

# Optimizing Intracellular Flow Cytometry:

Simultaneous Detection of Cytokines and  
Transcription Factors

Presented by  
Jurg Rohrer, PhD, BD Biosciences

# Outline

- Introduction
  - Cytokines
  - Transcription factors
- Basic concepts of intracellular flow cytometry
  - Optimization examples
- Treg/Th17 cell analysis
  - Considerations
  - Examples



# Cytokines

- Soluble polypeptides produced by most nucleated cells in the body
- Some potent producers include endothelial and epithelial cells and resident macrophages, especially near the interface with the external environment
- Critical to the development and functioning of both the innate and adaptive immune responses
- Promote cellular differentiation and proliferation
  - Example: IL-2 involved in T cell activation and maintenance of a Th1 response
- Work in either an autocrine or paracrine manner



# Th17 Cells

- A subset of CD4<sup>+</sup> T helper cells
- Developmentally distinct from Th1 and Th2 cells
- Immunity against bacterial and fungal infectious
- Play a key role in autoimmune diseases (tissue injury)
- Controlling Th17 activity could aid in the treatment of autoimmune diseases
- TGF- $\beta$ , IL-6, IL-21, IL-1 $\beta$ , and IL-23 appear to drive Th17 development
- Produce IL-17A, IL-17F; also IL-21, IL-22, IL-26, and less TNF and IL-6



# Transcription Factors

- Proteins that bind to specific DNA sequences
- Control the transfer of genetic information from DNA to RNA
- Regulators of gene expression
- A single transcription factor can bind hundreds of promoters



# Regulatory T Cells

- Tregs = CD4<sup>+</sup> T regulatory cells
- Comprise ~ 1–3% of human PBMCs and ~ 4–8% of mouse spleen
- Actively suppress T cell proliferation
- Play a crucial role in T cell homeostasis
- nTreg develop in the thymus, iTreg require TGFβ, IL-2 and RA
- FoxP3, a forkhead family transcription factor, is a specific marker for Tregs
- FoxP3 is necessary for the development and function of Tregs

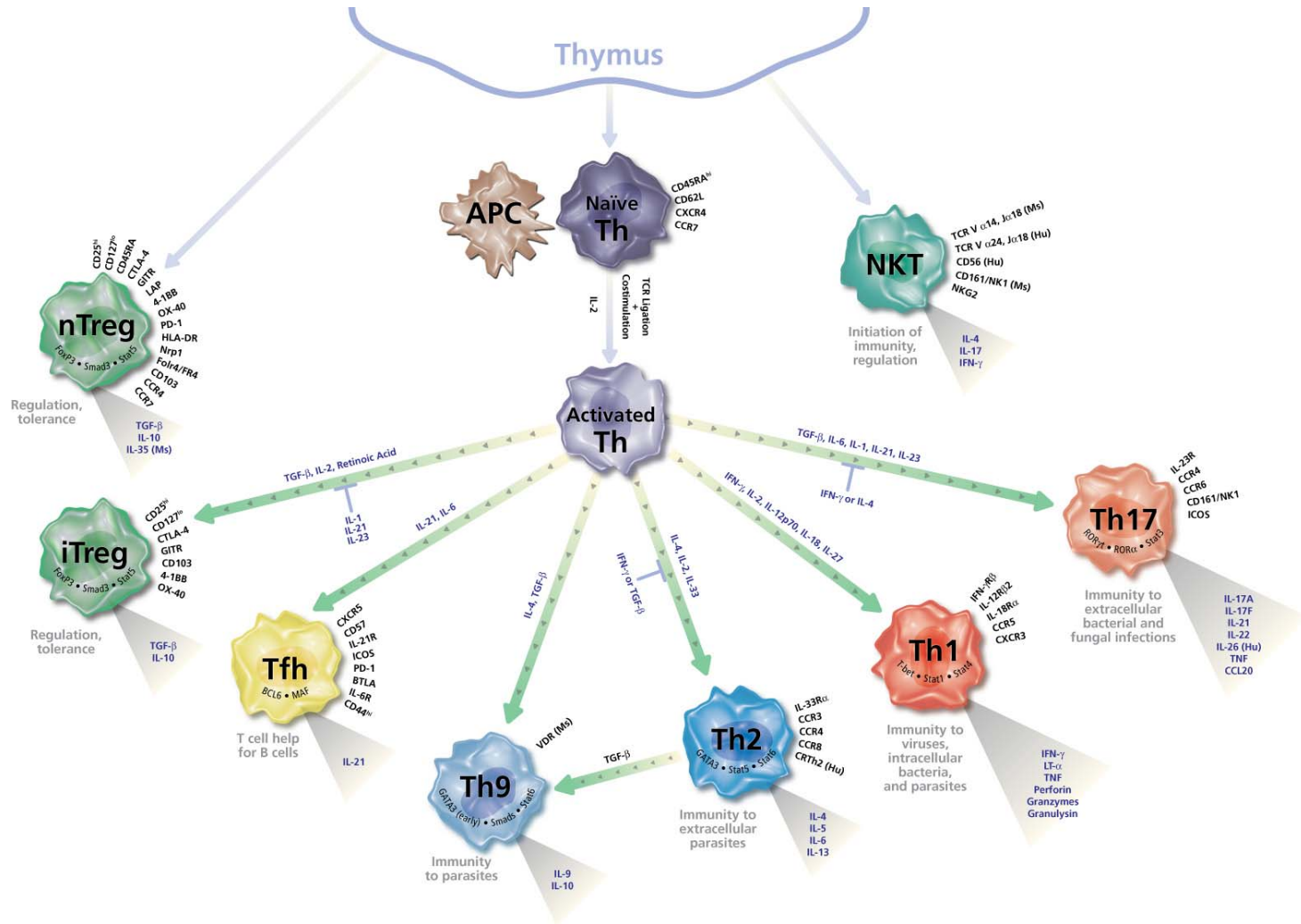


# Regulatory T Cells, *cont'd*

- Produce TGF $\beta$  and IL-10 and express high levels of CD25 and low levels of CD127
- Diminish immune responses against cancers, allogeneic transplants, and infectious pathogens
- Dampening Treg activity could improve anti-tumor responses and responses to vaccinations and chronic infections
- Deficiencies contribute to the development of autoimmune diseases
- Boosting Treg activity could be useful in the treatment of T cell induced diseases



# CD4<sup>+</sup> T Cell Differentiation





# What is Intracellular Flow Cytometry?

- Detection of:
  - Transcription factors
  - Intracellular signaling molecules
  - Cytokines
  - Structural proteins
  - Scaffold proteins
  - Pan and phospho-specific antigens



# Considerations for Intracellular Flow Cytometry

- Must permeabilize a cell to access cell contents
- If a cell is permeabilized, then contents could “leak” out and the protein of interest could be lost
- Therefore, cells are fixed first, followed by permeabilization
- To detect secreted proteins, they must be “trapped” within the cell prior to fixation and permeabilization to increase the likelihood of detection



# Considerations for Intracellular Flow, *cont'd*

- Protein transport inhibition
  - Monensin vs Brefeldin A (BD GolgiStop™ vs BD GolgiPlug™ inhibitor)
  - Optimal time for inhibition
  - Optimal concentration of inhibitor
- Fixation
  - Concentration (paraformaldehyde)
  - Time
  - Temperature
  - Compatibility with fluorochromes
  - Compatibility of cell surface markers



# Considerations for Intracellular Flow, *cont'd.*

- Permeabilization
  - Perm agent (saponin, methanol, Tween<sup>®</sup> 20, Triton X-100<sup>™</sup>)
  - Concentration
  - Time
  - Temperature
  - Compatibility with fluorochromes
  - Compatibility of cell surface markers
- Different locations in cells are more difficult to access
- Types of proteins being identified, single or in a complex?



