

Simultaneous correlation of cytokine production with Treg and Th17 cell proliferation

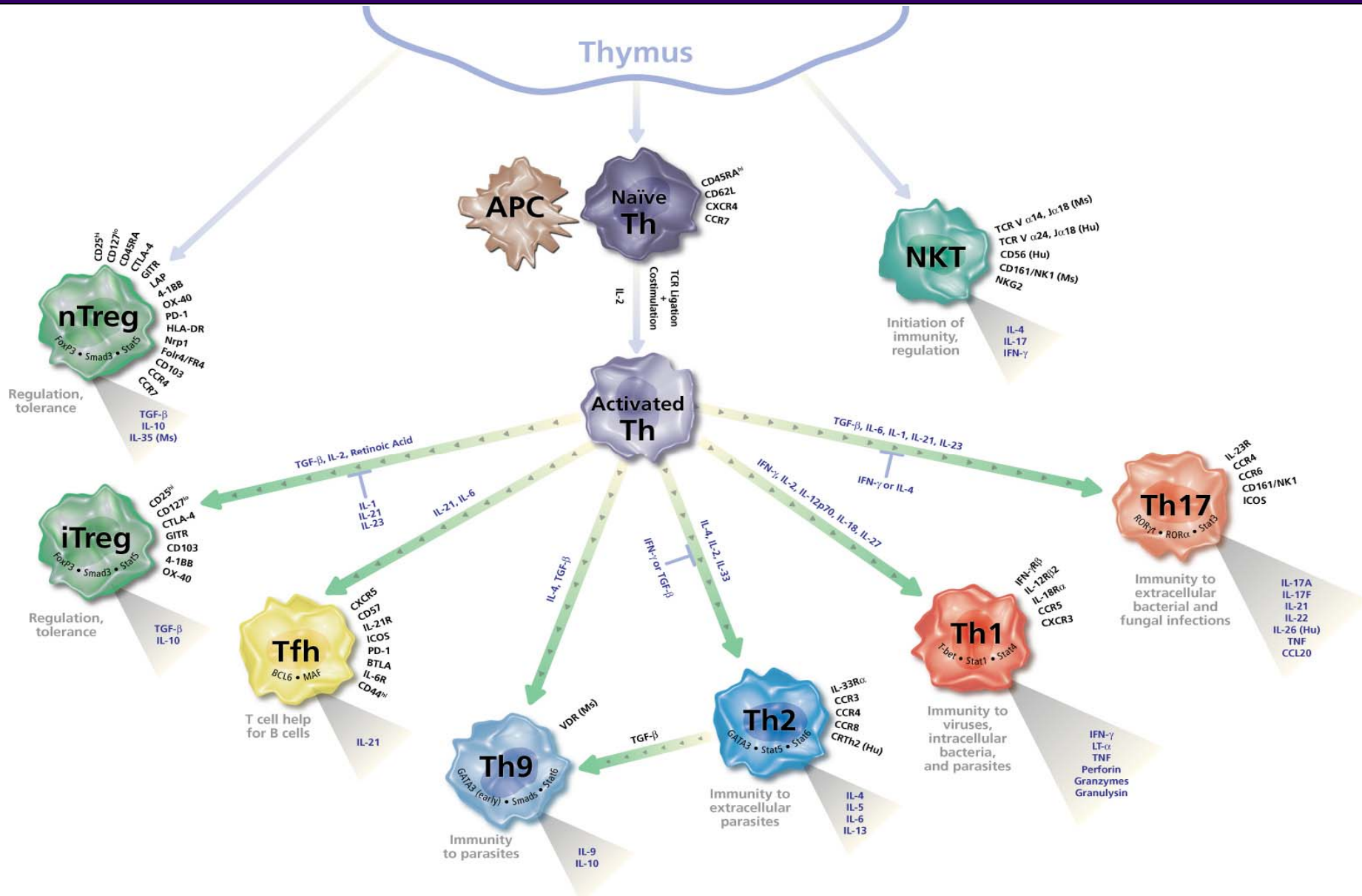
Jurg Rohrer, PhD
Director, R&D
BD Biosciences



Overview

- T helper (Th) cell overview
- Experimental setup
- Data analysis
- Conclusions

Introduction to Th biology

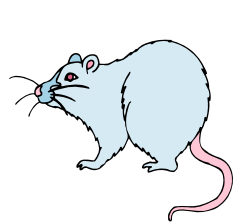


Experimental setup

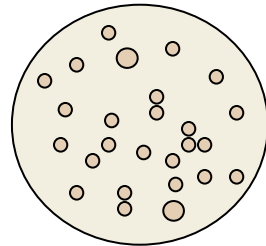
- Enrich Balb/c splenocytes by positive selection via CD4⁺ panning
- Load isolated cells with VFSE 1 μ M, 10 minutes
- Set up cultures as follows:
 - CD3/CD28
 - CD3/CD28/IL-6/IL-1 β
 - CD3/CD28/IL-6/IL-1 β /TGF- β
 - CD3/CD28/IL-6/IL-1 β /TGF- β /IL-23
- Harvest cells at 1, 2, 3, and 4 days
- Fix/perm and stain cells for IL-17A, Foxp3, IL-4, IL-2, and interferon- γ (IFN- γ)



Experimental setup continued



Harvest
Spleen



CD4 cells enriched
by panning

Cells loaded with
VFSE and washed

anti-CD3e+
anti-CD28

anti-CD3e+
anti-CD28+IL-1 β +IL-6

anti-CD3e+
anti-CD28+IL-1 β +IL-6+ TGF- β

anti-CD3e+
anti-CD28+IL-1 β +IL-6+ TGF- β + IL-23

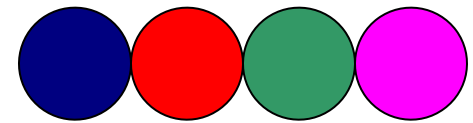
BD Cytofix/Cytoperm™ buffer
Cytokines

+ Monensin

Foxp3 Fix/Perm buffer
Foxp3 and some
cytokines

Harvest
Stimulate with PMA and
Ionomycin for 4–5 hours
+ or – Monensin

- Monensin



Day 1, 2, 3, 4

Supernatants: Analysis with BD™ CBA Flex Sets
Cells: BD™ Phosflow Fix and Perm buffer



VFSE histograms

Condition:

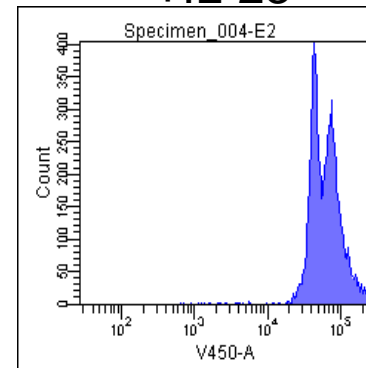
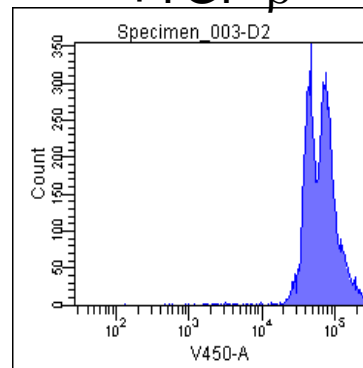
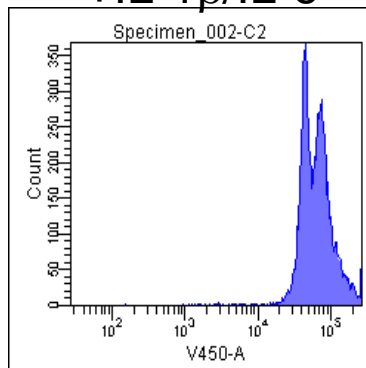
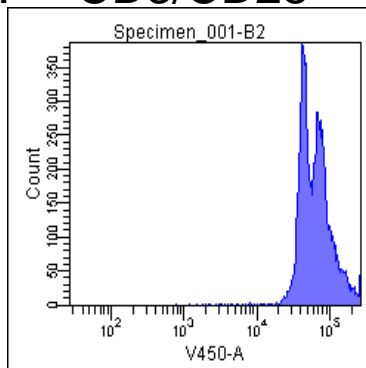
CD3/CD28

+IL-1 β /IL-6

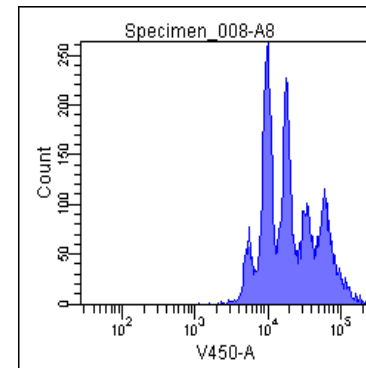
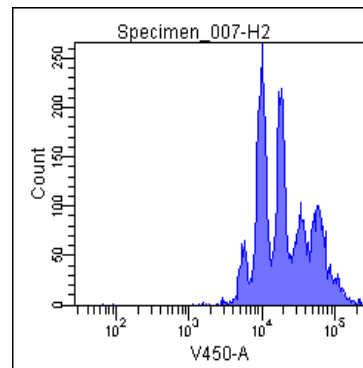
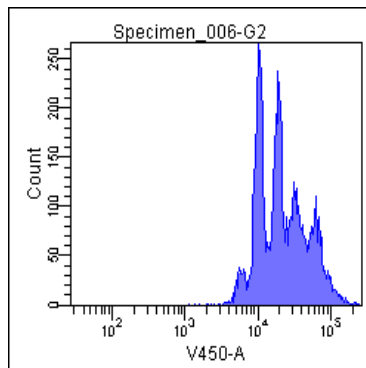
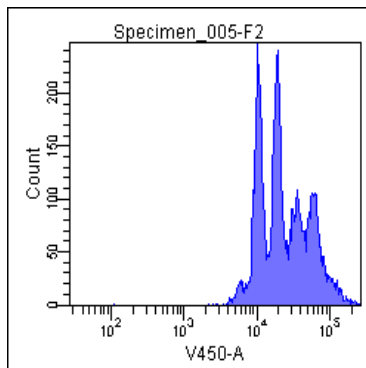
+TGF- β

