
Simplify Immunophenotyping Using Kits and Templates on the BD Accuri™ C6

Presented by:

Stacey Roys, Scientist, BD Biosciences

Lora Tirri, BD Accuri Product Manager, BD Biosciences



Helping all people
live healthy lives

Outline

- Background: Immunophenotyping
- Introduction to the BD Accuri™ C6 Flow Cytometer
- BD Kits and Templates
- BD Templates on the Web
- How to Create Your Own Templates
- Coming Soon....



Helping all people
live healthy lives

Outline

- **Background: Immunophenotyping**
- Introduction to the BD Accuri™ C6 flow cytometer
- BD Kits and Templates
- BD Templates on the Web
- How to Create Your Own Templates
- Coming Soon....



Helping all people
live healthy lives

CD4⁺ T-Cell Subsets

- **Th1**: Activate infected macrophages and help B cells; immunity to viruses, intracellular bacteria, and parasites
- **Th2**: Help B cells and switch antibody production; immunity to extracellular parasites
- **Th17**: Inflammation and autoimmunity; immunity to extracellular bacteria and fungi
- **Regulatory T cells (Tregs)**: Regulate the immune response (inhibition or downregulation)



Helping all people
live healthy lives

CD4 T-Cell Phenotypes/Subsetting

- Different CD4 T-cell phenotypes can be defined by cell surface markers, expression of transcription factors, and cytokine production.
- Many CD4 subsets cannot be detected using cell surface markers alone.
 - For example, identification of Th17 cells requires the detection of the IL-17 cytokine or ROR γ T transcription factor.
 - Detection of intracellular markers by flow cytometry requires special consideration.
 - Flow cytometry allows the study of multiple subsets within the same sample simultaneously.



Helping all people
live healthy lives

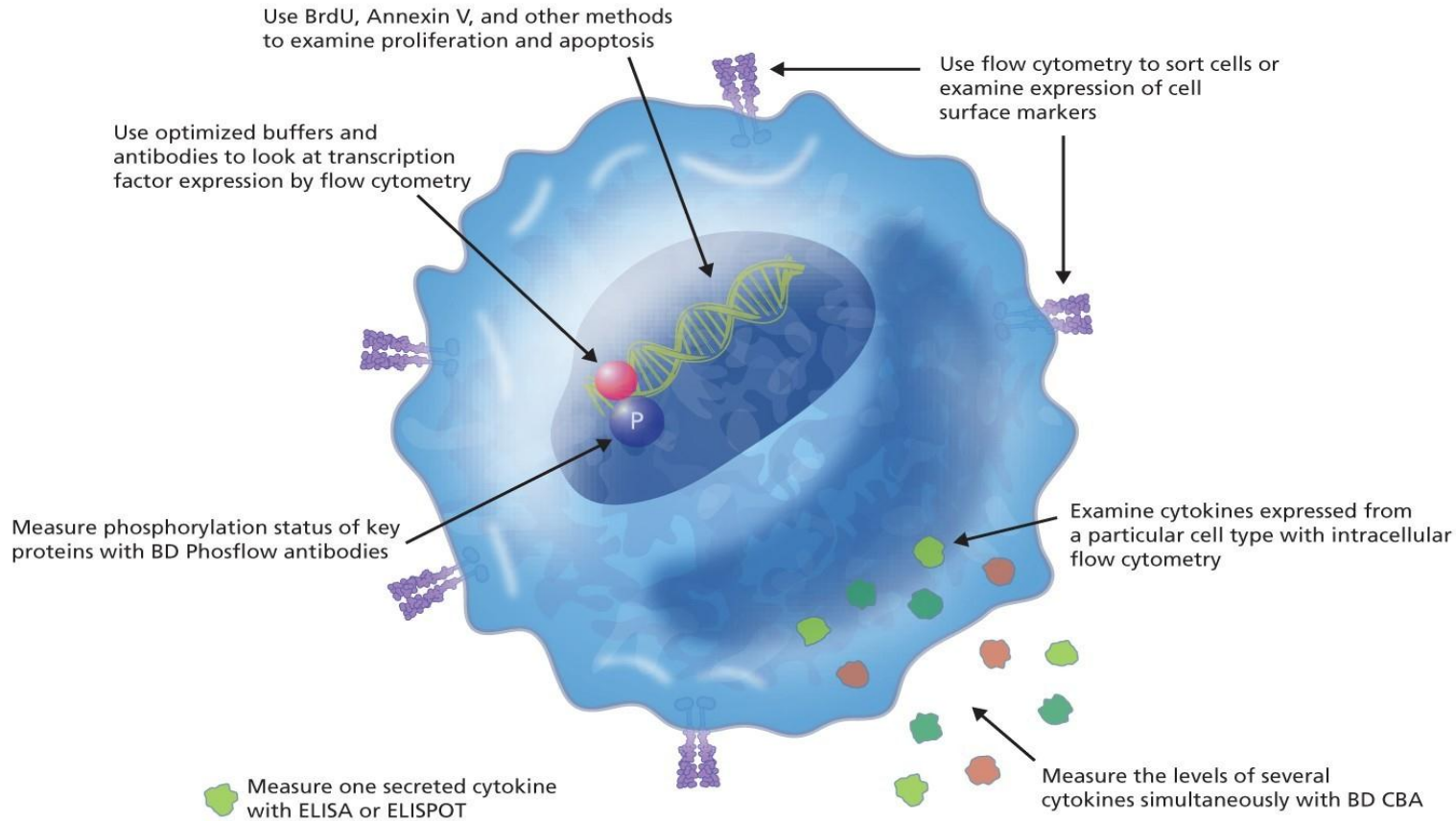
Key T-Cell Surface Markers and Cytokines

Marker	Significance
CD3	T-cell lineage marker, all T cells are CD3+
CD4	Marker for T helper cells
CD8	Marker for cytotoxic T cells
IFN- γ	Th1 signature cytokine
IL-4	Th2 signature cytokine
IL-17A	Th17 signature cytokine
CD45RA	Marker for naïve T cells
CD45RO	Marker for memory T cells
CD62L	Marker for memory cell activation
CD197	Marker for central/effector memory cells



Helping all people
live healthy lives

Flow Cytometry Allows the Detection of Multiple Parameters



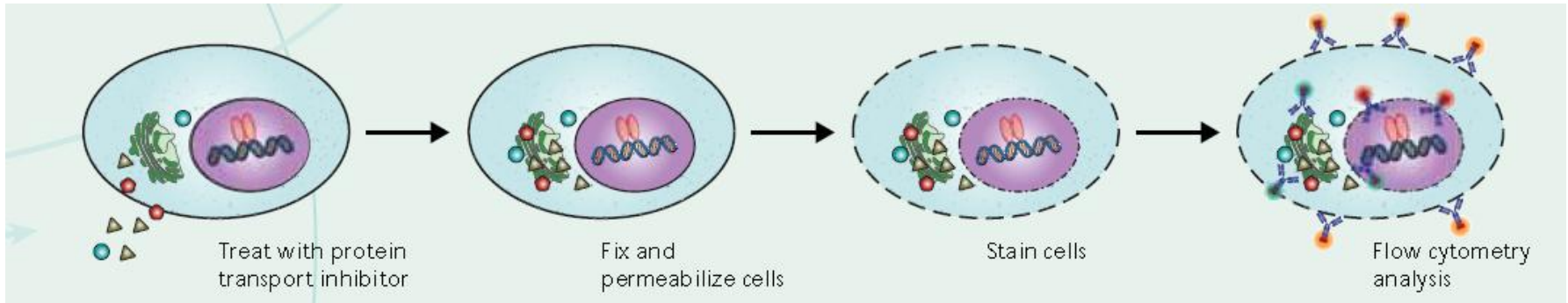
Methods to Detect Cytokines

Capability	BD CBA	ELISA	Intracellular Flow
Allows the detection of multiple cytokines in the same experiment			
Can obtain the phenotype of specific cells expressing the cytokine(s) of interest			
Can measure the quantity of the cytokine secreted			



Helping all people
live healthy lives

Detection of Cytokines by Intracellular Flow



- Cytokines are secreted proteins that need to be trapped inside the cell, fixed, and permeabilized for detection by intracellular flow.
- Cytokines can be trapped inside the cell using a protein transport inhibitor such as BD GolgiStop™ (Monensin) or BD GolgiPlug™ (Brefeldin A).
- BD kits contain the appropriate protein transport inhibitor, buffers, cocktails, and detailed protocols to simplify experiments and reduce variability across experiments and laboratories.



Helping all people
live healthy lives

Outline

- Background: Immunophenotyping
- **Introduction to the BD Accuri™ C6 flow cytometer**
- BD kits and templates
- BD templates on the web
- How to create your own templates
- Coming soon....



