



Analyzing Stem Cell Populations Using Flow Cytometry

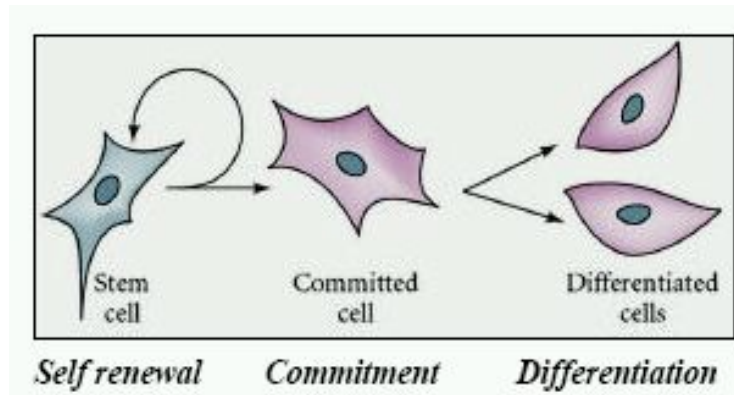
Robert Balderas
VP, Research & Development
BD Biosciences

*For Research Use Only. Not for use in diagnostic or therapeutic procedures.
Class I (1) Laser Products
23-9889-00*

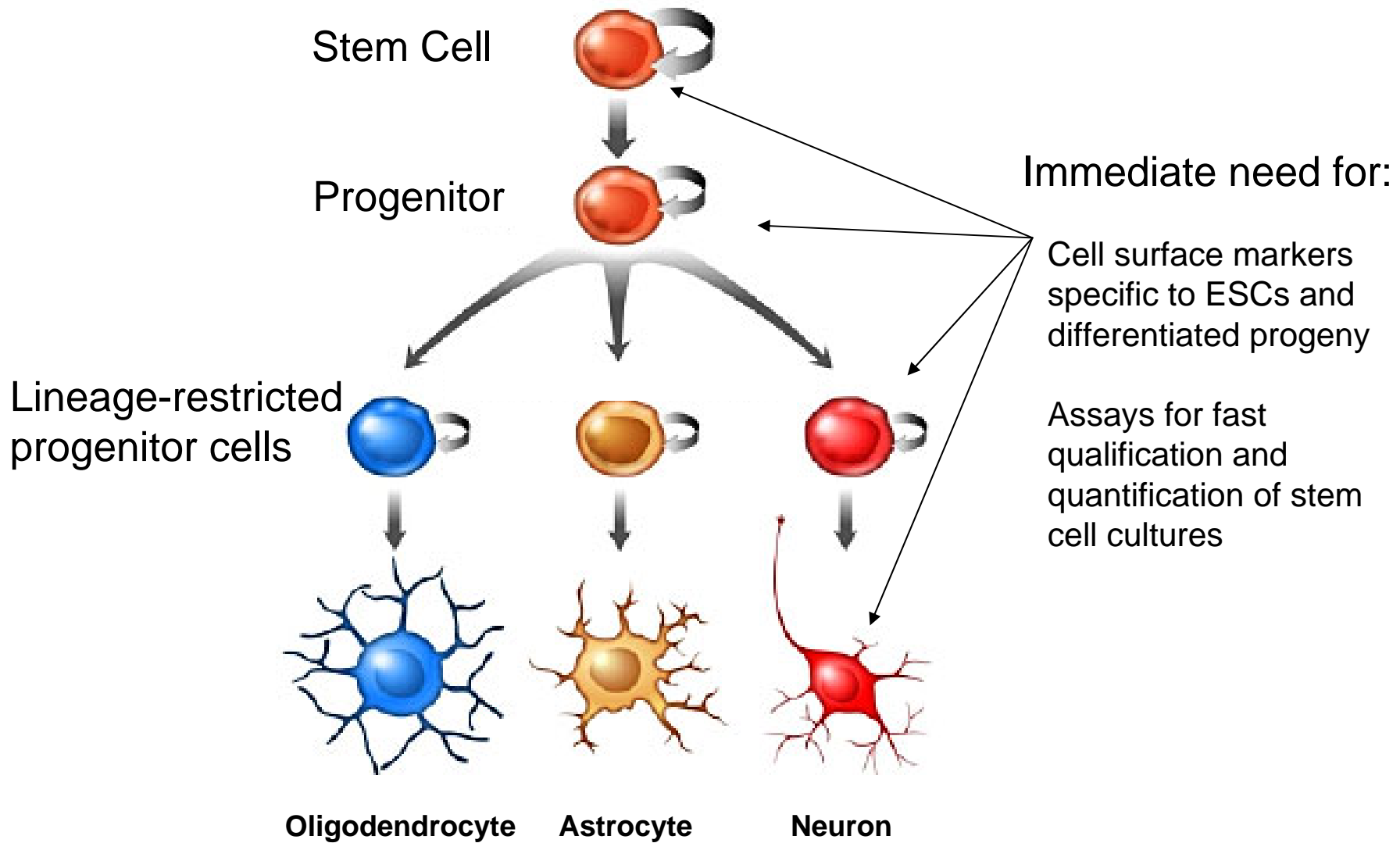
March 27, 2008



Stem Cell Research

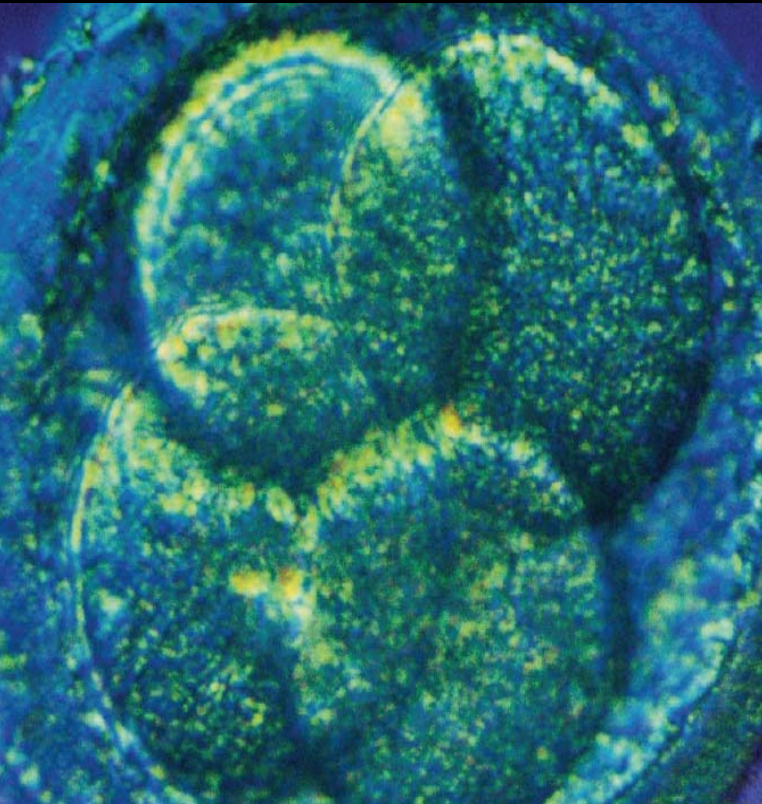


- **What drives cell differentiation in a specific direction?**