+IL-23

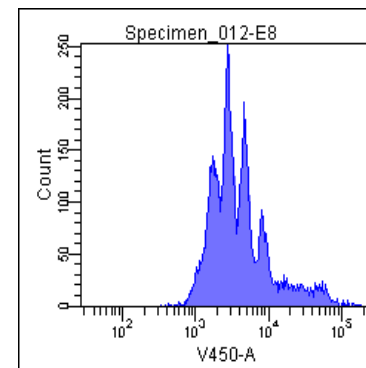
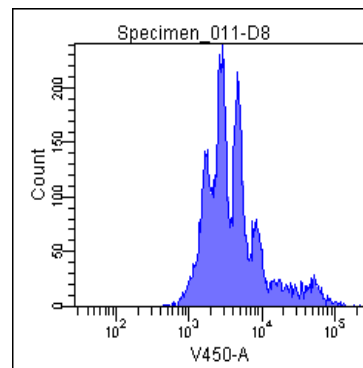
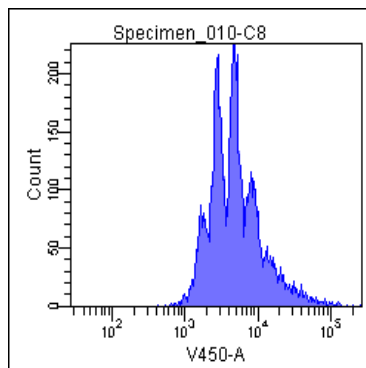
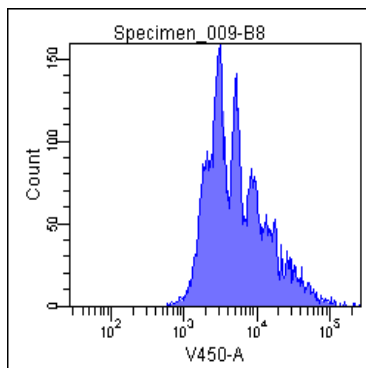
Day 1