Helping all people
live healthy lives

Introduction

The BD Accuri™ C6 Flow Cytometer System



An Affordable, Full-Featured, Easy-to-Use Flow Cytometer

Two lasers and six detectors



Helping all people
live healthy lives

Intuitive Software

Sample Grid

Cytometer Status

Fluidics Controls

Run Criteria

Real-Time Updates

The screenshot displays the BD CSampler software interface with the following components:

- Menu Bar:** File, Edit, Display, Instrument, About
- Navigation Tabs:** Manual Collect, Auto Collect, Analyze, Statistics, Batch Analysis
- Control Panel (Left):**
 - Plate Type: 24 tube rack
 - Eject Plate button
 - Plate Name: Click here to name plate
 - Panel 3 grid (A1-H12) with cell C1 highlighted.
 - Run Settings: Run with Limits, 10000 events in Lymphs.
 - Fluidics: Flow Rate 60 µL/min, Core Size 22 µm.
 - Threshold: Set Threshold, 250,000 on FSC-H.
 - Buttons: Backflush, Unclog, Wash, Agitate.
 - ADD to C01 button.
 - Last Run Summary: 0 Events, 170.136 Time, 4:31.7.
- Plots (Center):**
 - Plot 1: C01 Panel 3 (GATE) [No Gating] - Scatter plot of SSC-A vs FSC-A with a red gate for Lymphs (38.0%).
 - Plot 2: C01 Panel 3 (GATE) (Lymphs in all) - Scatter plot of FL2-A vs FL1-A with gates P2 (9.1%) and P3 (80.2%).
 - Plot 3: C01 Panel 3 (GATE) (Lymphs in all) - Histogram of FL4-A.
 - Plot 4: C01 Panel 3 (GATE) (Lymphs in all) - Scatter plot of FL1-A vs FL2-A.
 - Plot 5: Placeholder for a new plot.
 - Plot 6: Placeholder for a new plot.
 - Plot 7: C01 Panel 3 (GATE) (Lymphs in all) - Placeholder for a new plot.
- Statistics Tables (Bottom):**

Plot	Count	Volume (µL)	% of This Plot	% of All	Mean FL1-A	Mean FL2-A	CV FL1-A	CV FL2-A	Median FL1-A	Median FL2-A
Plot 2: C01 Panel 3 (Gated on (Lymphs in all))										
This Plot	64,614	297	100.00%	37.98%	33,627.31	4,263.52	77.45%	137.12%		
P2	5,896	297	9.12%	3.47%	2,400.14	15,828.54	28.98%	35.55%		
P3	51,839	297	80.23%	30.47%	41,321.67	3,420.89	32.62%	31.27%		
Plot 6: C01 Panel 3 (Gated on (Lymphs in all))										
This Plot	64,614	297	100.00%	37.98%	17,542.71	202.67%				
Plot 7: C01 Panel 3 (Gated on (Lymphs in all))										
This Plot	64,614	297	100.00%	37.98%	4,263.52	33,627.31	137.12%	77.45%		

Histograms

Dot Plots

Density Plots

Analysis and Gating Tools

Plot Statistics



Helping all people live healthy lives

Outline

- Background: Immunophenotyping
- Introduction to the BD Accuri™ C6 Flow Cytometer
- **BD Kits and Templates**
- BD Templates on the Web
- How to Create Your Own Templates
- Coming Soon....



Helping all people
live healthy lives

BD Kits and Templates

The BD Accuri™ C6 Personal Flow Cytometer
Conduct powerful assays in just a few easy steps.

The image illustrates the workflow for using the BD Accuri C6 Personal Flow Cytometer. It is divided into three numbered steps: 1. The cytometer unit, a compact white and red device. 2. Reagents and kits, including a blue box, a clear bottle, and two vials. 3. Data analysis, shown on a tablet displaying two flow cytometry plots. The first plot, titled 'Plot 1: HE2 (GATE) [No Gating]', shows a single population of cells at 100.0%. The second plot, titled 'Plot 1: All Concanavalin FITC + PI (GATE) [No Gating]', shows a population of cells at 91.5%. The background is a dark red with a grid pattern and red arrows indicating the flow from step 1 to 2, and then to 3.

Arguably, as easy as cell analysis is going to get.



Helping all people
live healthy lives

BD Immunology Kits

Application	Kit	Cat. No.
Naïve/Memory T Cell	CD45RA/CD45RO/ CD3/CD4	<u>340571</u>
	CD45RA/CD62L/CD3/CD4	<u>340977</u>
	Human Naïve/Memory T Cell Panel	<u>561438</u>
Intracellular T Cell	Th1/Th2/Th17 Phenotyping Kit	<u>560751</u>
	IFN- γ /CD69/CD8/CD3 Kit	<u>346048</u>
	IFN- γ /IL4 Kit	<u>340456</u>



Helping all people
live healthy lives

What is a Template?

The screenshot displays the BD Accuri C6 software interface. On the left, there is a 'Collect' panel with a 12x8 grid of sample wells (A1-H12) and a 'Run Settings' section. The 'Run Settings' section includes options for 'Run Unlimited' or 'Run with Limits', 'Fluidics' settings (Slow, Medium, Fast), and a 'Threshold' section. A large 'RUN' button is visible. The main area shows three plots: Plot 1: C09 (GATE) [No Gating], Plot 2: C09 (GATE) [No Gating], and Plot 3: C09 (GATE) (CD4+ in all). Below the plots, there are three prompts: 'Select plot type to make a new plot.' The bottom of the interface features three data tables for the plots.

Plot 1: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FSC-A	Mean SSC-A	CV FSC-A	CV SSC-A	Median FSC-A	Median SSC-A
All	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		
Lymphocytes	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		

Plot 2: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FL3 CD4 PerCP Cy5.5-A	Mean SSC-A	CV FL3 CD4 PerCP Cy5.5-A	CV SSC-A	Median FSC-A	Median SSC-A
All	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		
CD4+	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		

Plot 3: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FL1 CD45RA FITC-A	Mean FL4 CD197 AF 647-A	CV FL1 CD45RA FITC-A	CV FL4 CD197 AF 647-A	Median FSC-A	Median SSC-A
This Plot	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		
Central Memory	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		

A BD Accuri template is a predefined workspace that includes gates, labels, run criteria, and compensation settings for a specific assay.



Helping all people live healthy lives

Predefined Gating and Run Criteria for Ease of Use

BD Accuri™ C6

Collect
Analyze
Statistics
Batch Analysis

C09

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

C6 is connected and ready.

Plot 1: C09
GATE: [No Gating]

Plot 4: C09
GATE: [No Gating]

Plot 3: C09
GATE: [CD4+ in all]

Select plot type to make a new plot.

Run Settings

Run Unlimited

Run with Limits

Events: 0

In Ungated Sample

Min: 0 Sec

Do not collect events outside: Lymphocytes

Backflush Unclog

Fluidics

Slow Medium Fast

Flow Rate: 30 µL/min

Core Size: 10 µm

Custom

Flow Rate: 14 µL/min

Set Core Size

Core Size: 10 µm

Threshold

Set Threshold

80,000 on FSC-H

RUN

Set Color Compensation

Last Run: 0 Events, 0.00 Time, 0 Microliters, 0 Events / Sec, 0 Events / µL

Cumulative: All, Outside Lymphocytes

Delete Events: show warning

Data Capacity Used: 0% of 99,000,000 Events

Plot 1: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FSC-A	Mean SSC-A	CV FSC-A	CV SSC-A	Median FSC-A	Median SSC-A
All	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		
Lymphocytes	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		

Plot 4: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FL3 CD4 PerCP Cy5.5-A	Mean SSC-A	CV FL3 CD4 PerCP Cy5.5-A	CV SSC-A
All	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%
CD4+	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%

Plot 3: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FL1 CD45RA FITC-A	Mean FL4 CD197 AF 647-A	CV FL1 CD45RA FITC-A
Gated on (CD4+ in all)							
This Plot	0	0	100.00%	100.00%	0.00	0.00	0.00%
Central Memory	0	0	100.00%	100.00%	0.00	0.00	0.00%



Helping all people live healthy lives

Compensation

BD Accuri™ C6

Collect Analyze Statistics Batch Analysis

H12

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 Settings

Run Unlimited Run with Limits

Fluidics

Slow Medium Fast Custom

Flow Rate 86 µL/min
Core Size 22 µm

Flow Rate 14 µL/min
Set Core Size
Core Size 10 µm

Threshold

80,000 on FSC-H
None

Set Threshold

Backflush Unclog

RUN

Set Color Compensation

Last Run Cumulative

Events 0
Time 0:00:0
Microliters 0
Events / Sec 0
Events / µL 0

Delete Events show warning

All Outside Lymphs

Data Capacity Used
0% of 96,000,000 Events

Plot 1: A01 GATE [No Gating]

Plot 2: A01 GATE [Lymphs in all]

Plot 3: A01 GATE [CD4+ in (Lymphs in all...)]