# Considerations for Intracellular Flow, *cont'd.*

- Antibody staining
  - Order
  - Concentration
  - Time
  - Temperature
  - Fluorochromes
- Storage conditions
  - Buffer
  - Time
- Matching one antibody protocol with another antibody protocol



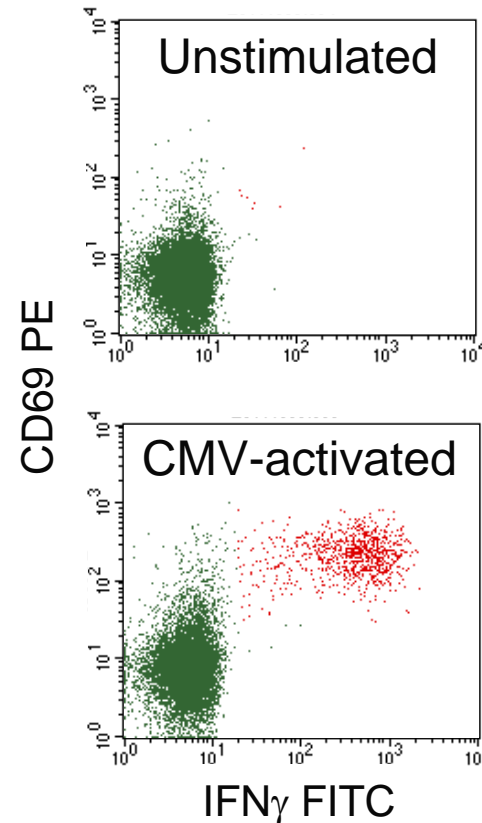
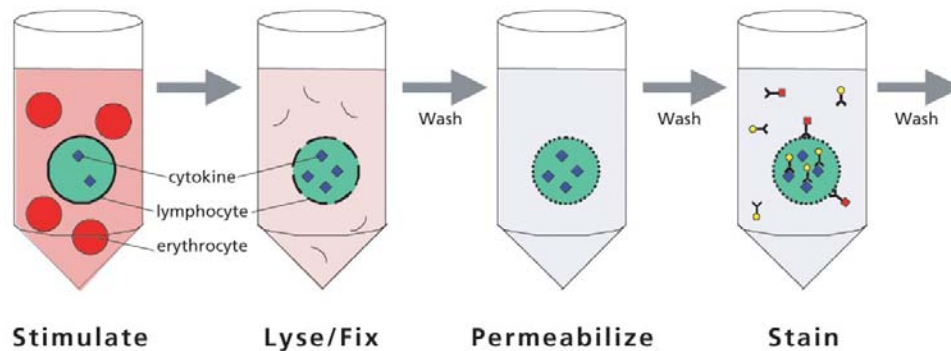
# Buffer Choices

- Fixation buffer
- BD Cytotfix/Cytoperm™ and BD™ Perm/Wash buffer
- BD Pharmingen™ FoxP3 buffer set (mouse or human)
- BD™ Phosflow Perm Buffer II
- BD™ Phosflow Perm Buffer III
- BD IntraSure™ kit
  
- BD FastImmune™ kits



# BD FastImmune™ Kits

- Optimized kits containing antibodies and buffers for simultaneous detection of cell surface markers and cytokines from whole blood



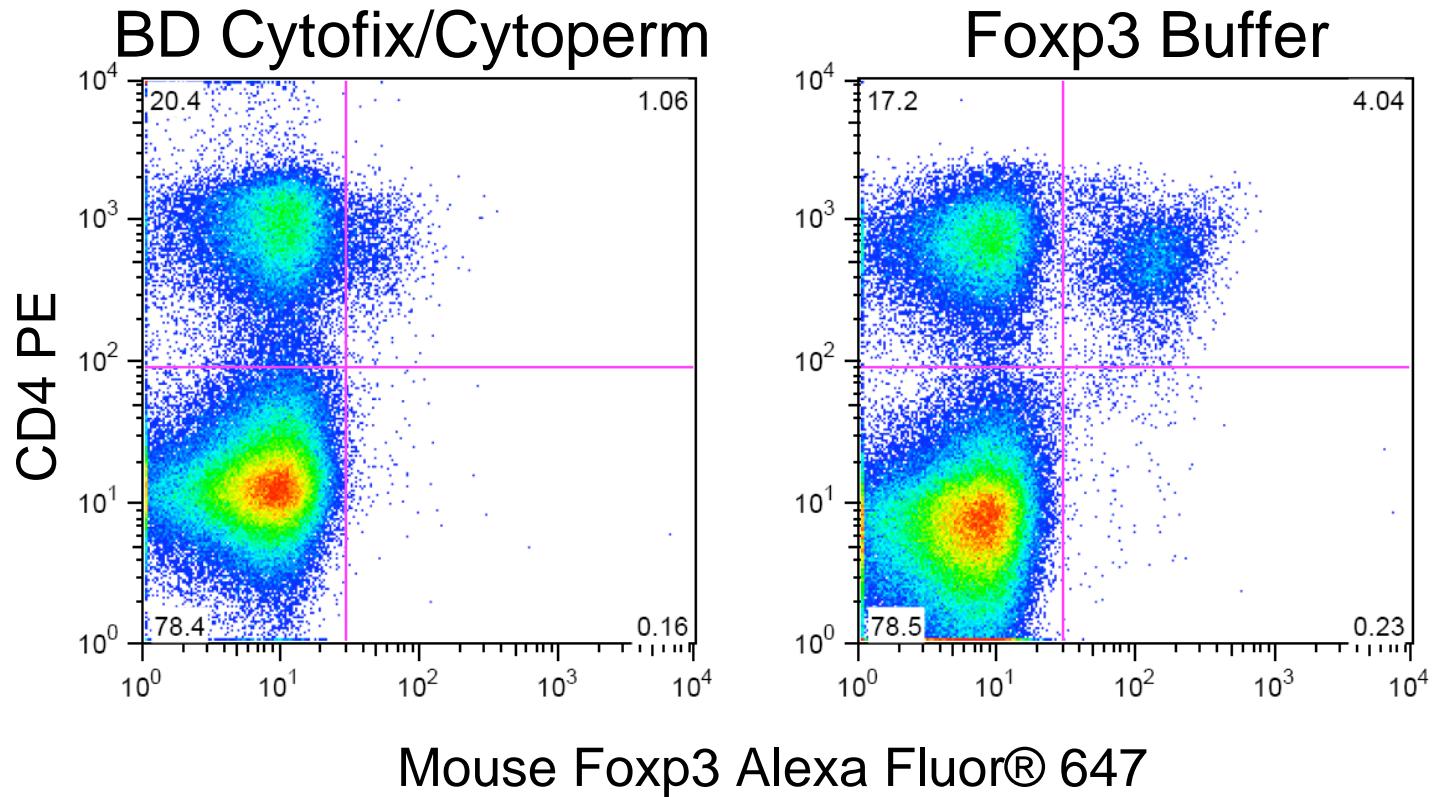
# Case Study

- The study of Treg and Th17 cells
- Requires the need to detect both FoxP3 and IL-17 in the same sample
- Unique protocols for both mouse and human FoxP3 staining
- Questions are:
  - How well does IL-17 staining work in the FoxP3 buffer system?
  - How well do other intracellular and surface markers work with the FoxP3 buffer system?
- Examples of FoxP3 optimization followed by addition of other markers