Day 2

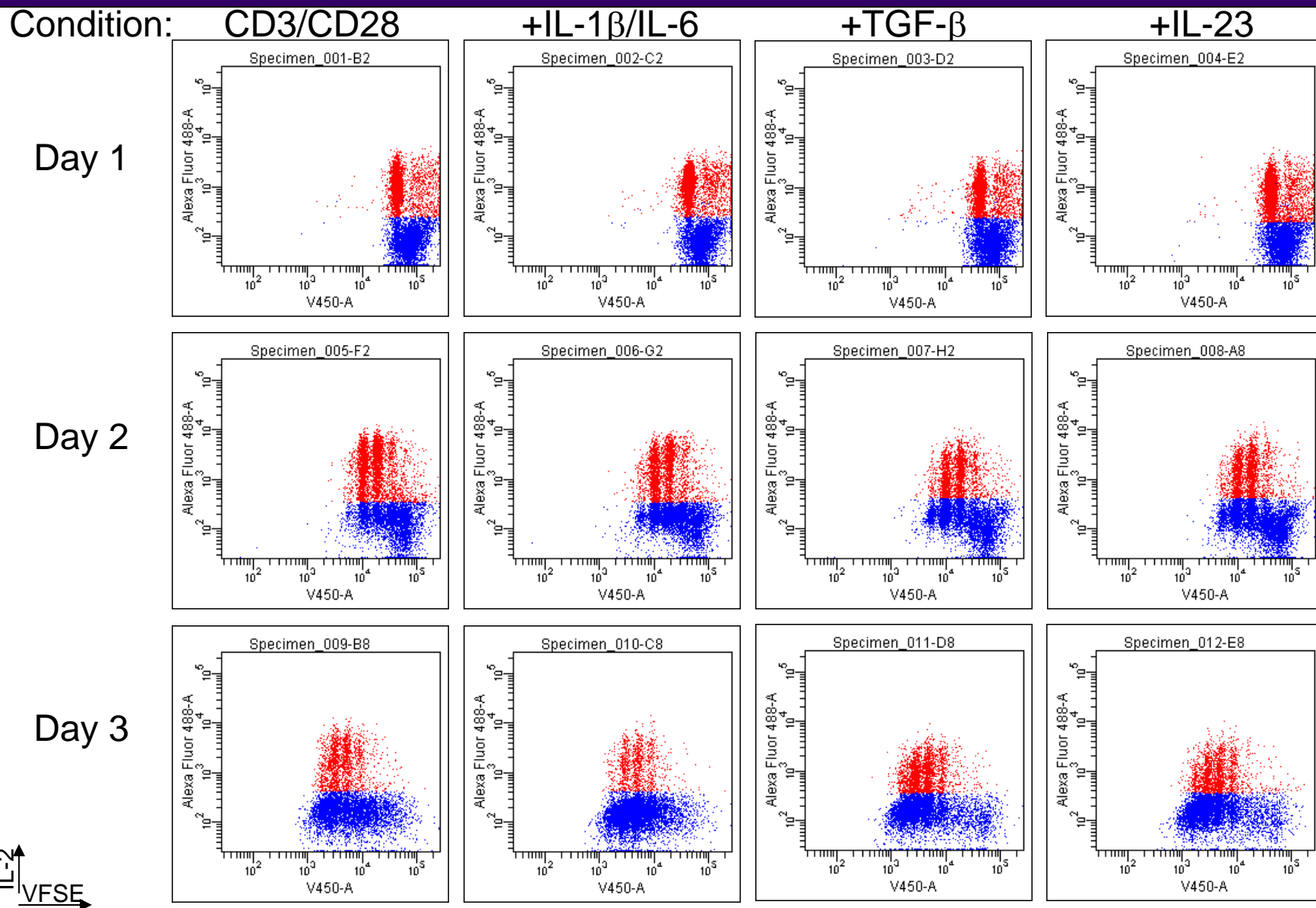


Day 3

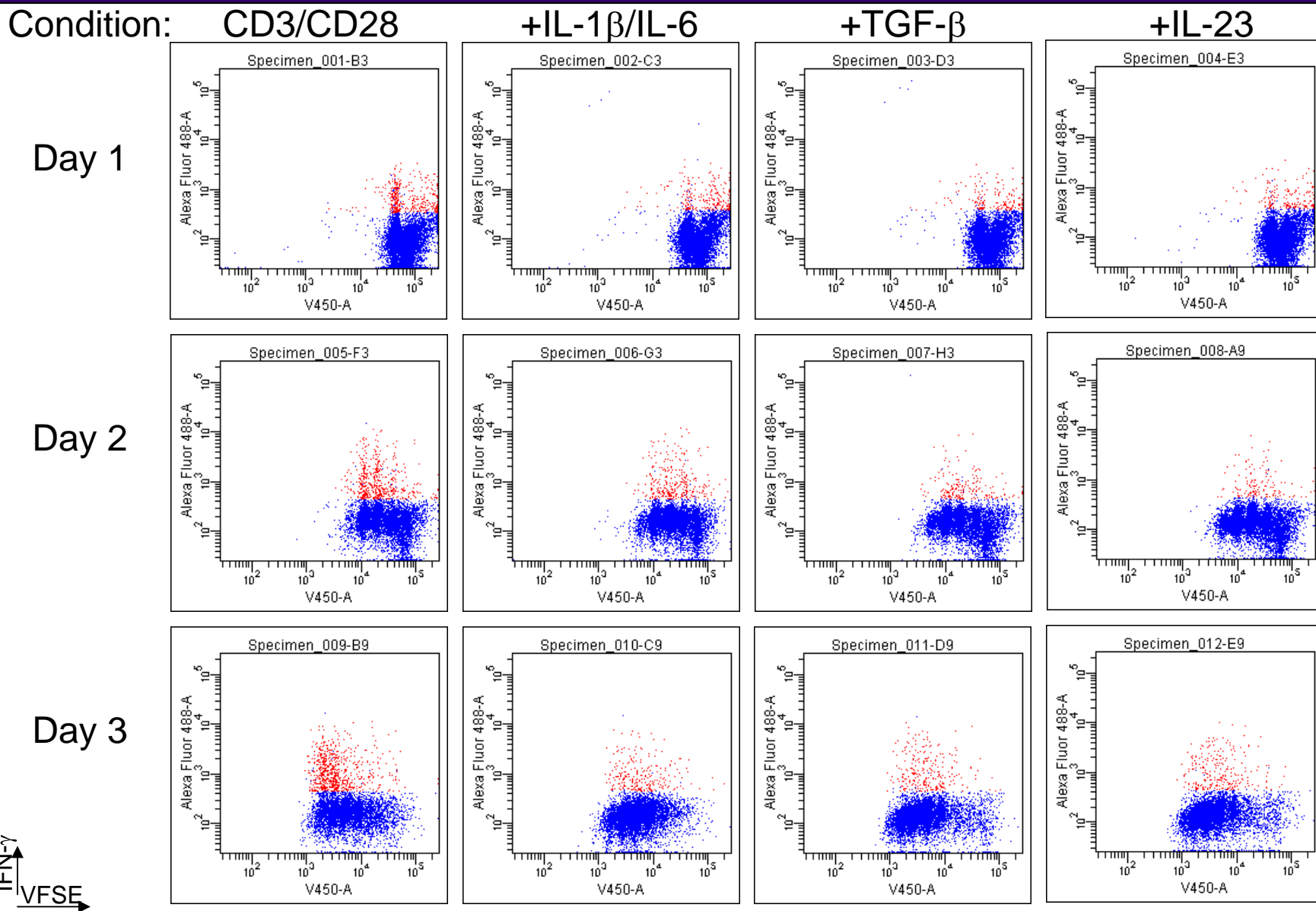


VFSE

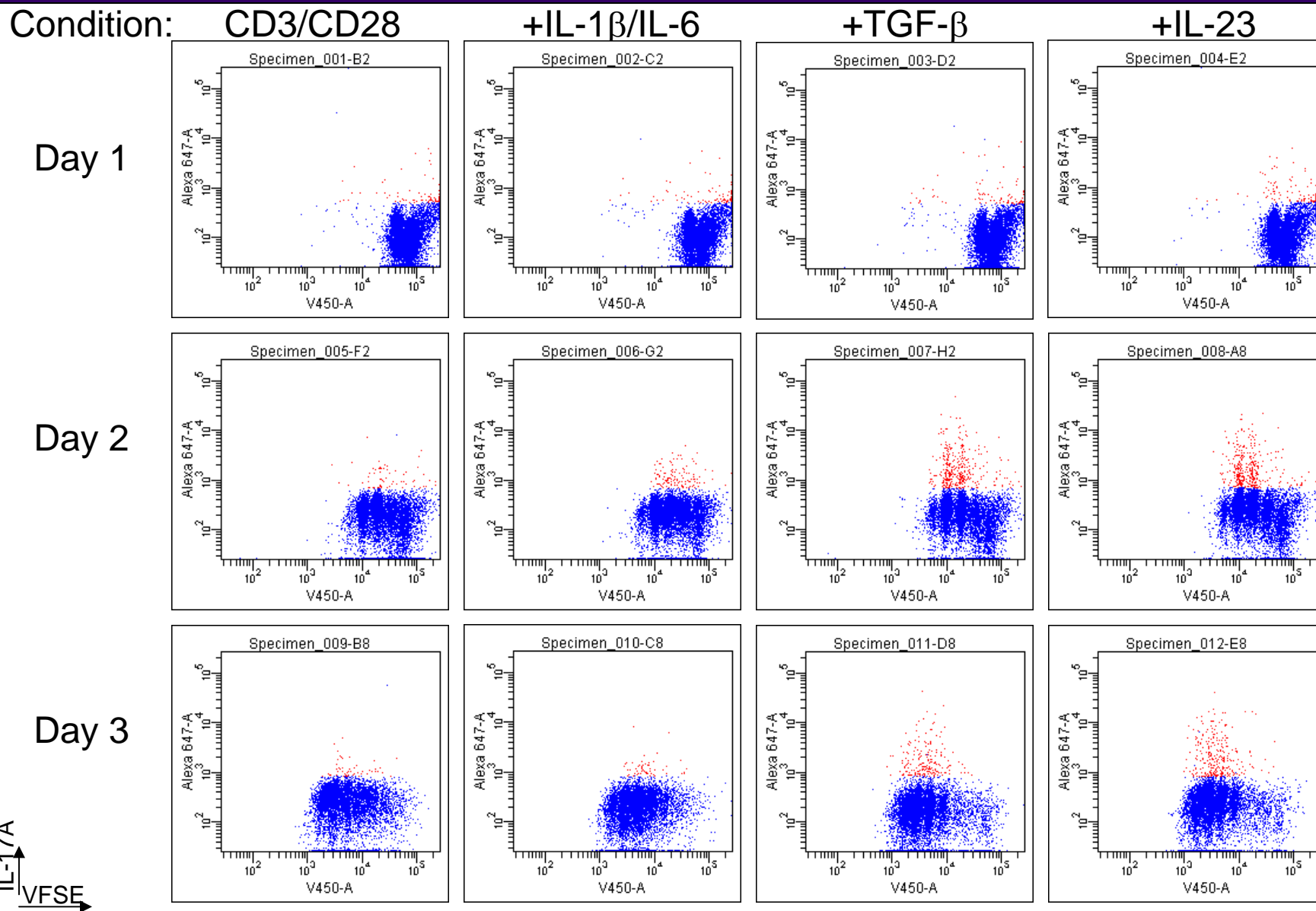
VFSE vs IL-2 data



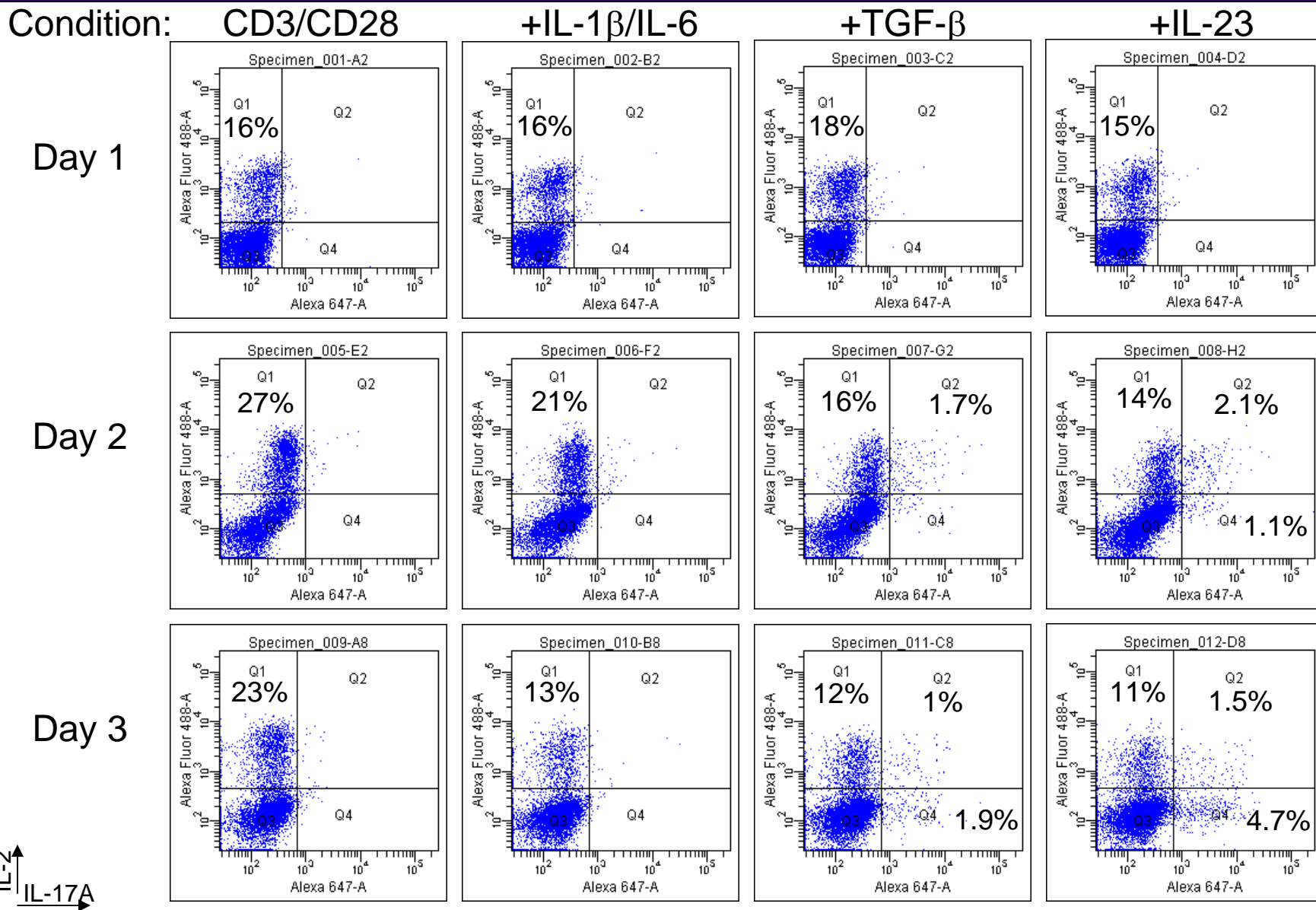
VFSE vs IFN- γ data



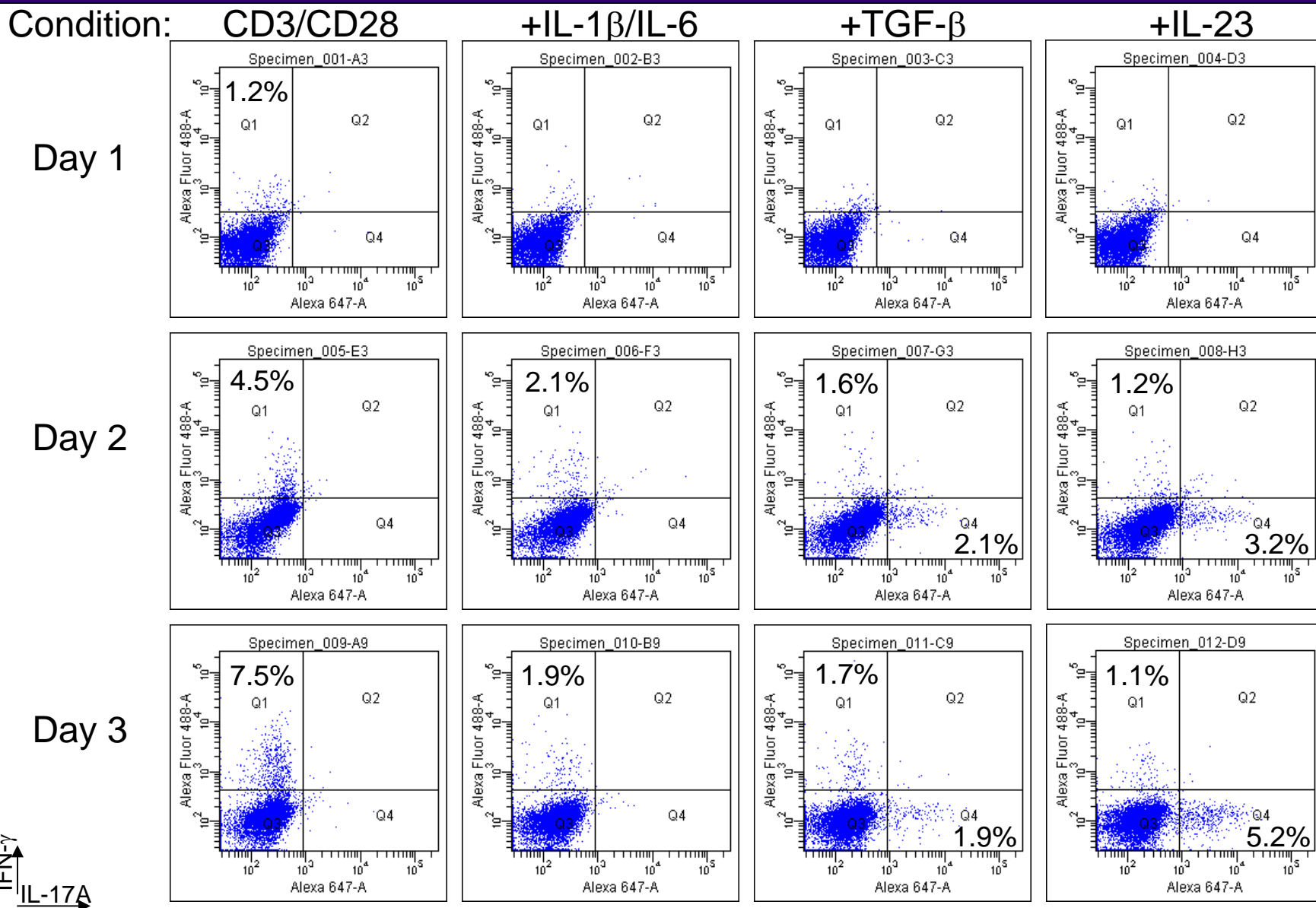
VFSE vs IL-17A data