Compensation Settings for A01

Correct FL1 by subtracting a percentage of:

FL2: 4.68 % FL3: 0.00 % FL4: 0.00 %

Correct FL2 by subtracting a percentage of:

FL1: 0.10 % FL3: 0.22 % FL4: 0.00 %

Correct FL3 by subtracting a percentage of:

FL1: 0.00 % FL2: 21.00 % FL4: 0.00 %

Correct FL4 by subtracting a percentage of:

FL1: 0.00 % FL2: 0.00 % FL3: 9.20 %

Apply to:

Reset all to 0.00% A01 All samples

Preview Apply & Close Cancel & Close

Select plot type to make a new plot.

	CV FSC-A	CV SSC-A	Median FSC-A	Median SSC-A
All	0	0	100.00%	100.00%
Lymphs	0	0	100.00%	100.00%

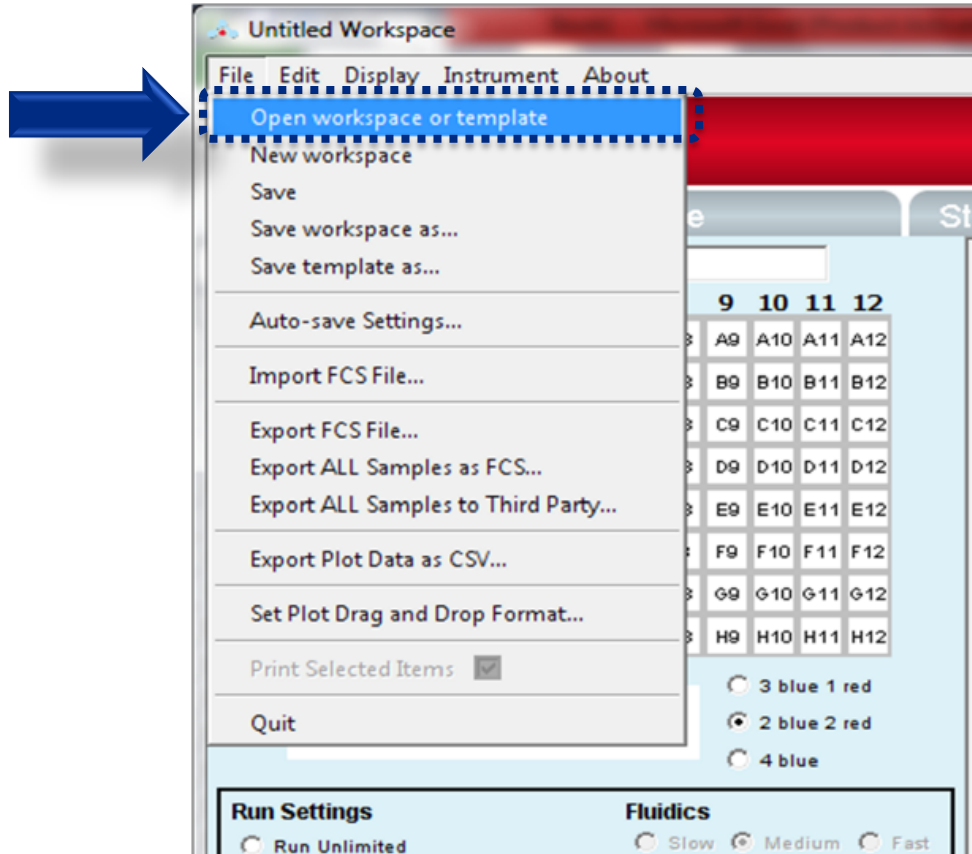
Plot 2: A01	Count	Volume (µL)	% of This Plot	% of All	Mean FL3 CD4 PerCP-Cy5.5-A	CV FL3 CD4 PerCP-Cy5.5-A	Median FL3 CD4
Gated on (Lymphs in all)							
This Plot	0	0	100.00%	100.00%	0.00	0.00%	0.00%
CD4+ (1,983.0 / 15,40...)	0	0	100.00%	100.00%	0.00	0.00%	0.00%

Plot 3: A01	Count	Volume (µL)	% of This Plot	% of All	Mean FL4 IL-4 APC-A	Mean FL2 IL-17A PE-A	CV FL4 IL-4 APC-A
Gated on (CD4+ in (Lymphs in all))							
This Plot	0	0	100.00%	100.00%	0.00	0.00	0.00%
Q1-UL	0	0	100.00%	100.00%	0.00	0.00	0.00%



Helping all people live healthy lives

How to Open a Template



Helping all people
live healthy lives

How to Open a Template

The screenshot displays the BD Accuri C6 software interface. The main window is titled 'Untitled Workspace' and has a menu bar with 'File', 'Edit', 'Display', 'Instrument', and 'About'. Below the menu bar is a red header with 'BD Accuri™ C6'. The interface is divided into several sections: 'Collect', 'Analyze', 'Statistics', and 'Batch Analysis'. The 'Collect' section shows a grid of sample wells (A1-A12, B1-B12, etc.) and a 'C6 Cytometer not connected.' message. The 'Analyze' section shows a plot titled 'Plot 1: A01 (GATE) [No Gating]'. The 'Statistics' section shows a table with columns for 'Count', 'Volume (µL)', '% of This Plot', '% of All', 'Mean FSC-A', 'Mean SSC-A', 'CV FSC-A', and 'CV SSC-A'. The 'Batch Analysis' section is currently empty.

An 'Open' dialog box is open, showing a list of templates. The 'Look In' field is set to 'Templates'. The list includes:

- Active caspase-3 PE_550914 template_24 tube rack
- Annexin V FITC_556570 template_24 tube rack
- Annexin V PE_559763 template
- Bacterial Viability_BD Cell Viab kit
- BD FastImmune_340456 template_24 tube rack
- BD FastImmune_346048 template

The 'File Name' field is set to 'Active caspase-3 PE_550914 template_24 tube rack.c6t'. The 'Files of Type' dropdown is set to 'Workspace and Template (*.c6, *.c6t, *.c6f, *.c6i)'. The 'Open' button is highlighted.

A blue arrow points from the 'Run Settings' section to the 'Open' dialog box. A dashed blue circle highlights the 'Open' dialog box.

Plot 1: A01	Count	Volume (µL)	% of This Plot	% of All	Mean FSC-A	Mean SSC-A	CV FSC-A	CV SSC-A
All	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%



Helping all people
live healthy lives

How to Use a Template: Example Using Human Naïve/Memory T Cell Panel (Cat. No. 561438)

BD Accuri™ C6

Collect Analyze Statistics Batch Analysis

C09

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 Settings

Run Unlimited

Run with Limits

0 events in Ungated Sample

0 Min 30 Sec

0 µL

Do not collect events outside Lymphocytes

Backflush Unclog

Fluidics

Slow Medium Fast

Flow Rate 35 µL/min
Core Size 16 µm

Custom

Flow Rate 14 µL/min
Set Core Size
Core Size 10 µm

Threshold

Set Threshold
80,000 on FSC-H
None

RUN

Set Color Compensation

Last Run

0 Events
0:00.0 Time
0 Microliters
0 Events / Sec
0 Events / µL

Cumulative

0
0:00.0
0
0
0

Delete Events show warning

All
 Outside Lymphocytes

Data Capacity Used
0% of 98,000,000 Events

Plot 1: C09 (GATE) (No Gating)

Select plot type to make a new plot.

Plot 4: C09 (GATE) (No Gating)

Select plot type to make a new plot.

Plot 3: C09 (GATE) (CD4+ in all)

Select plot type to make a new plot.

Plot 1: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FSC-A	Mean SSC-A	CV FSC-A	CV SSC-A	Median FSC-A	Median SSC-A
All	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		
Lymphocytes	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%		

Plot 4: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FL3 CD4 PerCP Cy5.5-A	Mean SSC-A	CV FL3 CD4 PerCP Cy5.5-A	CV SSC-A
All	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%
CD4+	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%

Plot 3: C09	Count	Volume (µL)	% of This Plot	% of All	Mean FL1 CD45RA FITC-A	Mean FL4 CD197 AF 647-A	CV FL1 CD45RA FITC-A	CV FL4 CD197 AF 647-A	CV FL1 CD45RA FITC-A
This Plot	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%	0.00%
Central Memory	0	0	100.00%	100.00%	0.00	0.00	0.00%	0.00%	0.00%

Panel:
CD4 PerCP-Cy™ 5.5
CD45RA FITC
CD197 Alexa
Fluor® 647



Helping all people
live healthy lives

How to Use a Template: Verify Settings, Collect, Adjust as Necessary

BD Accuri™ C6

Collect Analyze Statistics Batch Analysis

B01 Naive Memory panel 561438

	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 Settings: Run with Limits, 0 events, Ungated Sample, 0 Min 30 Sec, Do not collect events outside Lymphocytes, Backflush, Unclog.