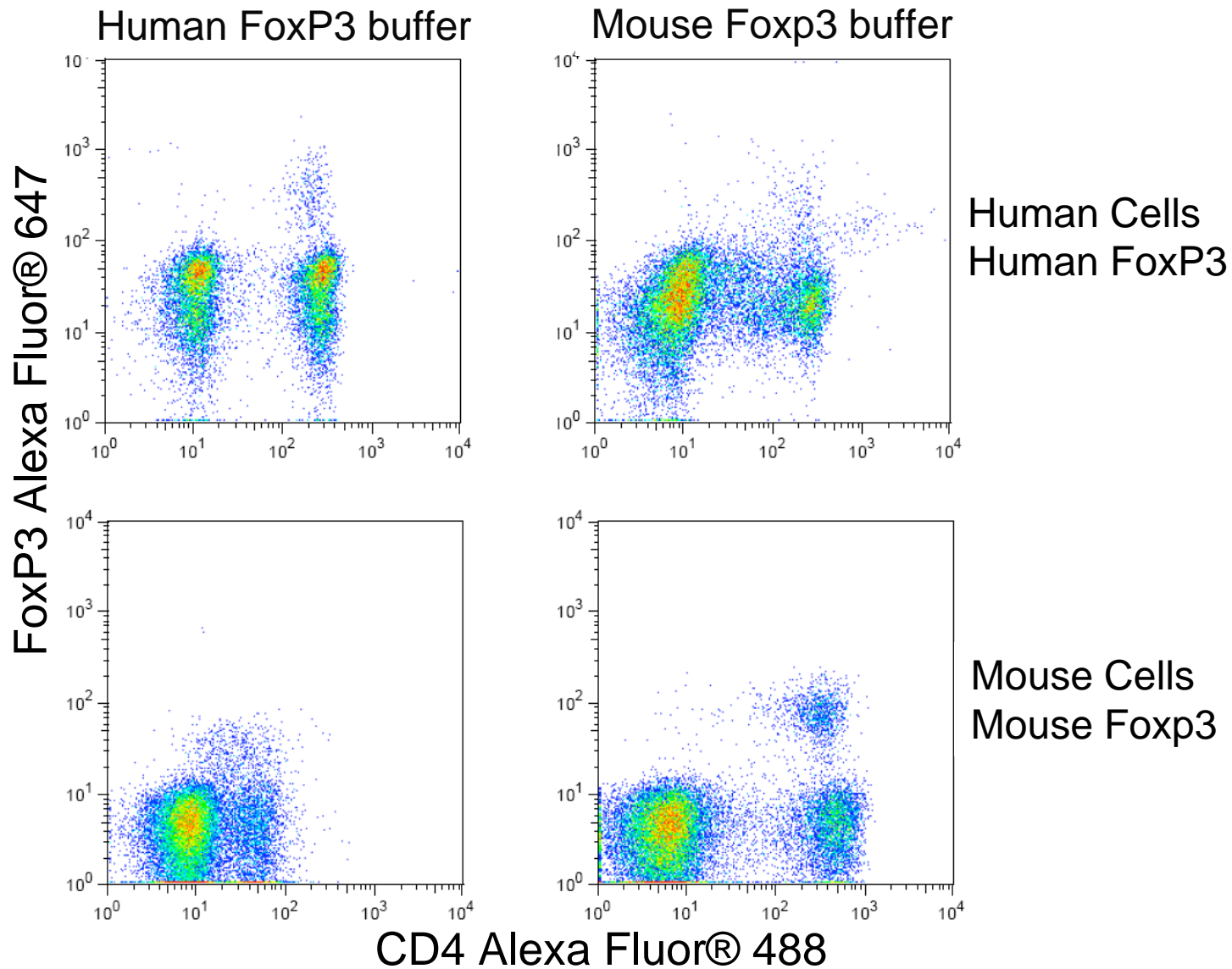




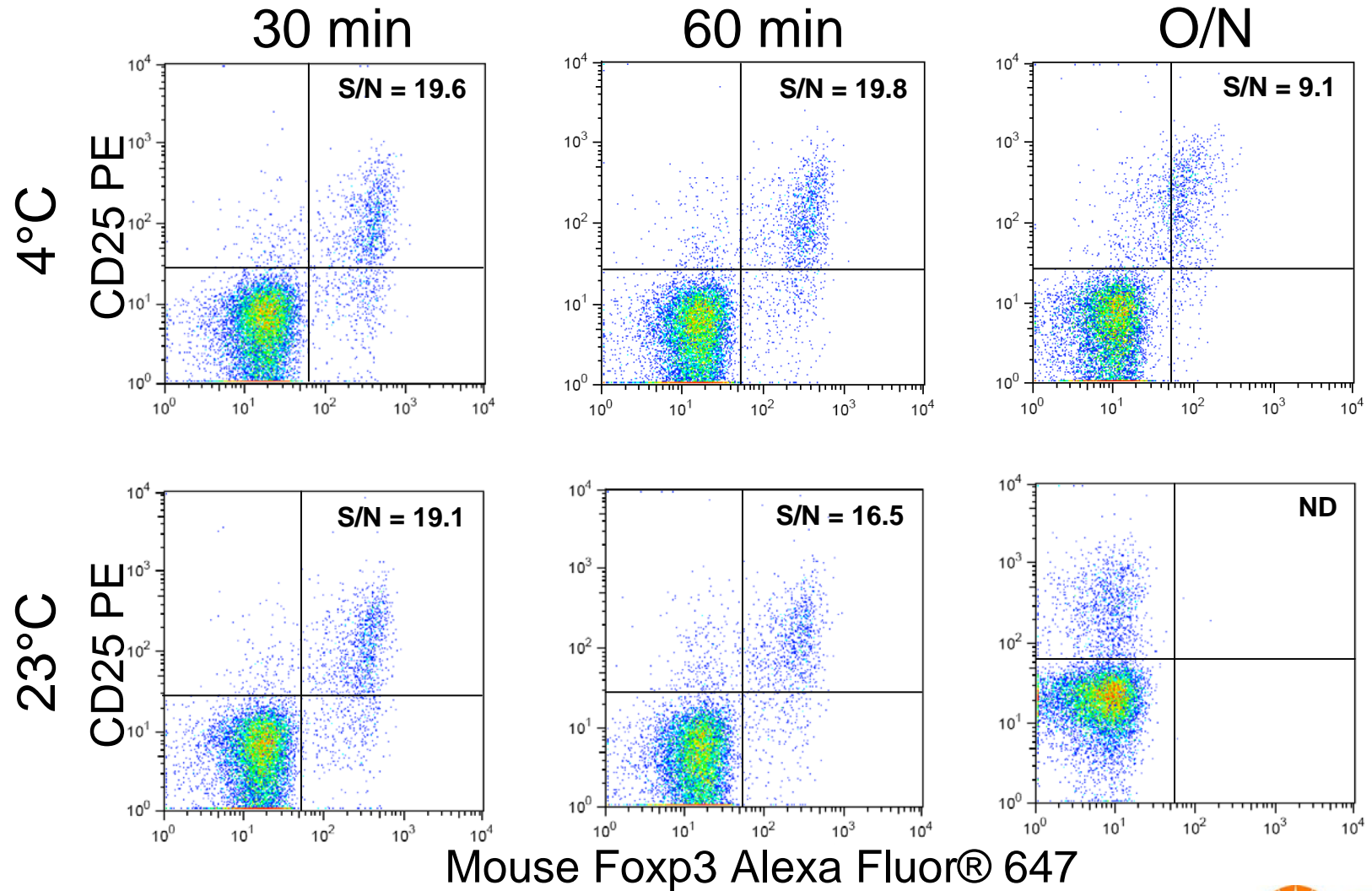
# Effect of BD Cytotfix/Cytoperm Buffer on Mouse Foxp3 Staining