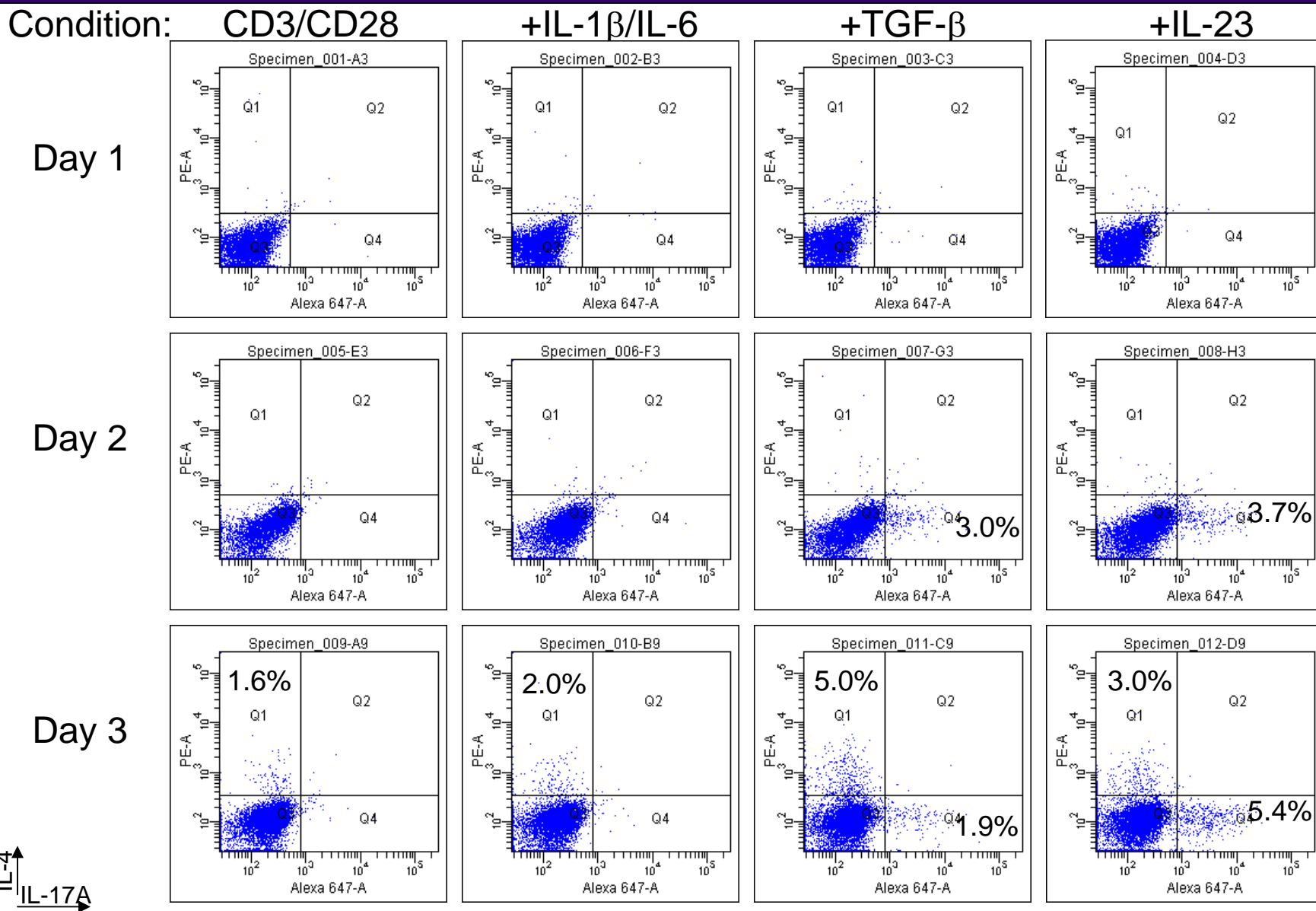
Co-expression of IL-17A vs IL-2



Co-expression of IL-17A vs IFN- γ

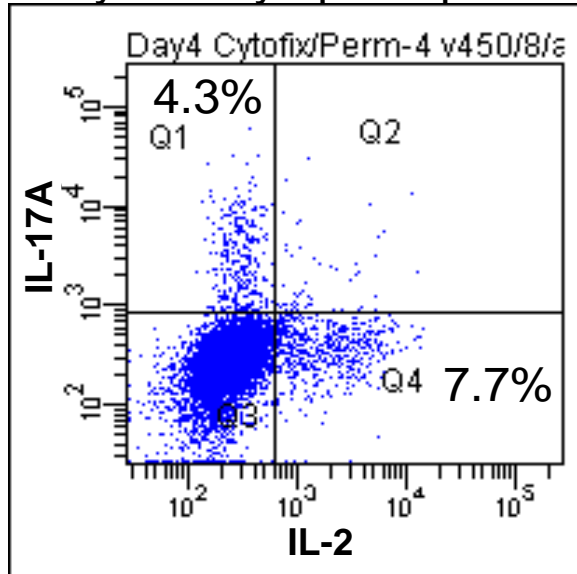


Co-expression of IL-17A vs IL-4

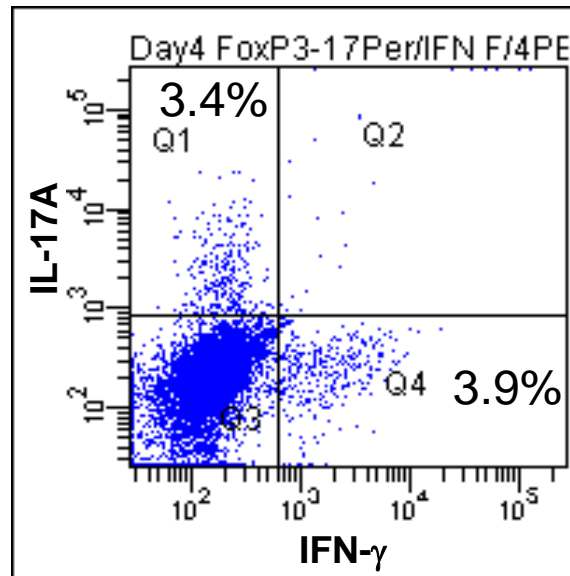
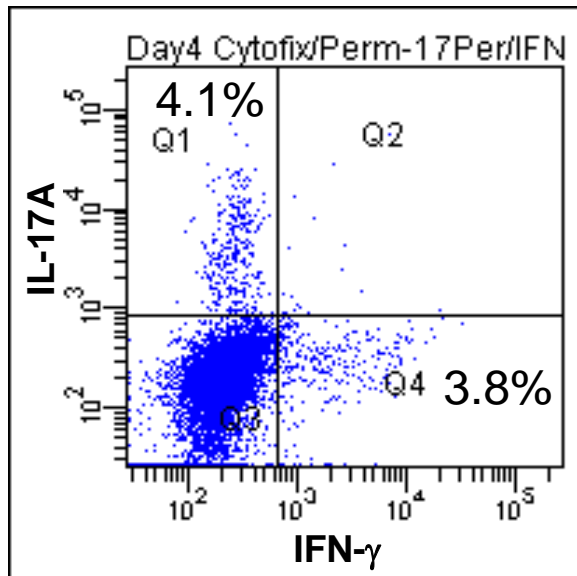
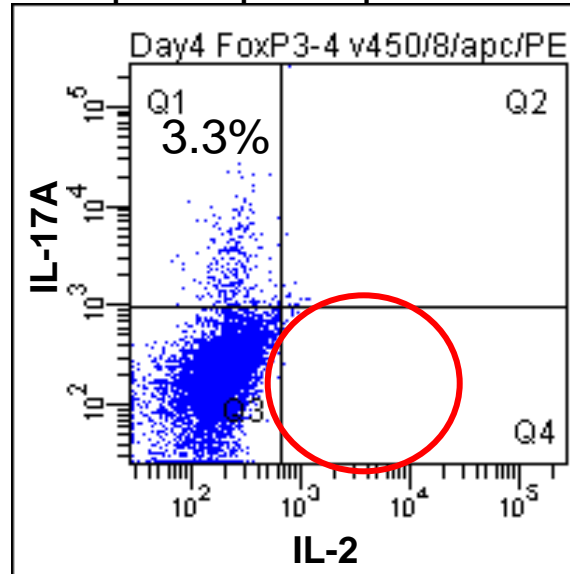