Fluidics: Slow Medium Fast, Flow Rate 06 µL/min, Core Size 22 µm, Custom, Flow Rate 14 µL/min, Core Size 10 µm.

Threshold: Set Threshold, 80,000 on FSC-H, None.

ADD to B01

Set Color Compensation

Last Run: 0 Events, 0.00.0 Time, 2:01.5, 0 Microliters, 133 Events / Sec, 743 Events / µL, 680. Delete Events, show warning, All, Outside Lymphocytes, Data Capacity Used <1% of 96,000,000 Events.

Plot 1: B01 Naive Memory panel 561...
GATE [No Gating]

SSC-A vs FSC-A plot showing Lymphocytes gate at 21.3%.

Plot 4: B01 Naive Memory panel 561...
GATE [No Gating]

SSC-A vs FL3 CD4 PerCP Cy5.5-A plot showing CD4+ gate at 5.0%.

Plot 3: B01 Naive Memory panel 561...
GATE [CD4+ in all]

FL1 CD45RA FITC-A vs FL4 CD197 AF 647-A plot showing Central Memory (25.7%), Naive (37.9%), Effector Memory (36.0%), and Q1-LR (0.4%) populations.

Select plot type to make a new plot.

Plot 4: B01 Naive Memory panel 561438	Count	Volume (µL)	% of This Plot	% of All	Mean FL3 CD4 PerCP Cy5.5-A	Mean SSC-A	CV FL3 CD4 P
All	90,411	133	100.00%	100.00%	2,241.89	420,307.18	
CD4+	4,514	133	4.99%	4.99%	18,112.84	87,837.91	

Plot 3: B01 Naive Memory panel 561438 Gated on (CD4+ in all)	Count	Volume (µL)	% of This Plot	% of All	Mean FL1 CD45RA FITC-A	Mean FL4 CD197 AF 647-A	CV I
This Plot	4,514	133	100.00%	4.99%	8,411.61	5,388.40	
Central Memory	1,161	133	25.72%	1.28%	1,191.51	5,723.84	
Naive	1,710	133	37.88%	1.89%	20,352.13	8,850.47	
Effector Memory	1,624	133	35.98%	1.80%	924.83	1,541.82	
Q1-LR	19	133	0.42%	0.02%	14,874.37	2,087.11	



Helping all people live healthy lives

Ready to Acquire Data

BD Accuri™ C6

Collect Analyze Statistics Batch Analysis

B01 Naive Memory panel 561438

	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 Settings

Run Unlimited
 Run with Limits

2000 events
in CD4+

Do not collect events outside
Lymphocytes

Backflush Unclog

Fluidics

Slow Medium Fast

Flow Rate 86 µL/min
Core Size 22 µm

Custom

Flow Rate 14 µL/min
Core Size 10 µm

Set Core Size

Threshold

Set Threshold

350,000 on FSC-H
None

ADD to B01

Set Color Compensation

Last Run

0	Events	90,411
0:00.0	Time	2:01.5
0	Microliters	133
0	Events / Sec	7.43
0	Events / µL	680

Cumulative

Delete Events show warning

All
 Outside Lymphocytes

Data Capacity Used
<1% of 96,000,000 Events

Select plot type to make a new plot.

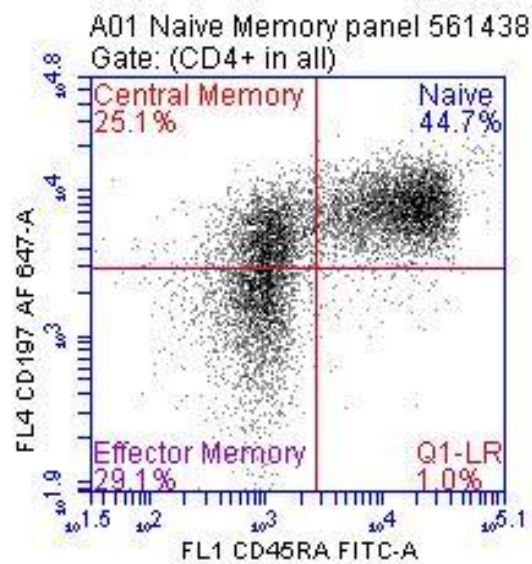
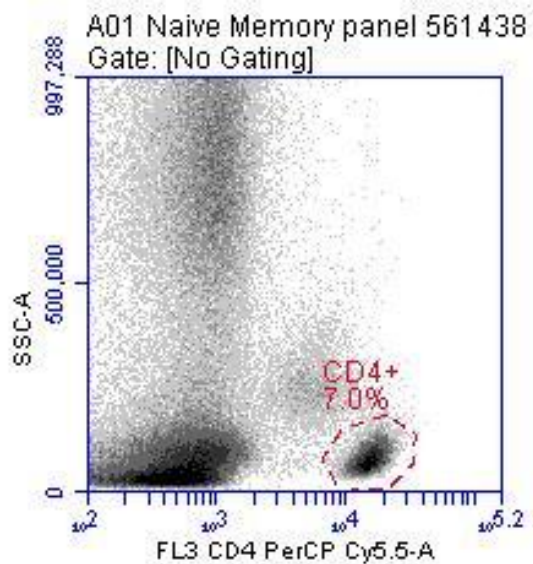
Select plot type to make a new plot.

Select plot type to make a new plot.



Helping all people
live healthy lives

Naïve/Memory T-Cell Data



Helping all people
live healthy lives

How to Further Optimize a Template: BD Multitest CD45RA/CD45RO/CD3/CD4 (340571)

The screenshot displays the BD Accuri C6 software interface. At the top, there is a menu bar (File, Edit, Display, Instrument, About) and the software name 'BD Accuri™ C6'. Below this are tabs for 'Collect', 'Analyze', 'Statistics', and 'Batch Analysis'. The 'Analyze' tab is active, showing a grid of 12 sample wells (A1-H12) with 'D6' selected in well D6. A status bar indicates 'C6 is connected and ready.' Below the grid are 'Run Settings' and 'Fluidics' sections. The 'Run Settings' section includes options for 'Run Unlimited' or 'Run with Limits', 'Flow Rate' (66 µL/min), 'Core Size' (22 µm), and 'Threshold' (Set Threshold). The 'Fluidics' section includes options for 'Slow', 'Medium', or 'Fast' flow rate, 'Flow Rate' (14 µL/min), and 'Core Size' (10 µm). The 'Run' button is highlighted. Below the 'Run' button is a 'Set Color Compensation' button. The main area shows four flow cytometry plots: Plot 1 (SSC-A vs FSC-A) with 'Lymphocytes 6.5%' gate; Plot 2 (SSC-A vs FL3 CD3 PerCP-A) with 'CD3+ 3.7%' gate; Plot 5 (FL4 CD4 APC-A vs FL1 CD45RA FITC-A) with 'Naive 11.5%' gate; and Plot 6 (FL4 CD4 APC-A vs FL2 CD45RO PE-A) with 'Memory 49.1%' gate. Below the plots are two instructions: 'Select plot type to make a new plot.' and 'Select plot type to make a new plot.' with icons for histogram, scatter plot, and contour plot. At the bottom, there are three data tables for the plots.

Plot 1: D06 Multitest 340571 4 color	Count	Volume (µL)	% of This Plot	% of All	Mean FSC-A	Mean SSC-A	CV FSC-A	CV SSC-A	Median F
All	130,063	87	100.00%	100.00%	608,351.90	110,126.28	117.06%	214.07%	
Lymphocytes	8,517	87	6.55%	6.55%	1,013,316.42	84,868.57	23.65%	39.44%	


Plot 2: D06 Multitest 340571 4 color	Count	Volume (µL)	% of This Plot	% of All	Mean FL3 CD3 PerCP-A	Mean SSC-A	CV FL3 CD3 PerCP-A
All	130,063	87	100.00%	100.00%	982.33	110,126.28	279.65%
CD3+	4,787	87	3.68%	3.68%	11,875.01	82,745.39	34.32%

Plot 5: D06 Multitest 340571 4 color	Count	Volume (µL)	% of This Plot	% of All	Mean FL1 CD45RA FITC-A	Mean FL4 CD4 APC-A	CV FL1 CD45
This Plot	4,787	87	100.00%	3.68%	2,554.49	20,614.70	
Naive	550	87	11.49%	0.42%	5,731.86	29,650.44	

Panel:
CD3 PerCP
CD4 APC
CD45RA FITC
CD45RO PE

How to Optimize a Template: Zoom Tool

File Edit Display Instrument About

BD CSampler™ 

Manual Collect Auto Collect Analyze Statistics Batch Analysis

Plate Type: 24 tube rack

Plate Name:

A01 Multitest 340571 4 color

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

Run Unlimited Run with Limits

0 events in Ungated Sample Min 30 Sec

Do not collect events outside Lymphocytes

Backflush Unclog Wash Agitate

ADD to A01

Set Color Compensation

Last Run Cumulative Delete Events show warning

0 Events 130,063
0:00:00 Time 1:19.2
0 Microliters 87
0 Events / Sec 1,641
0 Events / μ L 1,495

All Outside Lymphocytes

Data Capacity Used 1% of 24,000,000 Events

Plot 1: A01 Multitest 340571 4 color (GATE) [No Gating]

Plot 2: A01 Multitest 340571 4 color (GATE) [No Gating]

Plot 5: A01 Multitest 340571 4 color (GATE) (CD3+ in all)

Plot 6: A01 Multitest 340571 4 color (GATE) (CD3+ in all)

Plot 7: A01 Multitest 340571 4 color (GATE) [No Gating]

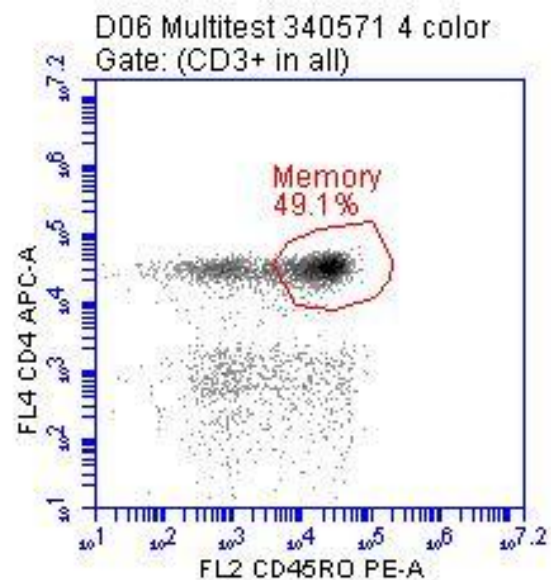
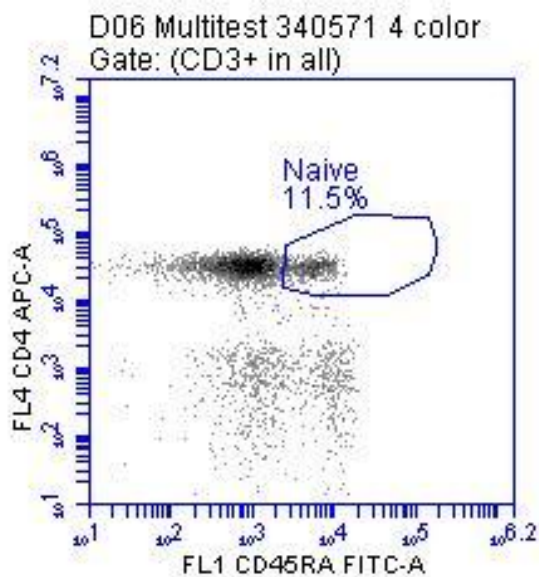
Select plot type to make a new plot.

Plot 3: A01 Multitest 340571 4 color	Count	Volume (μ L)	% of This Plot	% of All	Mean FL3 CD3 PerCP-A	Mean FL4 CD4 APC-A	CV FL3 CD3 PerCP-A
All	130,063	87	100.00%	100.00%	982.33	1,172.21	27.1
CD3+CD4+	3,495	87	2.69%	2.69%	12,114.68	29,970.84	10.1



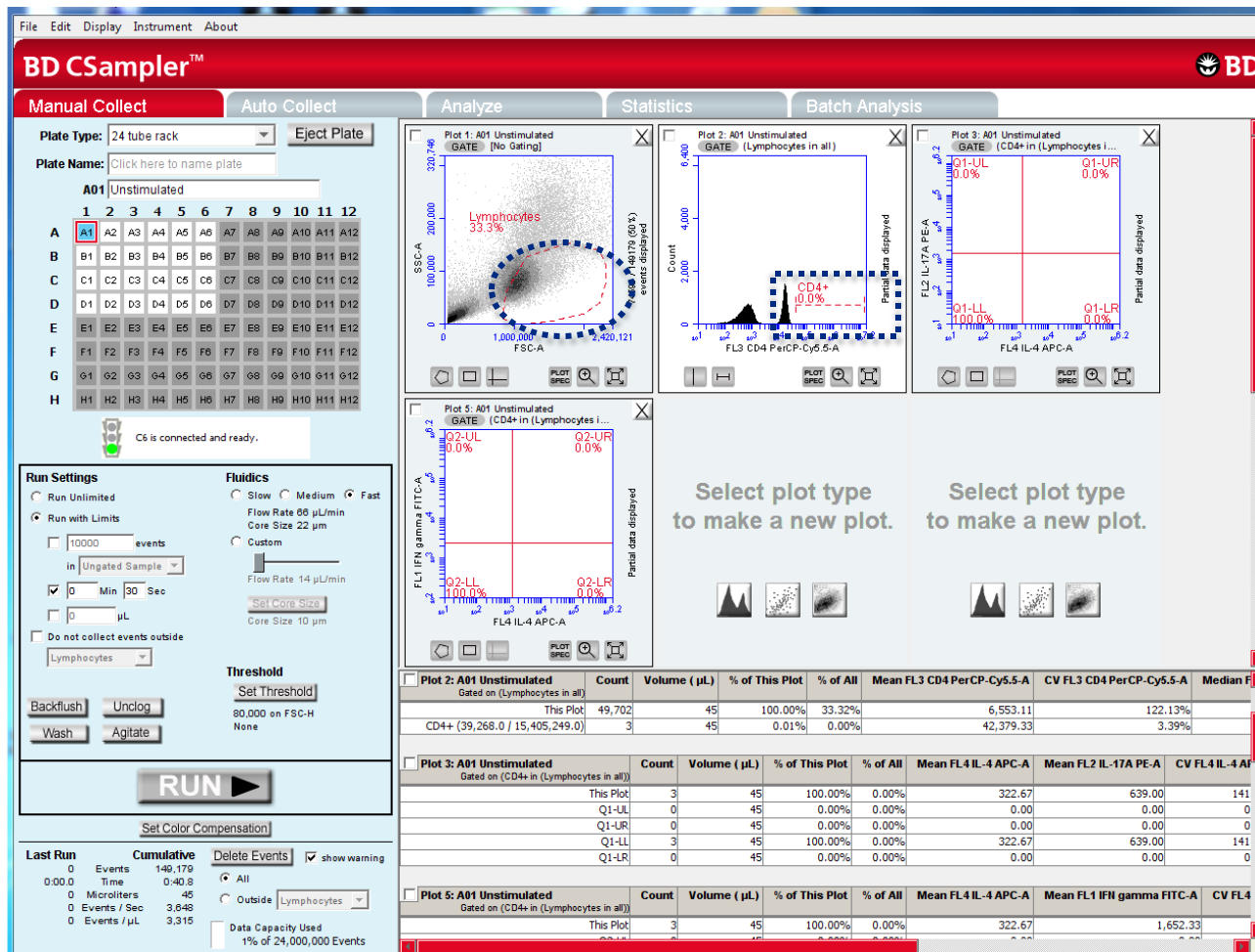
Helping all people live healthy lives

BD Multitest CD45RA/CD45RO/CD3/CD4 Data



Helping all people
live healthy lives

How to Optimize a Template for the BD CSampler™ Accessory: Th1/Th2/Th17 Phenotyping Kit (560751)



Panel:
CD4 PerCP-Cy5.5
IL-17A PE
IFN-γ FITC
IL-4 APC



Helping all people
live healthy lives

How to Optimize a Template: BD CSampler

File Edit Display Instrument About

BD CSampler™

Manual Collect Auto Collect Analyze Statistics Batch Analysis

Plate Type: 24 tube rack Eject Plate

Plate Name: Click here to name plate

A01 Unstimulated

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

C6 is connected and ready.

Run Settings

Run Unlimited Run with Limits

10000 events in Ungated Sample

0 Min 30 Sec

0 µL

Do not collect events outside Lymphocytes

Fluidics

Slow Medium Fast

Flow Rate 66 µL/min Core Size 22 µm

Flow Rate 14 µL/min

Set Core Size

Core Size 10 µm

Set Threshold

300,000 on FSC-H

None

Backflush Unclog Wash Agitate

RUN

Set Color Compensation

Last Run

0 Events 110.572 Cumulative
0.00 Time 0-40.8
0 Microliters 45
0 Events / Sec 2.704
0 Events / µL 2.457

Delete Events show warning

All Outside Lymphocytes

Data Capacity Used <1% of 24,000,000 Events

Plot 1: A01 Unstimulated GATE (No Gating)

Plot 2: A01 Unstimulated GATE (Lymphocytes in all)

Plot 3: A01 Unstimulated GATE (CD4+ in Lymphocytes i...)

Plot 5: A01 Unstimulated GATE (CD4+ in Lymphocytes i...)