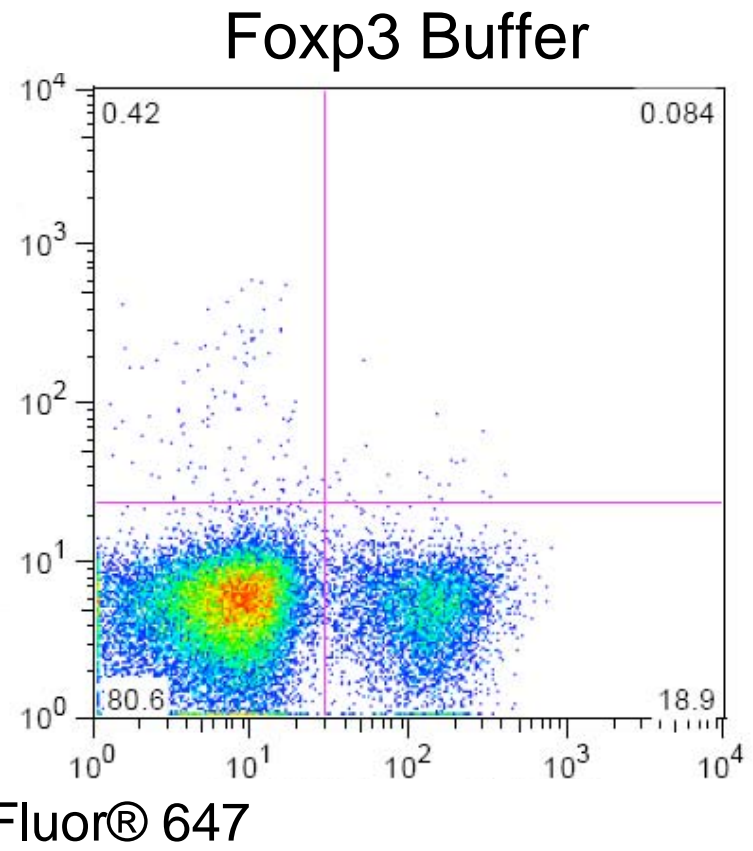
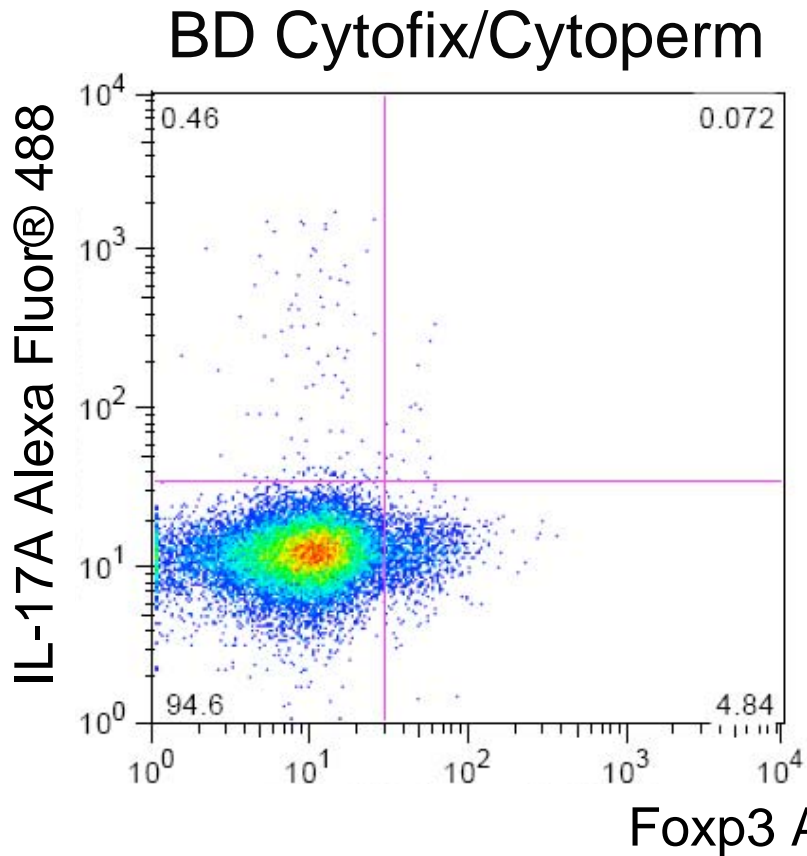
# Effect of Human FoxP3 Buffer System on Mouse Foxp3 Staining



# Effect of Fixation Time and Temperature on Mouse Foxp3 Staining



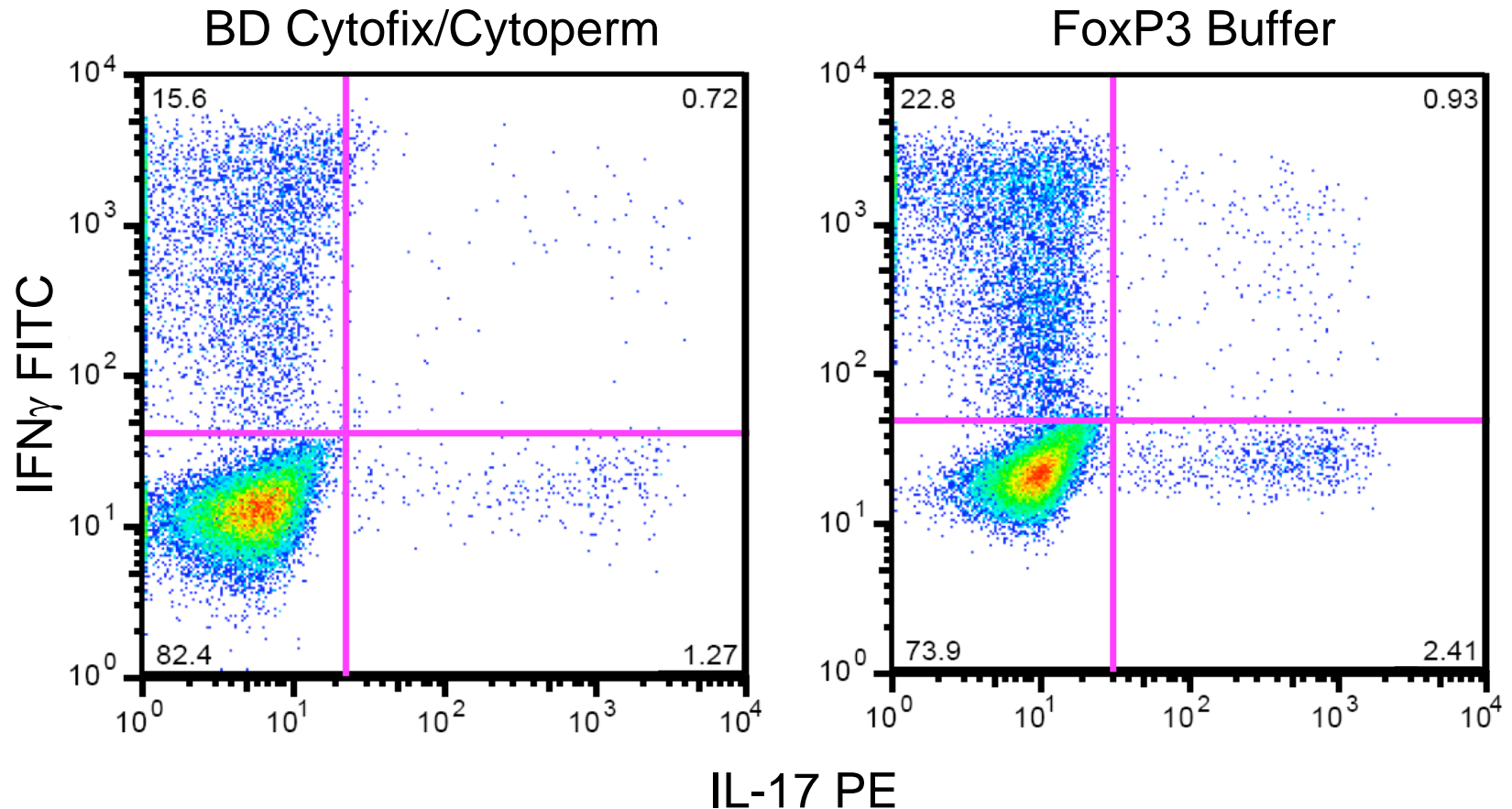
# Effect of FoxP3 Buffer on Mouse IL-17 Staining