Comparison of two fix/perm protocols

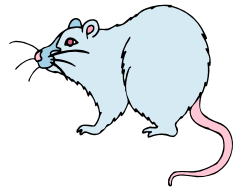
BD Cytfix/Cytoperm protocol



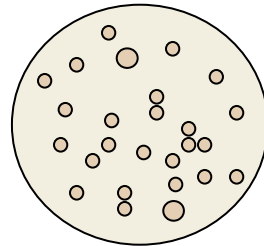
Foxp3 fix/perm protocol



Experimental setup



Harvest
Spleen



CD4 cells enriched
by panning

Cells loaded with
VFSE and washed

anti-CD3e+
anti-CD28

anti-CD3e+
anti-CD28+IL-1 β +IL-6

anti-CD3e+
anti-CD28+IL-1 β +IL-6+ TGF- β

anti-CD3e+
anti-CD28+IL-1 β +IL-6+ TGF- β + IL-23

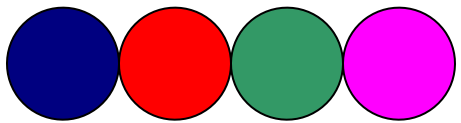
BD Cytofix/Cytoperm™ buffer
Cytokines

+ Monensin

Foxp3 Fix/Perm buffer
Foxp3 and some
cytokines

Harvest
Stimulate with PMA and
Ionomycin for 4–5 hours
+ or – Monensin

- Monensin

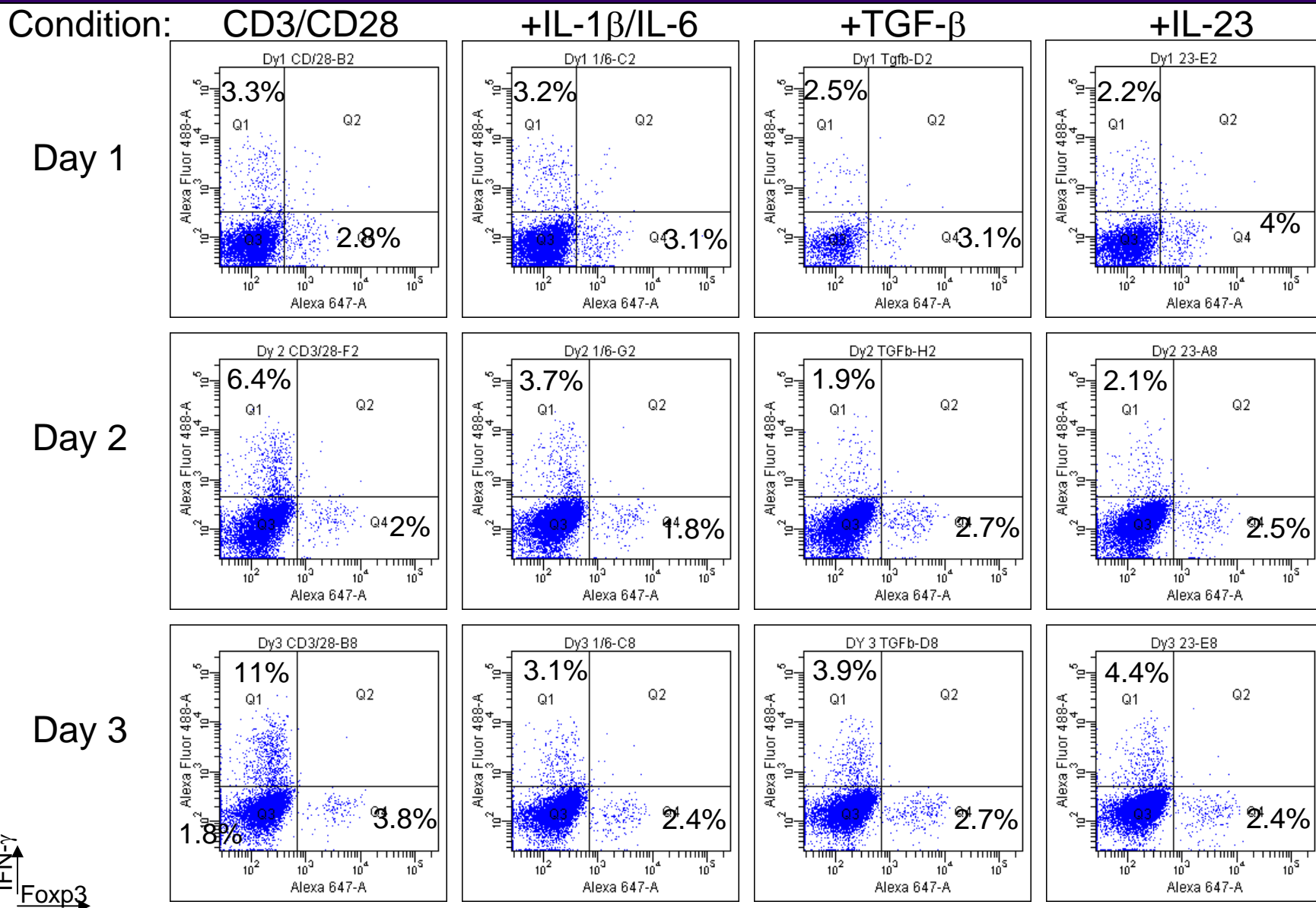


Day 1, 2, 3, 4

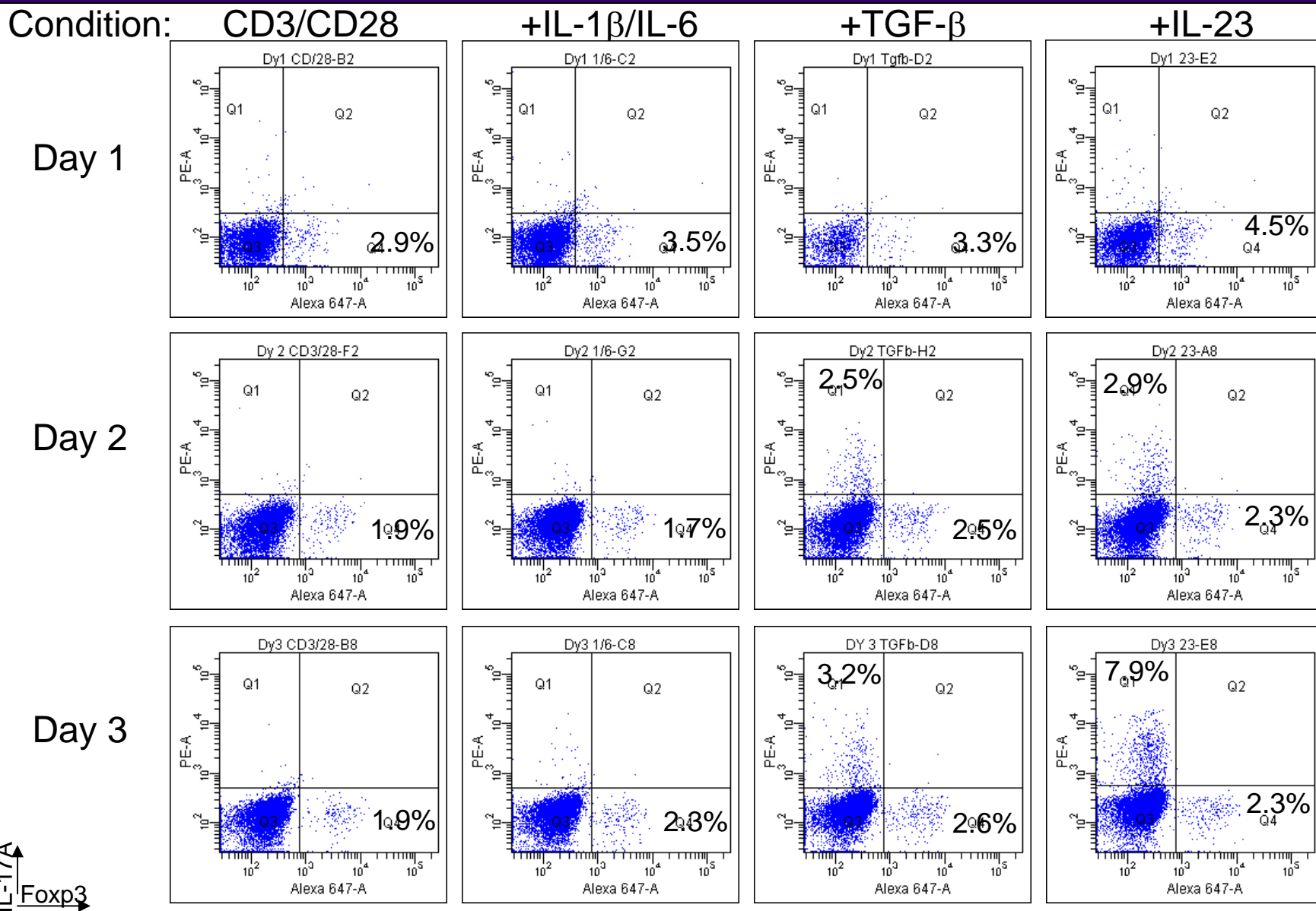
Supernatants: Analysis with BD™ CBA Flex Sets
Cells: BD™ Phosflow Fix and Perm buffer