Select plot type to make a new plot.

Select plot type to make a new plot.

Plot 2: A01 Unstimulated Gated on (Lymphocytes in all)	Count	Volume (µL)	% of This Plot	% of All	Mean FL3 CD4 PerCP-Cy5.5-A	CV FL3 CD4 PerCP-Cy5.5-A	Median F
This Plot	71,783	45	100.00%	64.92%	7,180.20	162.29%	
CD4+ (3,308.0 / 15,405,249.0)	29,667	45	41.33%	26.83%	16,580.21	80.40%	

Plot 3: A01 Unstimulated Gated on (CD4+ in Lymphocytes in all)	Count	Volume (µL)	% of This Plot	% of All	Mean FL4 IL-4 APC-A	Mean FL2 IL-17A PE-A	CV FL4 IL-4 AP
This Plot	29,667	45	100.00%	26.83%	582.95	360.59	13,796
Q1-UL	6	45	0.02%	0.01%	183.17	8,644.17	73
Q1-UR	1	45	0.00%	0.00%	1,646,264.00	1,606.00	0
Q1-LL	29,644	45	99.92%	26.81%	62.99	358.85	286
Q1-LR	16	45	0.05%	0.01%	861,232.50	401.56	386


Plot 5: A01 Unstimulated Gated on (CD4+ in Lymphocytes in all)	Count	Volume (µL)	% of This Plot	% of All	Mean FL4 IL-4 APC-A	Mean FL1 IFN gamma FITC-A	CV FL4
This Plot	29,667	45	100.00%	26.83%	582.95	932.94	



Helping all people live healthy lives

How to Optimize a Template: BD CSampler

File Edit Display Instrument About

BD CSampler™ 


Manual Collect **Auto Collect** Analyze Statistics Batch Analysis

Plate Type: 24 tube rack

Plate Name:

Ctrl-click to view sample settings. [Select All](#) [Deselect All](#)

A	1	2	3	4	5	6	7	8	9	10	11	12
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

 CS is connected and ready.

Run Limits

in Min S

Do not collect events outside less than

Wash Settings

Agitate Plate

None every Well(s) Run Horizontally

None every Min Run Vertically

Fluidics

Flow Rate: 14 µL/min
Pore Size: 10 µm

Set Threshold

Delete events on less than (Minimum=10)

Well	Sample Name	Rename FL1	Rename FL2	Rename FL3	Rename FL4	Notes
A01	Unstimulated	FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A02		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A03		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A04		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A05		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A06		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B01		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B02		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B03		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B04		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B05		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B06		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C01		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C02		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C03		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C04		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C05		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C06		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D01		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D02		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D03		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D04		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D05		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D06		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	



Helping all people live healthy lives

How to Optimize a Template: BD CSampler

File Edit Display Instrument About

BD CSampler™

Manual Collect **Auto Collect** Analyze Statistics Batch Analysis

Plate Type: 24 tube rack Eject Plate

Plate Name: [Click here to name plate](#)

Ctrl-click to view sample settings. Select All Deselect All

1 2 3 4 5 6 7 8 9 10 11 12

Well	Sample Name	Rename FL1	Rename FL2	Rename FL3	Rename FL4	Notes
A01	Unstimulated	FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A02	PMA+Ionomycin	FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A03	FITC single	FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A04	APC single	FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A05	PE single	FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
A06	PerCpCy5.5 single	FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B01		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B02		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B03		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B04		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B05		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
B06		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C01		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C02		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C03		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C04		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C05		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
C06		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D01		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D02		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D03		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D04		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D05		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	
D06		FL1 IFN gamma FITC	FL2 IL-17A PE	FL3 CD4 PerCP-Cy5.5	FL4 IL-4 APC	

C6 is connected and ready.

Run Limits

0000 events

in CD4+ in (Lymph...)

0 μL

Do not collect events outside Lymphocytes

Fluidics

Flow Rate: 86 μL/min
Core Size: 22 μm

Slow Medium Fast

Set Threshold

Delete events on (Minimum=10)

FSC-H less than 900000

less than 0

Wash Settings

None

Apply Settings Remove Settings

Agitate Plate

None every 1 Well(s) Run Horizontally

None every 1 Min Run Vertically

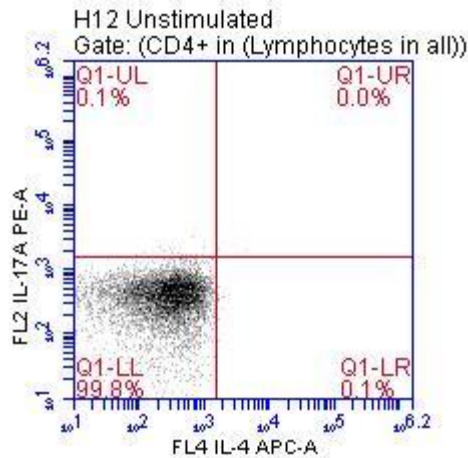
OPEN RUN DISPLAY



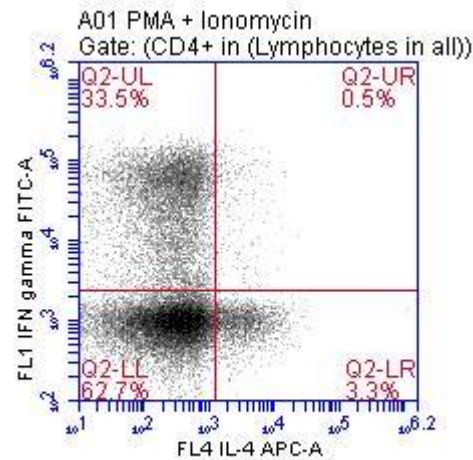
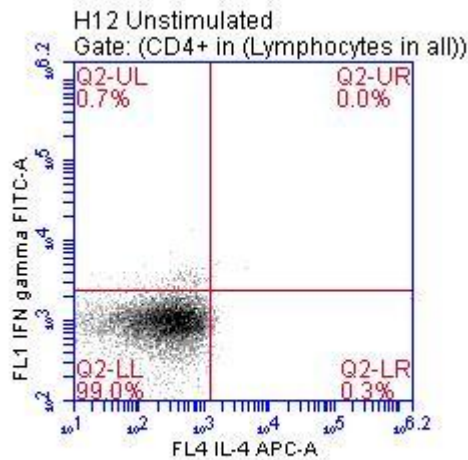
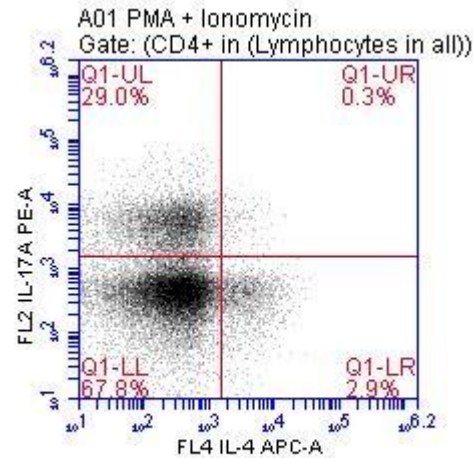
Helping all people
live healthy lives

BD Pharmingen™ Human Th1/Th2/Th17 Phenotyping Kit Data

Unstimulated



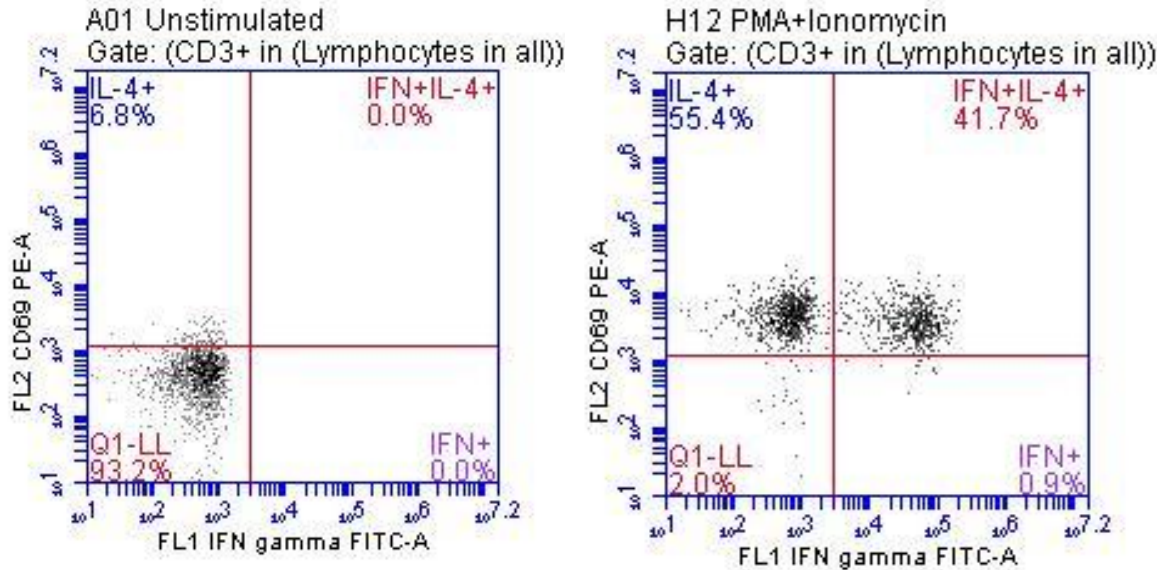
Stimulated



Helping all people
live healthy lives

Application-Specific Templates: BD FastImmune™ Kits (346048 and 340456)

- Detection of intracellular T-cell cytokines using preconfigured kits. Available with or without CD69 (activation marker).