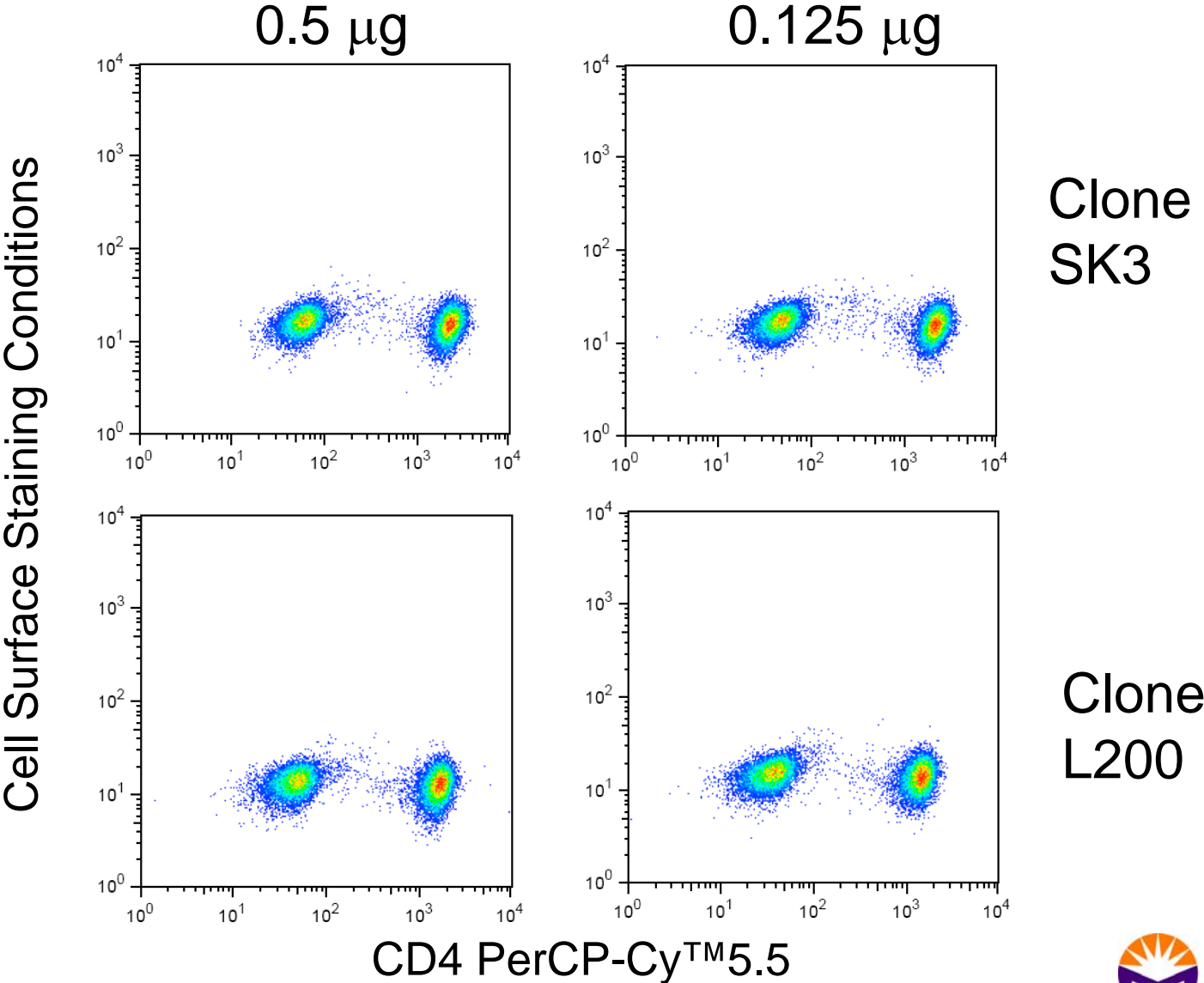
Gated on CD4<sup>+</sup> lymphocytes



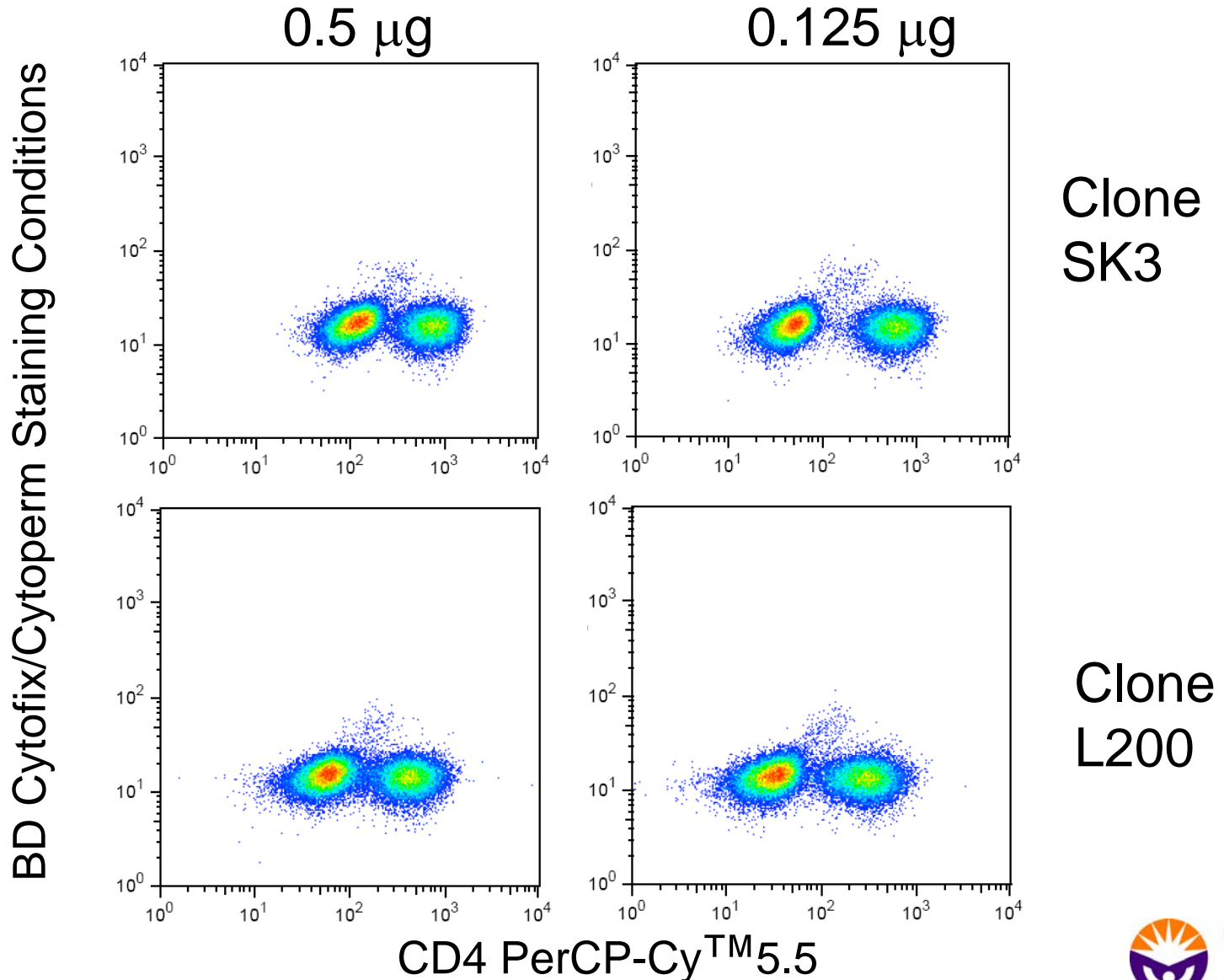
# Effect of FoxP3 Buffer on Human Cytokine Staining



