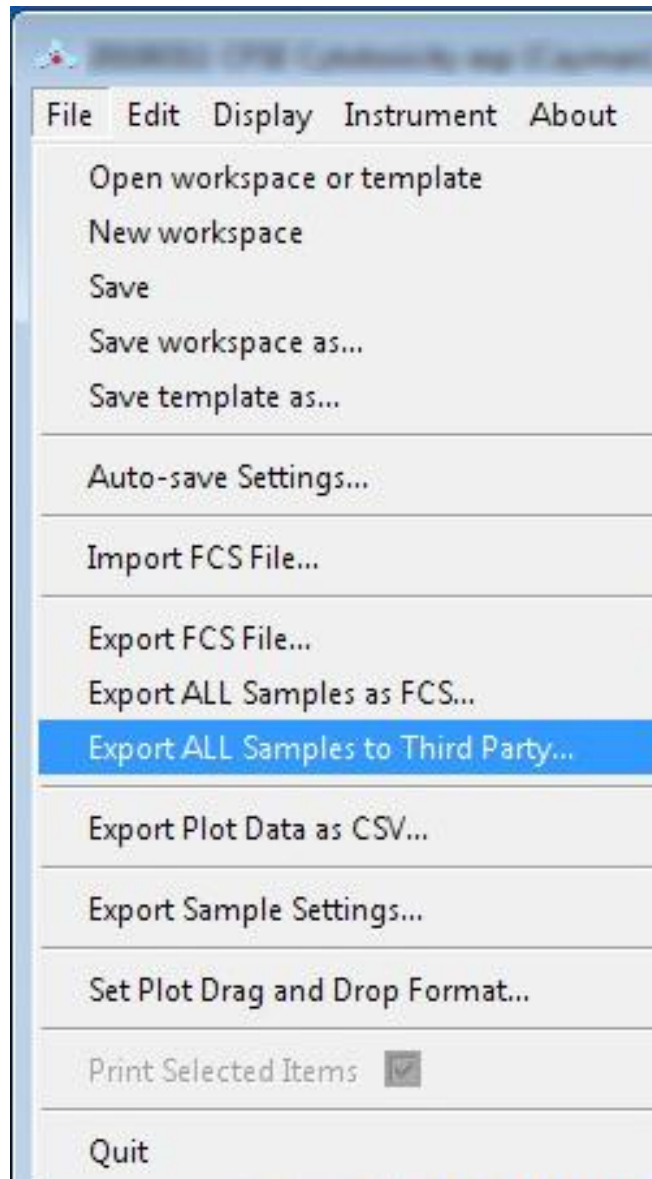
Histogram, Dot Plot, and Density Plot Display Area

Analysis and Gating Tools

Plot Statistics



Export Data to Third Party



Resources: Technical Bulletins

- Display of BD Accuri™ C6 Software-Generated FCS 3.0 Files Using FlowJo™ Version 8.8.7 or 9.4.11 for the Mac®
- Display of BD Accuri™ C6 Software-Generated FCS 3.0 Files Using FlowJo™ for PC and Mac® v7.6

www.bdbiosciences.com/support/resources/accuri



Acknowledgments

Stacey Roys

For Research Use Only. Not for use in diagnostic or therapeutic procedures.

FlowJo is a trademark of Tree Star, Inc.

Class 1 Laser Product.

BD, BD Logo and all other trademarks are property of Becton, Dickinson and Company. © 2013 BD