Panel:
CD8 PerCP-Cy5.5
CD3 APC
IFN- γ FITC
CD69 PE



Helping all people
live healthy lives

Where to Find the Templates

- www.bdbiosciences.com/go/templates



Category	Product Information Sheet	Brand	Kit	Cat. No.	Template
Cell Biology	Apoptosis Kits and Templates	BD Pharmingen™	Annexin V FITC Apoptosis Detection Kit II	556570	Download
		BD Pharmingen™	Annexin V PE Apoptosis Detection Kit I	559763	Download
		BD™	MitoScreen (JC-1) Kit	551302	Download
		BD Pharmingen™	Caspase-3 PE Assay Kit	550914	Download
	Cell Cycle and DNA Kits and Templates	BD Cycletest™ Plus	DNA Reagent Kit	340242	Download
		BD Pharmingen™	FITC BrdU Flow Kit	559619	Download
		BD Pharmingen™	APC BrdU Flow Kit	552598	Download
Immunology	Naïve/Memory T- Cell Kits and Templates	BD Multitest™	CD45RA/CD45RO/CD3/CD4	340571	Download
		BD Multitest™	CD45RA/CD62L/CD3/CD4	340977	Download
		BD Pharmingen™	Human Naïve/Memory T Cell Panel	561438	Download
	Intracellular T-Cell Kits and Templates	BD Pharmingen™	Th1/Th2/Th17 Phenotyping Kit	560751	Download
		BD FastImmune™	IFN-γ/CD69/CD8/CD3 Kit	346048	Download
		BD FastImmune™	IFN-γ/IL4 Kit	340456	Download

BD Templates Available for Free on the Web

EXISTING

6, 8-Peak Bead
BD™ CBA
Water Quality

NEW

Proliferation: BrdU
Apoptosis: Annexin V
Apoptosis: Caspase-3
DNA Analysis
BD™ MitoScreen
Stem Cell Flow Kits
T-Cell Cytokines
Naïve/memory T Cells



Helping all people
live healthy lives

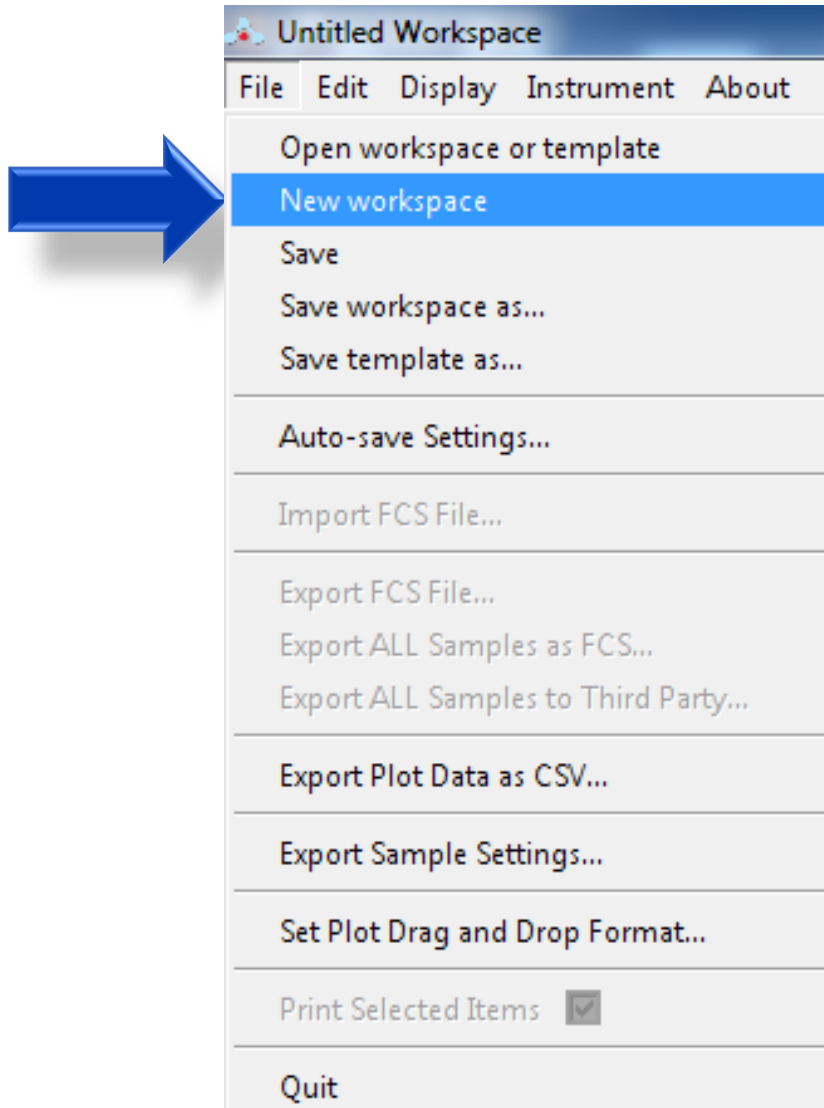
Outline

- Background: Immunophenotyping
- Introduction to the BD Accuri™ C6 Flow Cytometer
- BD Kits and Templates
- BD Templates on the Web
- **How to Create Your Own Templates**
- Coming Soon....



Helping all people
live healthy lives

How to Create Your Own Templates

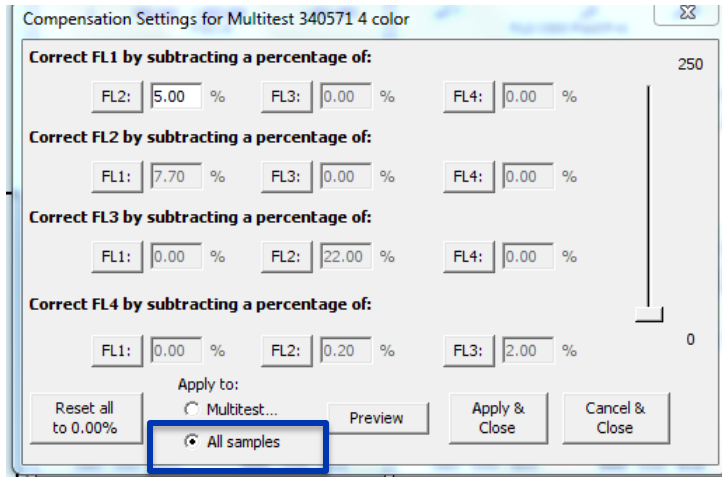


1. Open a new workspace or existing workspace.



Helping all people
live healthy lives

How to Create Your Own Templates



Compensation Settings for Multitest 340571 4 color

Correct FL1 by subtracting a percentage of: 250

FL2: 5.00 % FL3: 0.00 % FL4: 0.00 %

Correct FL2 by subtracting a percentage of:

FL1: 7.70 % FL3: 0.00 % FL4: 0.00 %

Correct FL3 by subtracting a percentage of:

FL1: 0.00 % FL2: 22.00 % FL4: 0.00 %

Correct FL4 by subtracting a percentage of:

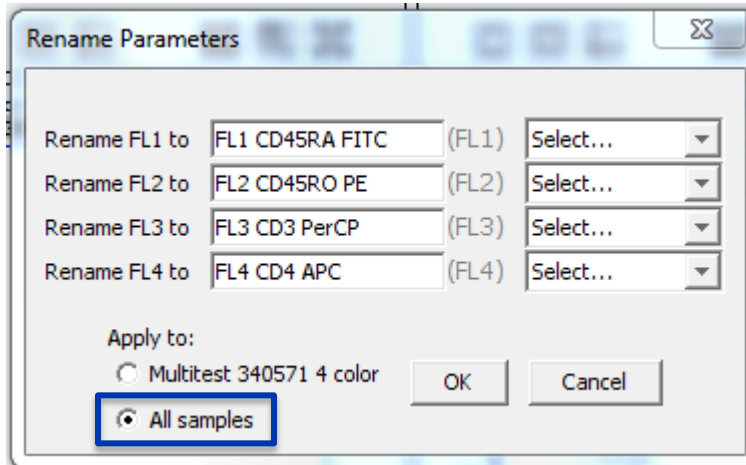
FL1: 0.00 % FL2: 0.20 % FL3: 2.00 %

Apply to:

Multitest... All samples

Reset all to 0.00% Preview Apply & Close Cancel & Close

2. Collect samples, optimize plots and run criteria to desired settings.
3. If desired, input compensation, parameter names, and threshold settings and click Apply to: All samples.