# Optimizing Cell Surface Staining – Surface Stain



# Optimizing Cell Surface Staining – BD Cytofix/Cytoperm Stain



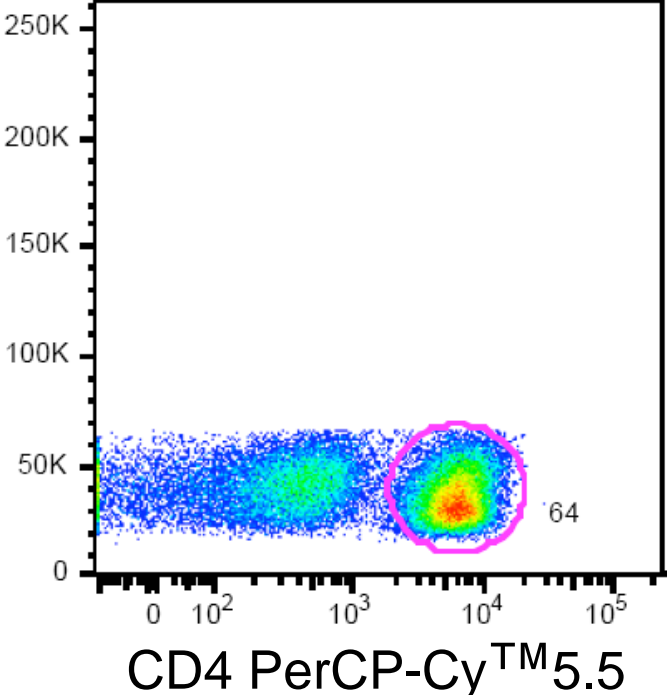
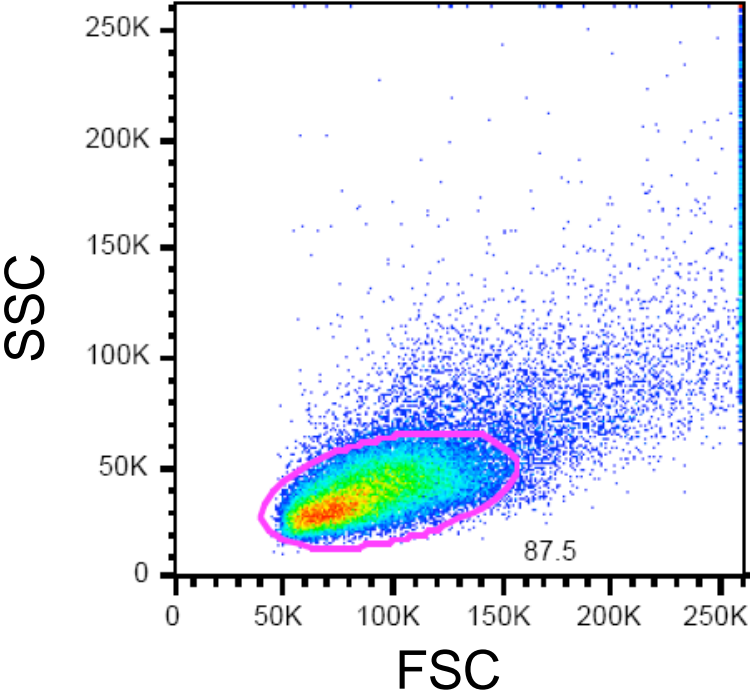
## Example: Simultaneous detection of human FoxP3, IL-17, IL-4, and IFN $\gamma$ in CD4<sup>+</sup> T cells.

- Freshly isolated PBMC
- Either stimulated or not
  - PMA/Ionomycin with GolgiStop™
  - 5 hours 37°C
- Fix (2 ways) and stored O/N in stain buffer
- Perm (2 ways) and stain 40 minutes
  - CD4 PerCP-Cy5.5
  - FoxP3 V450
  - IL-17 Alexa Fluor® 647
  - IFN $\gamma$  FITC
  - IL-4 PE
- Acquire and analyze