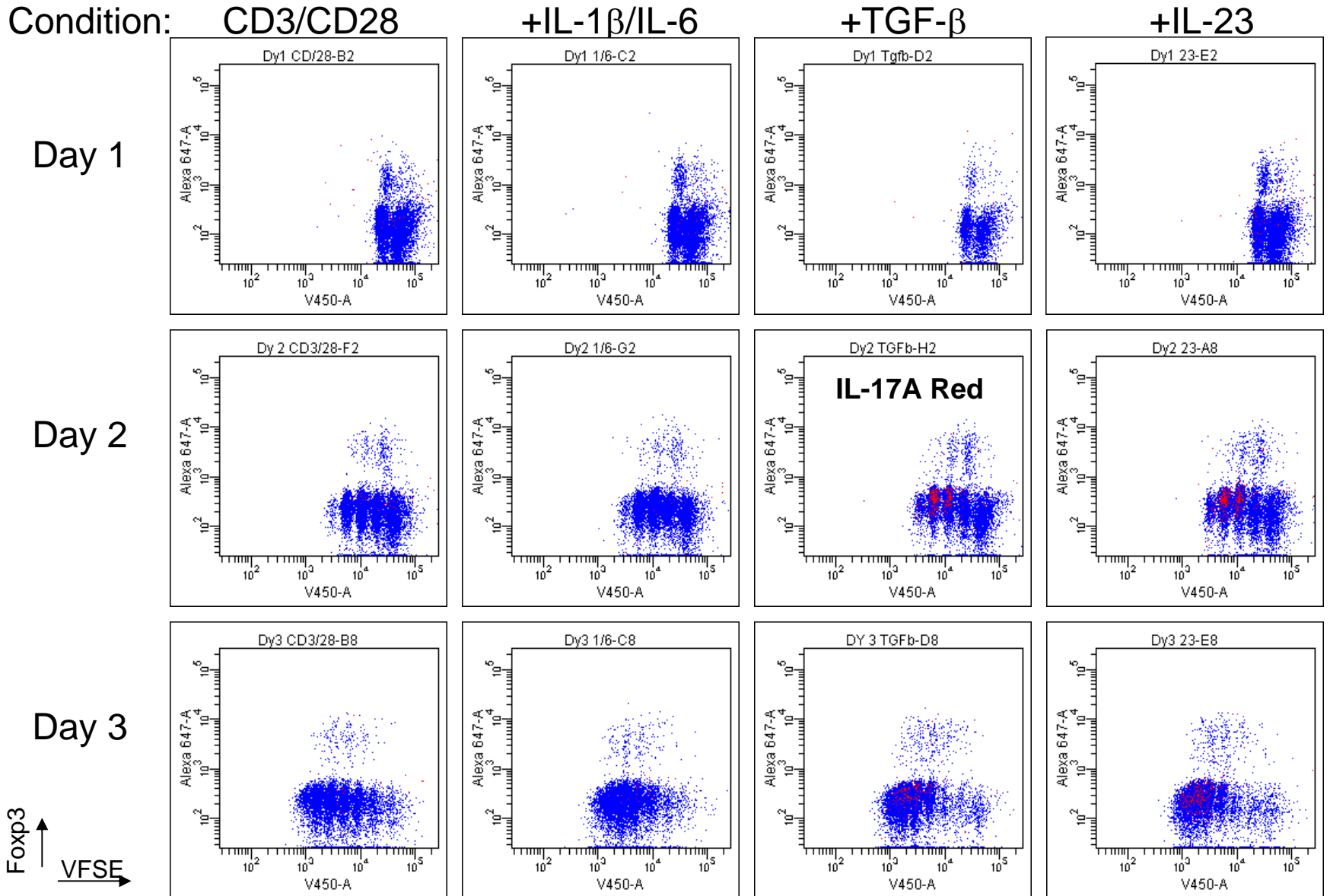
Co-expression of *Foxp3* vs *IFN- γ*



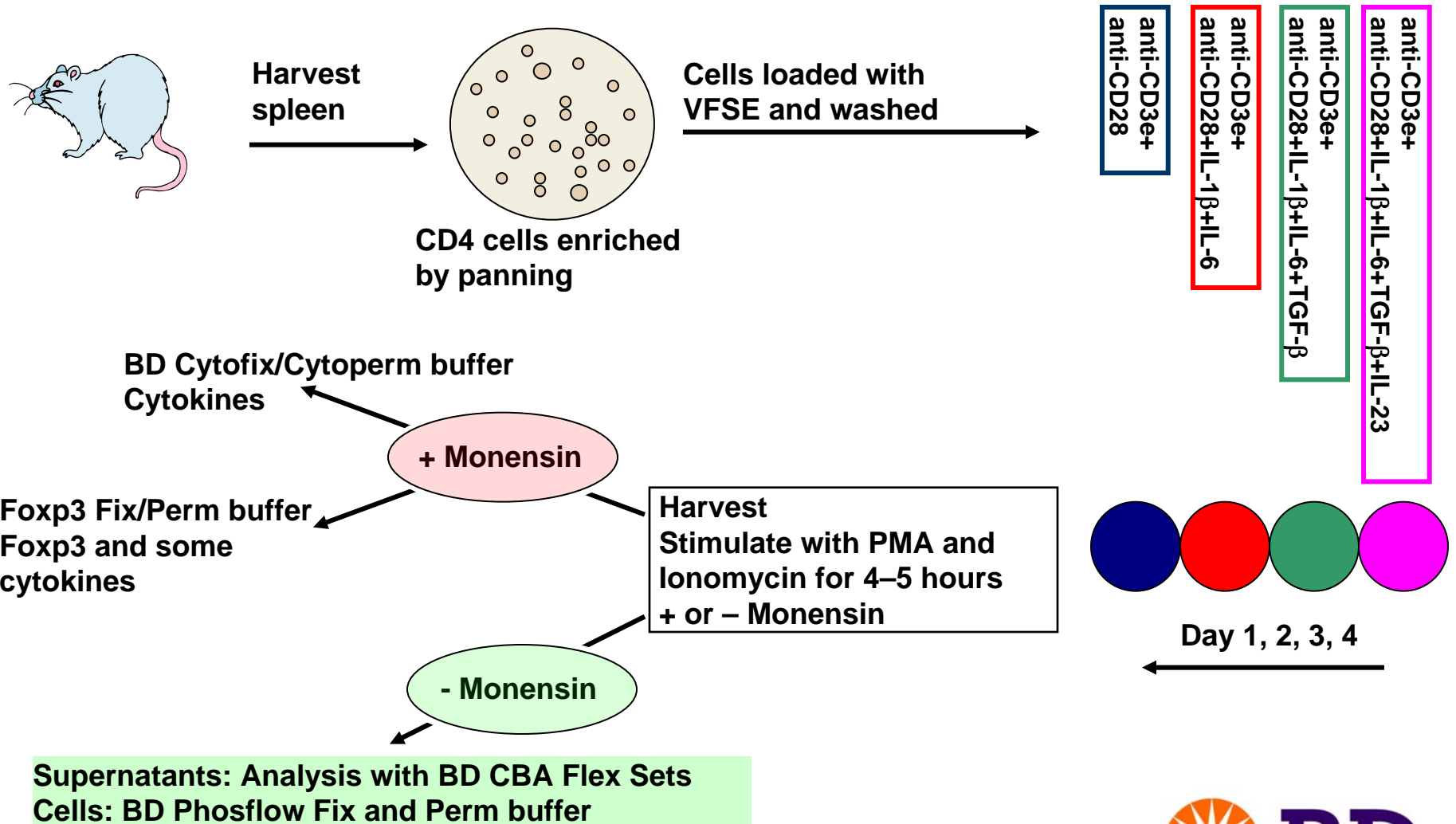
Co-expression of Foxp3 vs IL-17A



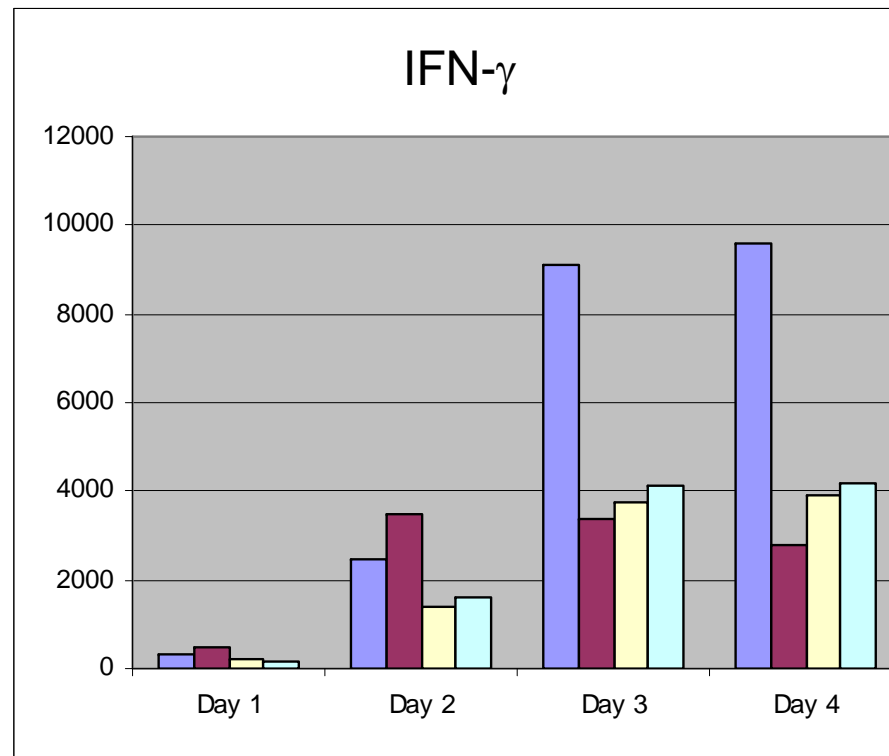
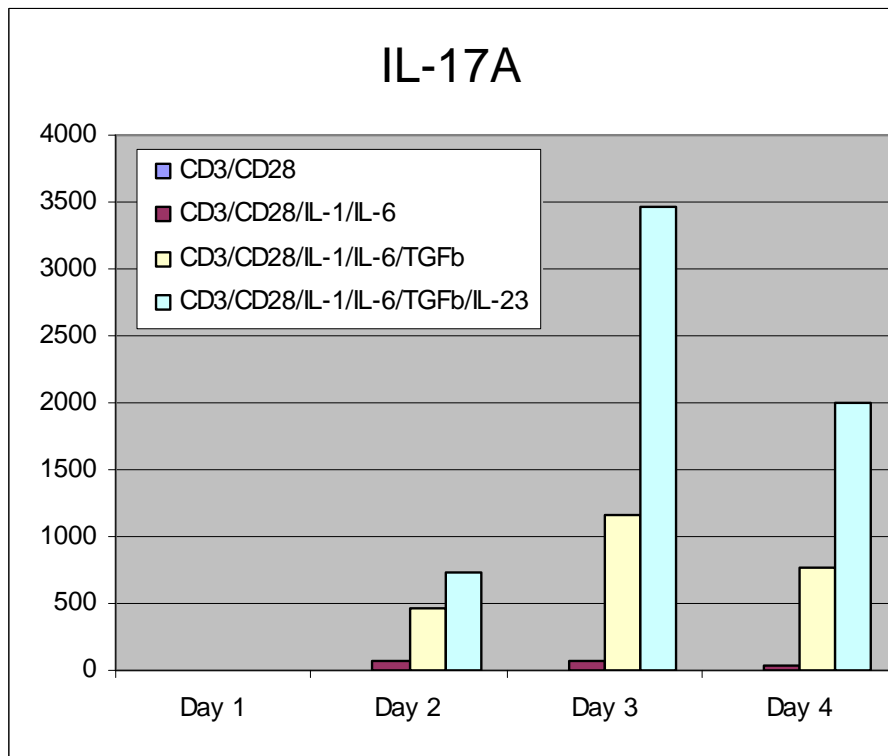
Proliferation of Treg and Th17 cells