Rename Parameters

Rename FL1 to FL1 CD45RA FITC (FL1) Select...

Rename FL2 to FL2 CD45RO PE (FL2) Select...

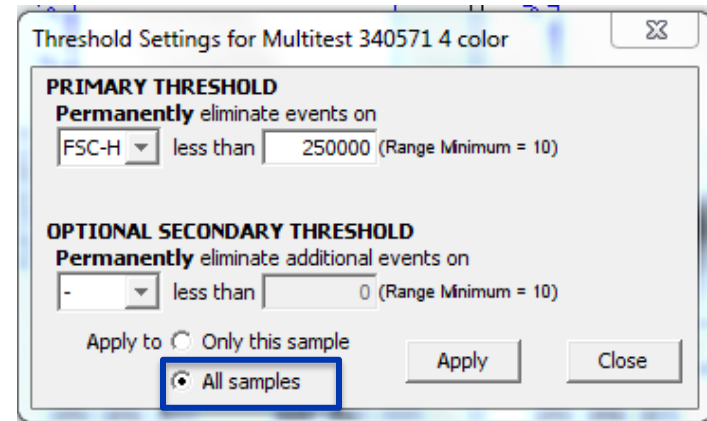
Rename FL3 to FL3 CD3 PerCP (FL3) Select...

Rename FL4 to FL4 CD4 APC (FL4) Select...

Apply to:

Multitest 340571 4 color All samples

OK Cancel



Threshold Settings for Multitest 340571 4 color

PRIMARY THRESHOLD
Permanently eliminate events on
FSC-H less than 250000 (Range Minimum = 10)

OPTIONAL SECONDARY THRESHOLD
Permanently eliminate additional events on
- less than 0 (Range Minimum = 10)

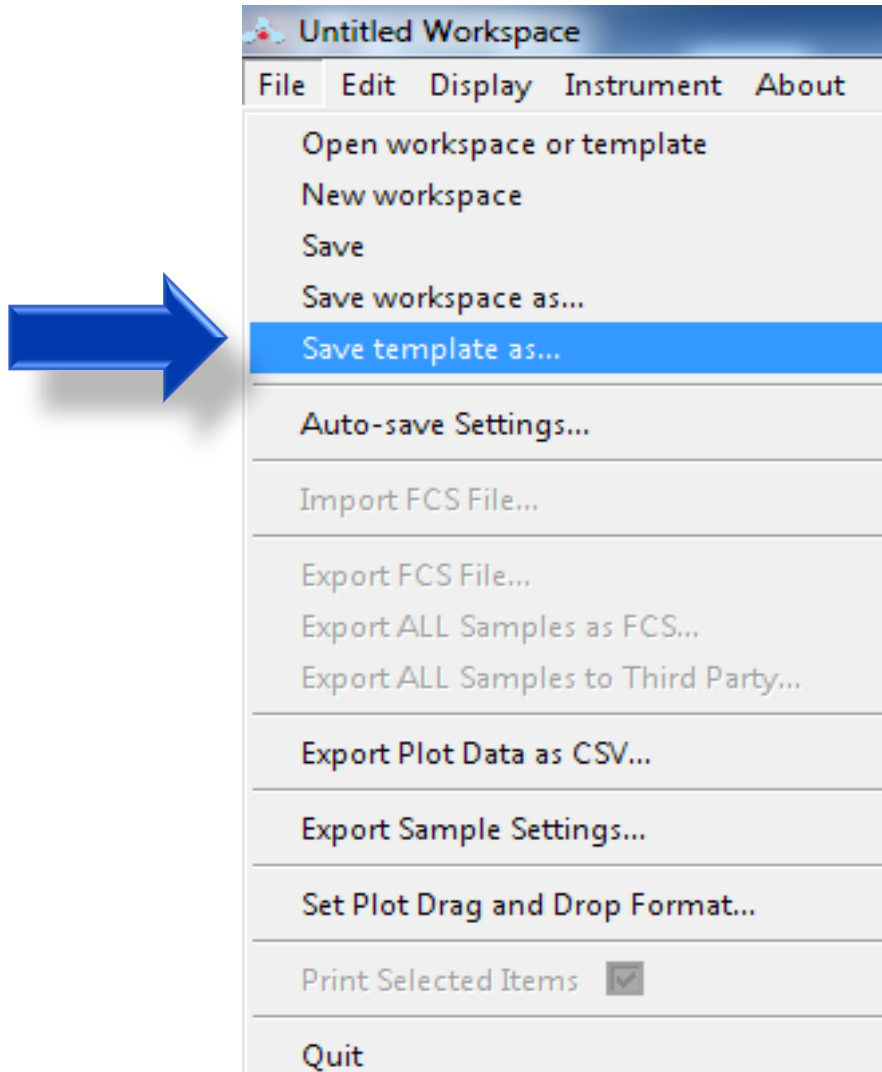
Apply to Only this sample All samples

Apply Close



Helping all people
live healthy lives

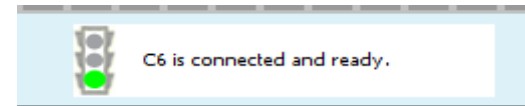
How to Create Your Own Templates



4. Select File > Save Template As.

5. Save the template with a new name.

Note: Creating templates works best at the cytometer workstation, when the BD Accuri C6 is powered on.



Helping all people
live healthy lives

Important Points to Remember

- Customize templates to fit your sample type and workflow.
- Verify all settings, including thresholds and compensation.
- When using the BD CSampler, copy desired settings to the Auto Collect tab.
- For best results, create templates when the software is connected to the cytometer.



Helping all people
live healthy lives

Outline

- Background: Immunophenotyping
- Introduction to the BD Accuri™ C6 Flow Cytometer
- BD Kits and Templates
- BD Templates on the Web
- How to Create Your Own Templates
- **Coming Soon....**



Helping all people
live healthy lives

BD Templates Coming Soon to the Web.....

COMING SOON

NEW

Tregs
Mouse Lymphocytes
Rat Lymphocytes
Mouse Tregs
Mouse T-Cell
Cytokines
Counting and Viability
of Bacteria
Mesenchymal Stem
Cell Kit

Proliferation: BrdU
Apoptosis: Annexin V
Apoptosis: Caspase-3
DNA Analysis
BD MitoScreen
Stem Cell Flow Kits
T-Cell Cytokines
Naïve/memory T Cells

EXISTING

6, 8-Peak Bead
BD CBA
Water Quality



Helping all people
live healthy lives

Special Promotional Offer

For a limited time, take advantage of **big savings** and value-added packages for the BD Accuri C6 flow cytometer with our reagent kit discount.

Get a 40% discount on reagent kits now through June 2014

New to BD Accuri?

Buy a BD Accuri C6 now and receive an immediate **10% discount** off the purchase price. You'll also be eligible to receive **40% off the list price of all BD Pharmingen reagents** you purchase for use on the BD Accuri C6 for the next 2 years.

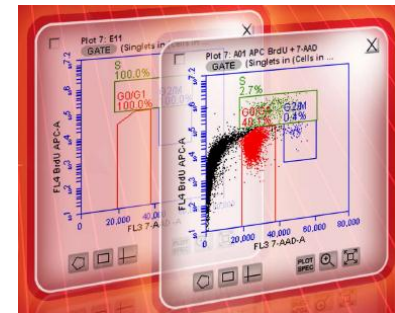
Offer expires June 30th, 2014.

Now you can simplify your workflow while dramatically reducing the cost of use of your personal flow cytometer.

www.bdbiosciences.com/go/templates

Summary

- The **BD Accuri C6** is making it even easier to apply the power of flow cytometry to your research with **free software templates** and **ready-to-go reagent kits** specific to your studies.
- Find out more about how the BD Accuri C6 puts the power of 4-color cell analysis within reach by visiting www.bdbiosciences.com/go/templates



Flow Cytometry Within Reach.



Helping all people
live healthy lives

For more information...

If you have further questions:

Contact Technical Support (US) at:

877-232-8995, Prompt 3, 2

or email: ResearchApplications@bd.com

Please visit our BD Accuri resources site at:

www.bdbiosciences.com/resources/accuri.

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

Class 1 Laser Product.

Alexa Fluor® is a registered trademark of Life Technologies Corporation.

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

BD, BD Logo and all other trademarks are property of Becton, Dickinson and Company. © 2014 BD



Helping all people
live healthy lives