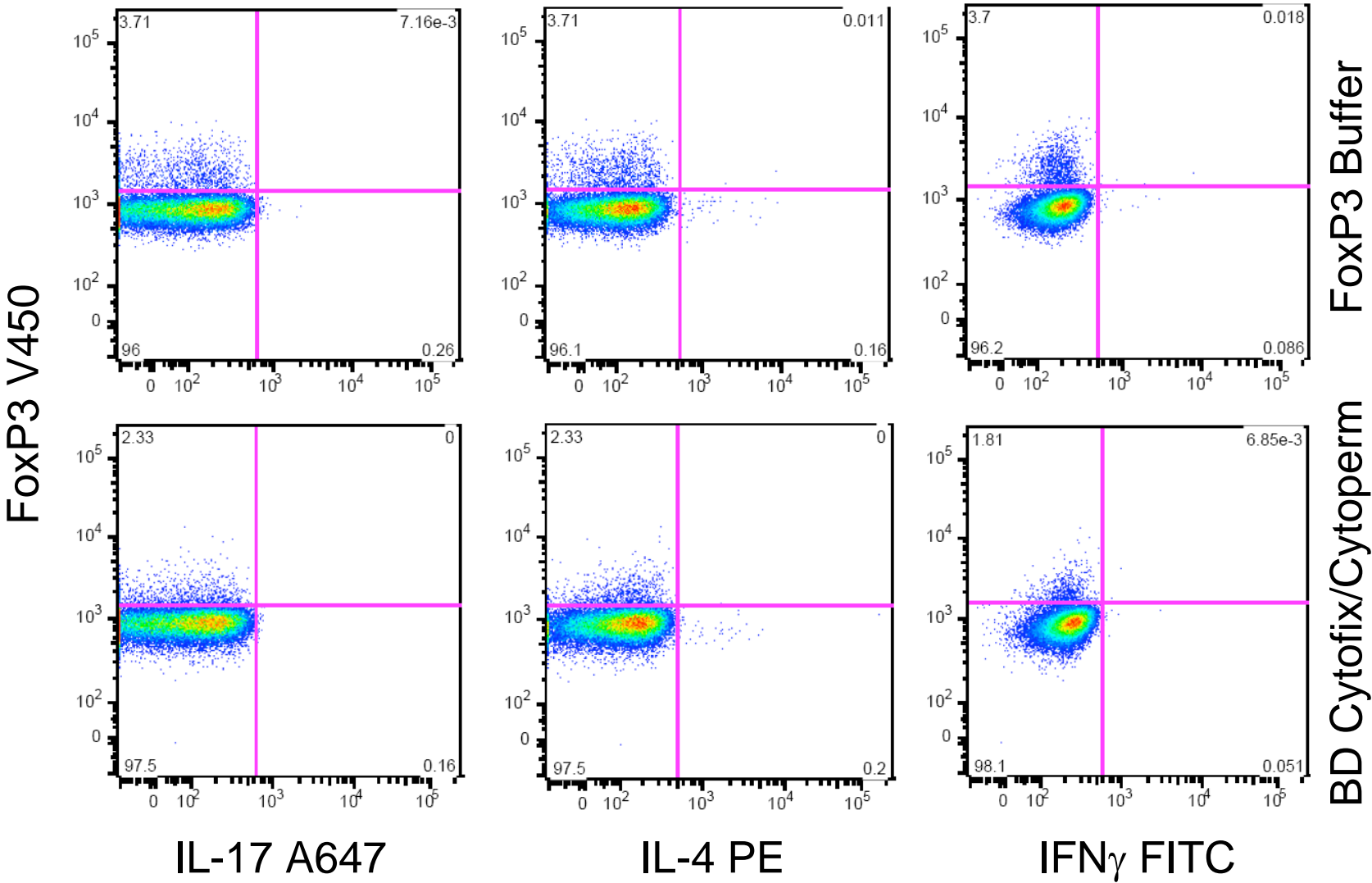




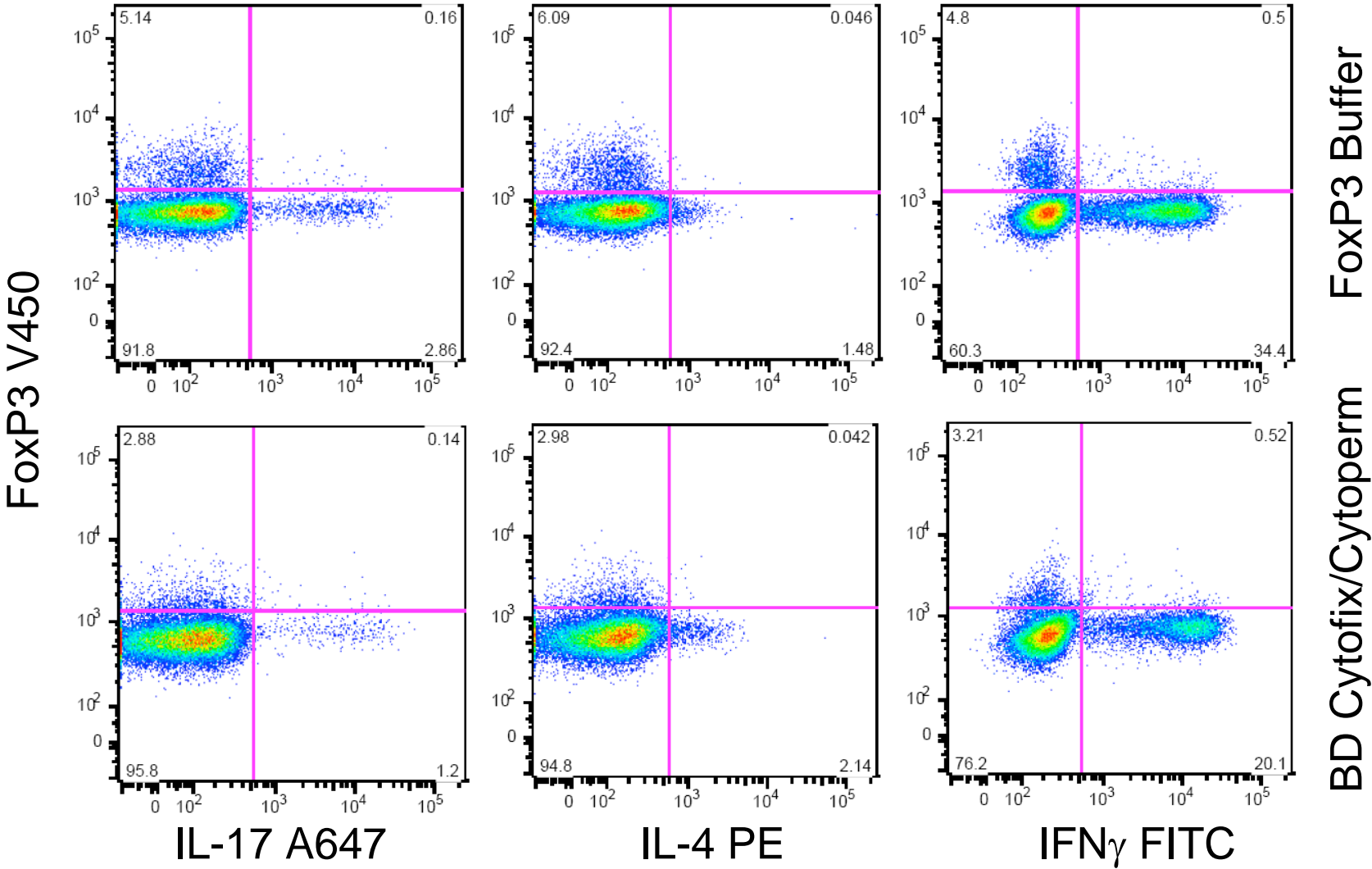
# Setting the CD4+ gate



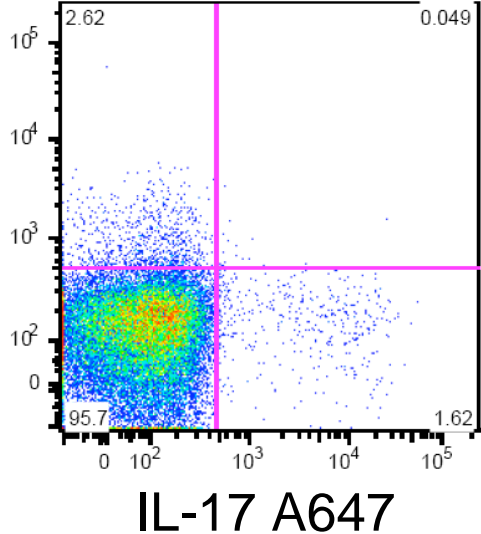
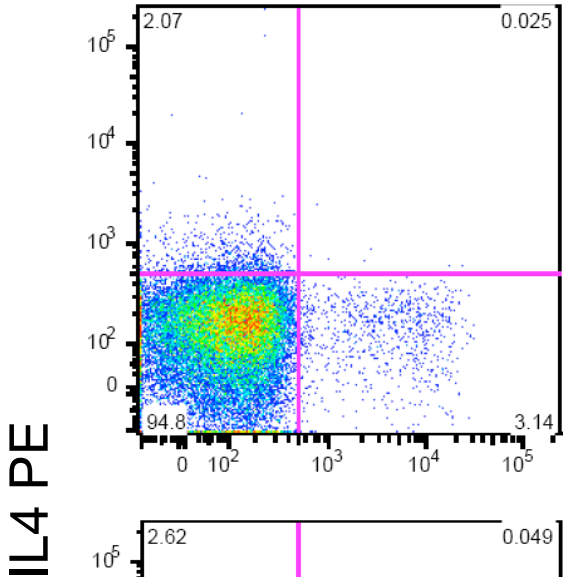
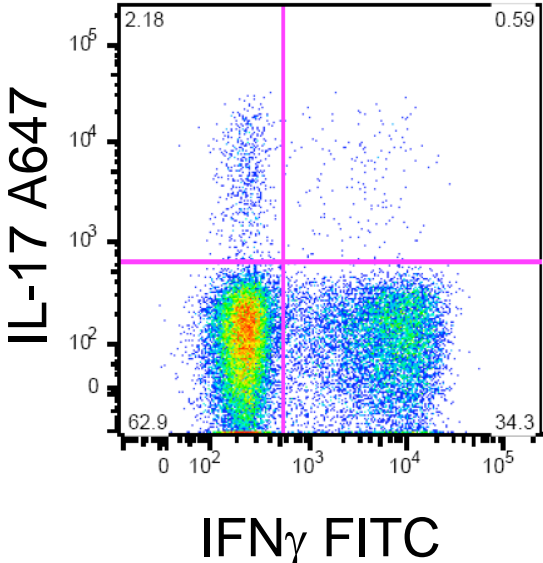
# Unstimulated PBMC