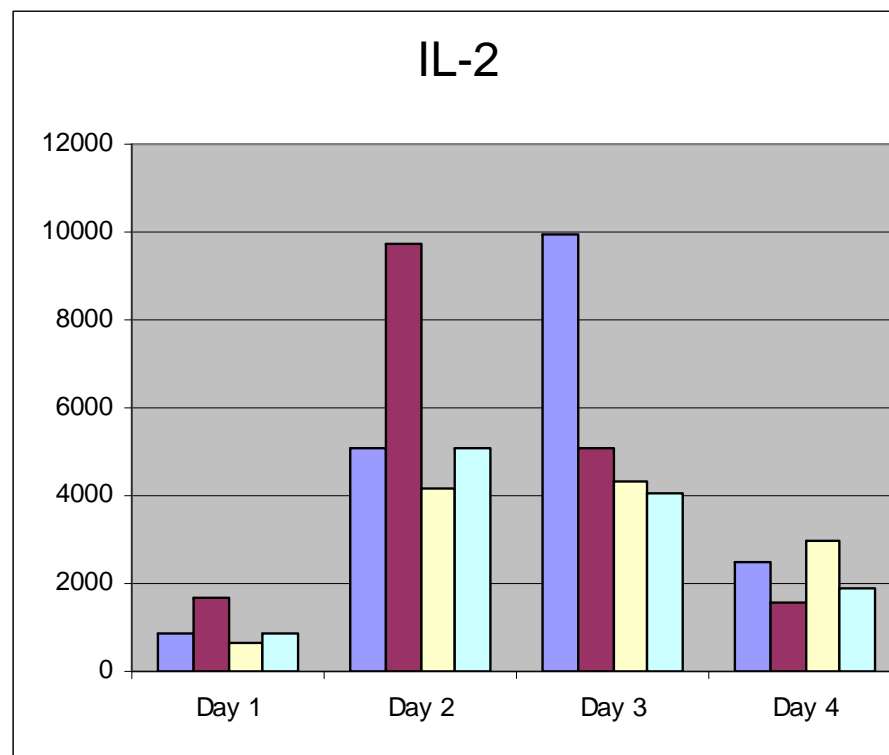
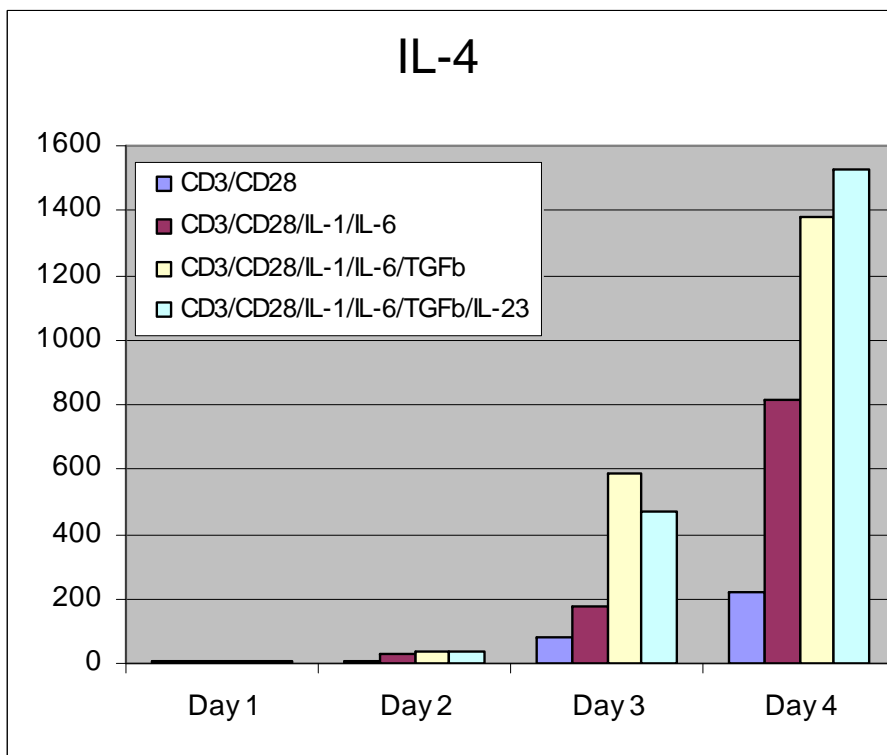
Experimental setup