# Stimulated PBMC



# Stimulated PBMC, *cont'd.*



FoxP3 Buffer

BD Cytotfix/Cytoperm

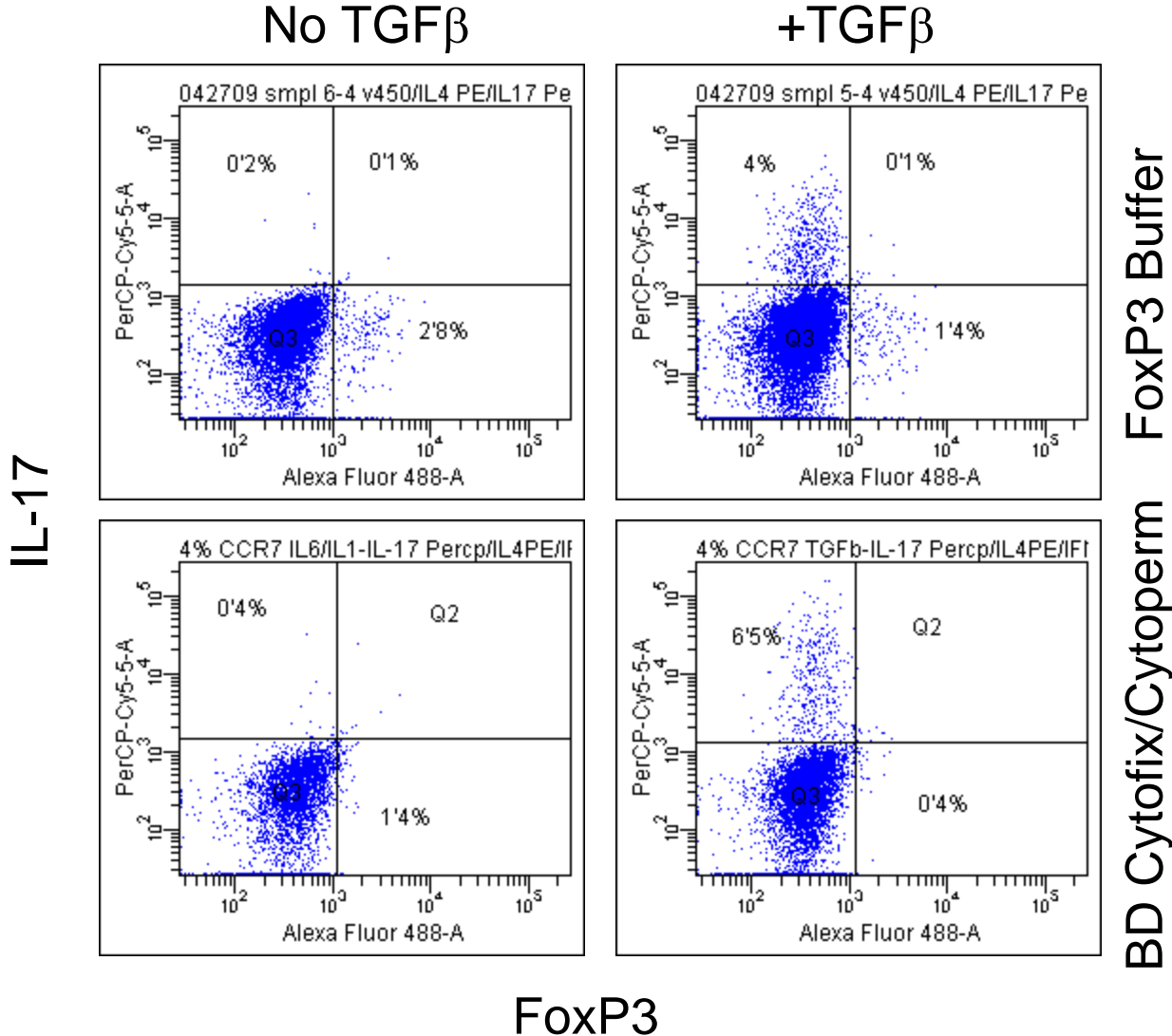


## Example: Requirement of TGF $\beta$ for the differentiation of mouse Th17 CD4<sup>+</sup> T cells.

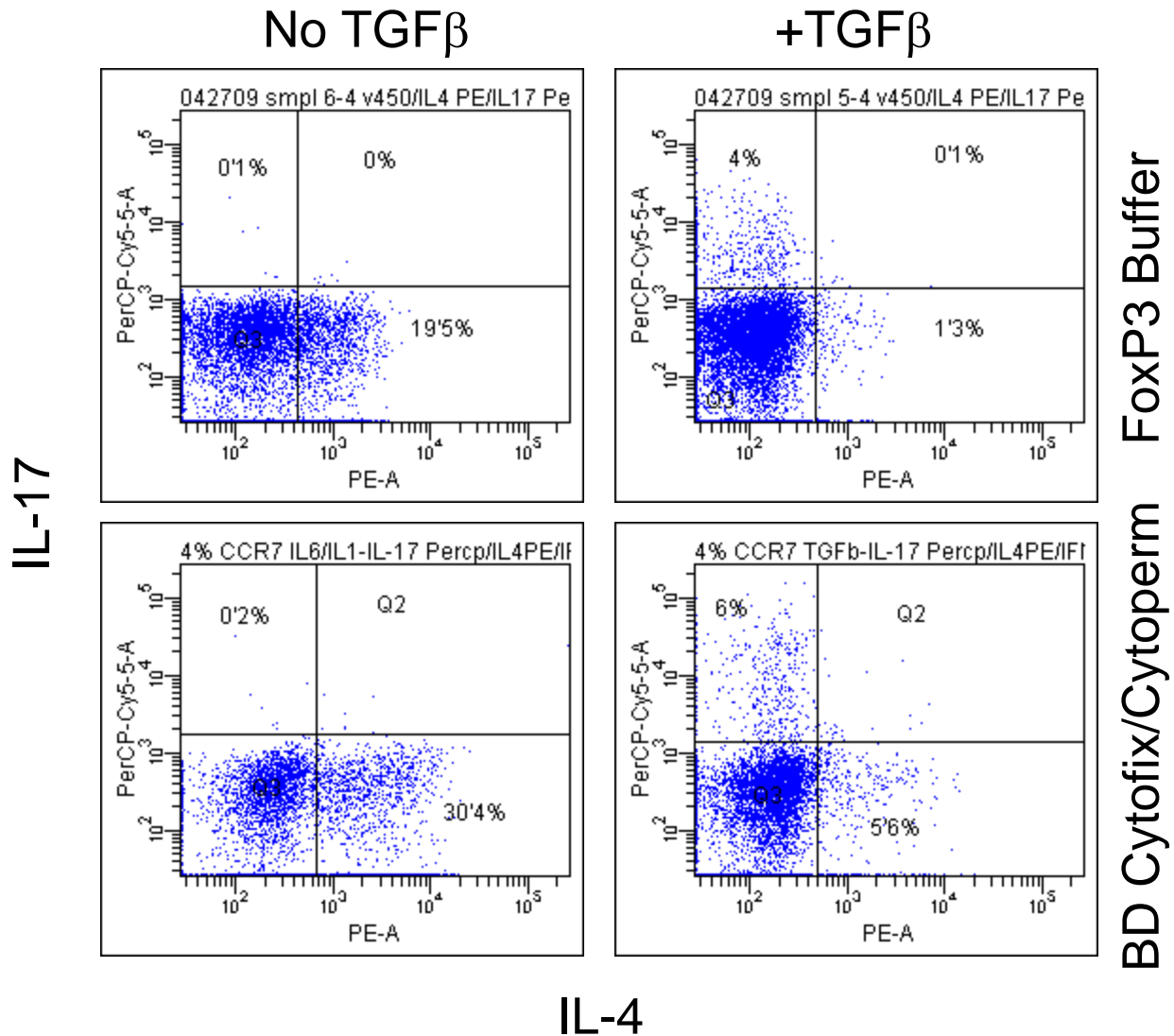
- Freshly isolated spleen
- Purify CD4<sup>+</sup> T cells by panning
- Polarize T cells on anti-CD3 coated plates in the presence of CD28, IL-6 and IL1 $\beta$  either with or without TGF $\beta$
- After 4 days harvest the cells and stimulate with PMA/Ionomycin with GolgiStop<sup>TM</sup> for 5 hours
- Fix (2 ways) and store O/N in stain buffer
- Perm (2 ways) and stain 40 minutes
  - CD4 V450
  - FoxP3 Alexa Fluor<sup>®</sup> 488
  - IL-17 PerCP-Cy<sup>TM</sup>5.5
  - IL-4 PE
- Acquire and analyze



# Differentiated CD4+ T cells



# Differentiated CD4<sup>+</sup> T cells, *cont'd.*



IL-17

IL-4

FoxP3 Buffer  
BD Cytotfix/Cytoperm



# Summary

- Determine marker combination(s) for your experiment
- Pair the brightest dye with dimmest marker
- Determine optimal buffers for your antibodies
- Begin cross testing antibodies in different buffers
  - Typically optimize conditions for intracellular staining first and then determine what works best for your chosen cell surface markers
  - Understand what compromises can be made
- Once optimal conditions have been determined for your particular needs, proceed with experiments





# Acknowledgements

- Xiao-Wei Wu
- Ai-Li Wei
- Li Li
- Ravi Hingorani
- Jeanne Elia
- Christopher Boyce



If you have further questions:

**Contact your US Reagent Sales Rep**  
or e-mail: [ResearchApplications@bd.com](mailto:ResearchApplications@bd.com)

Alexa Fluor® is a registered trademark of Molecular Probes, Inc.

Cy™ is a trademark of Amersham Biosciences.

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