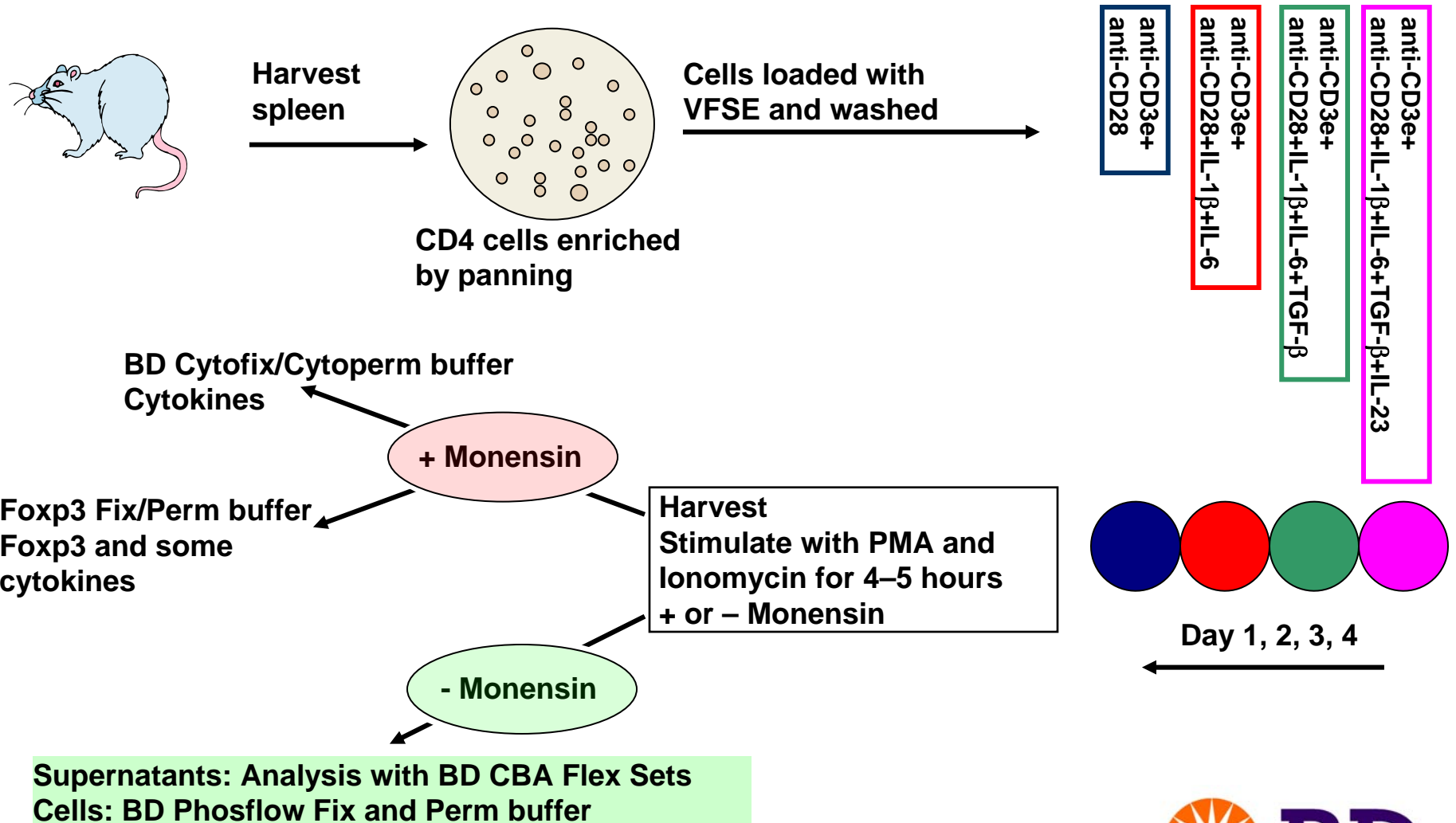
IL-17A and IFN- γ production



IL-4 and IL-2 production



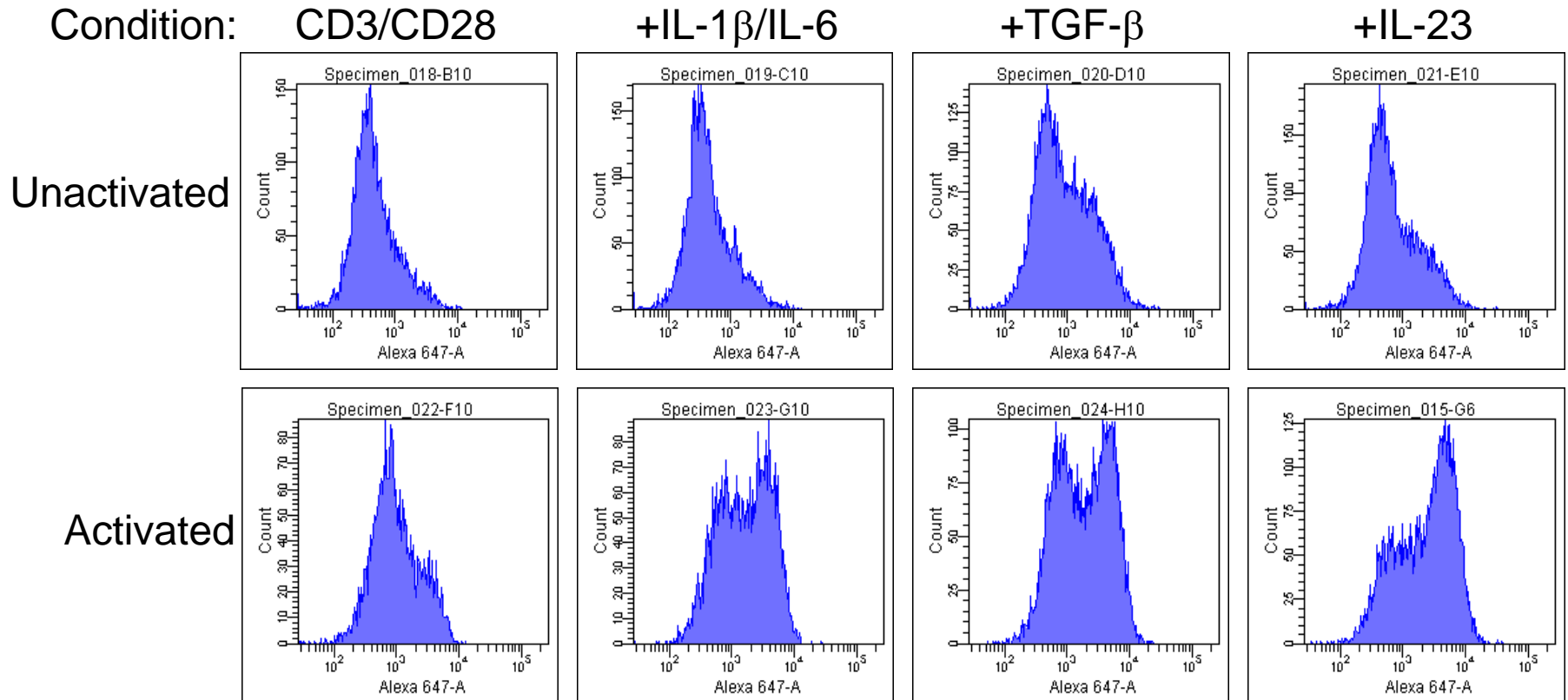
Experimental setup



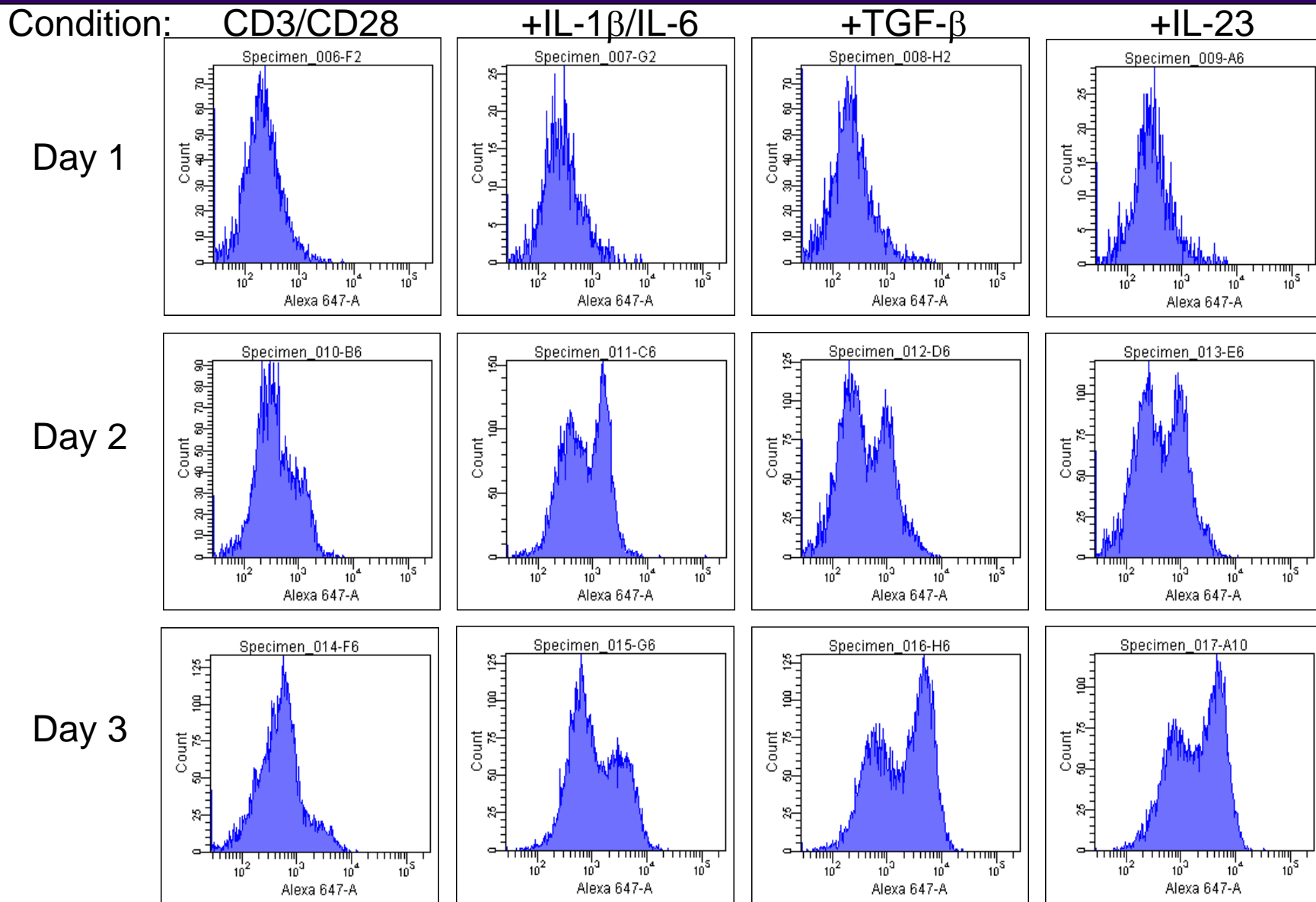
pStat5 detection on day 4

Unactivated: Cells were cultured, harvested, and stained with phosphospecific Stat5 antibody.

Activated: Cells were cultured and activated with PMA/Ionomycin for 5 hours and then stained with phosphospecific Stat5 antibody.



pStat5 in activated cells over time



pStat5 in proliferating cells

Condition:

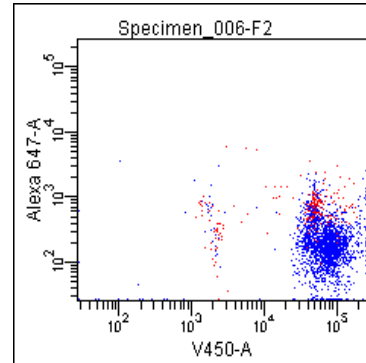
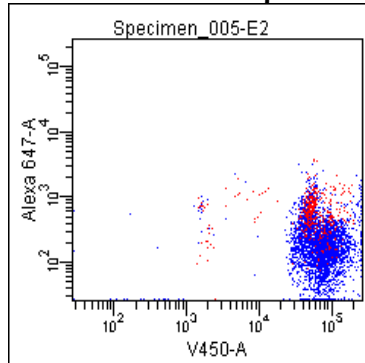
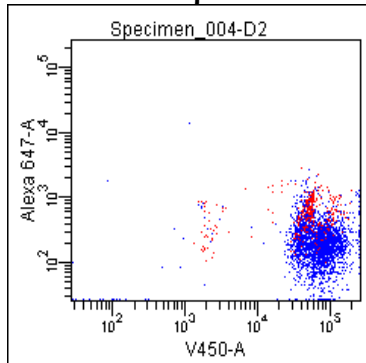
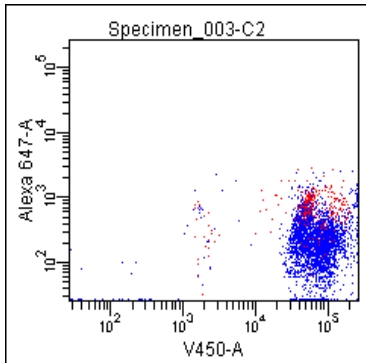
CD3/CD28

+IL-1 β /IL-6

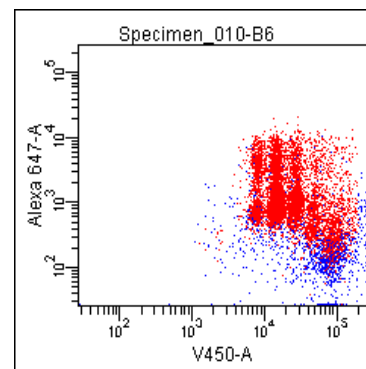
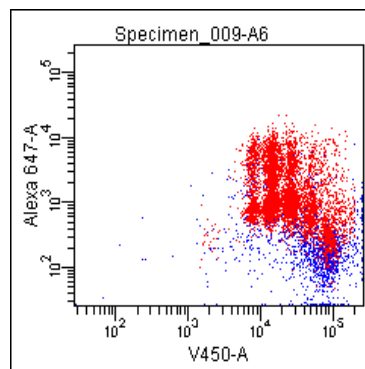
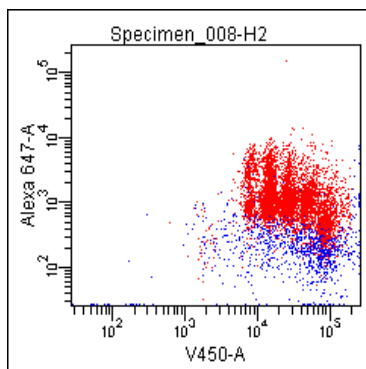
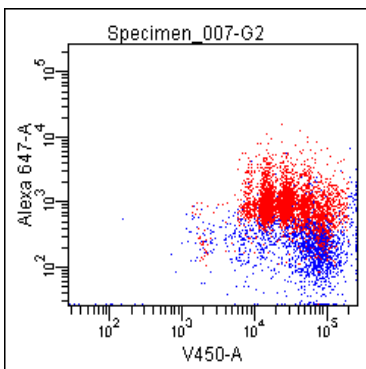
+TGF- β

+IL-23

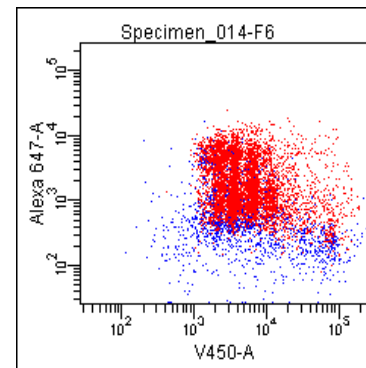
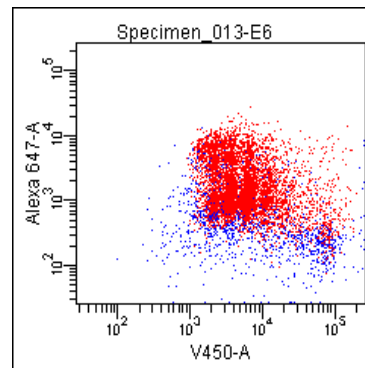
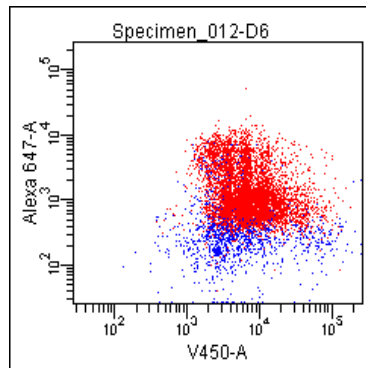
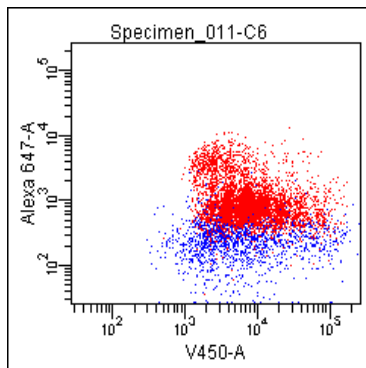
Day 1



Day 2



Day 3



pStat5
VFSE

Conclusions

- Cells proliferated equally well under all four polarization conditions
- Invitro cultures showed that TGF- β was important for polarization of CD4 cells towards Th17
- Initial cultures show co-expression of IL-2 and IL-17 that later become independent of each other
- Detection of secreted cytokines (by CBA) correlated with the intracellular staining
- Cytokine production by proliferating cells resulted in increased phosphorylation of the signal transducer Stat5



Acknowledgments

- Jeanne Elia
- Xiao-Wei Wu
- Ravi Hingorani

- Jacob Rabenstein