 - How are stem cells programmed?
 - What are the environmental factors?
- **What are the characteristic features of a cell in a particular state of differentiation?**
 - Which biomarkers are expressed?
 - What is the function behind expression patterns?
- **What are the required factors for keeping a cell in a particular differentiation state?**



Stem Cell Subpopulations

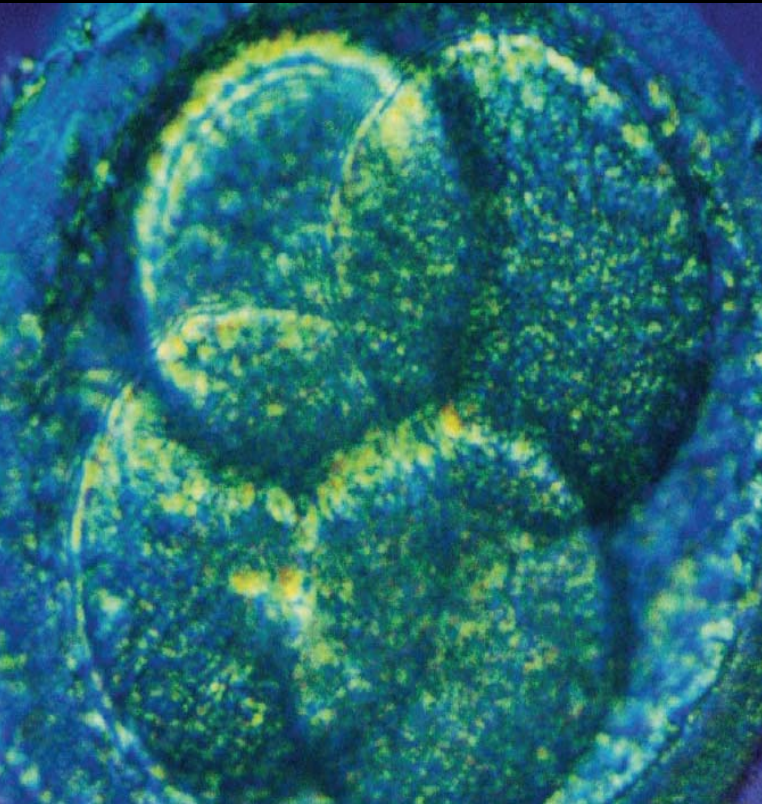
Determining How Stem Cells Differentiate



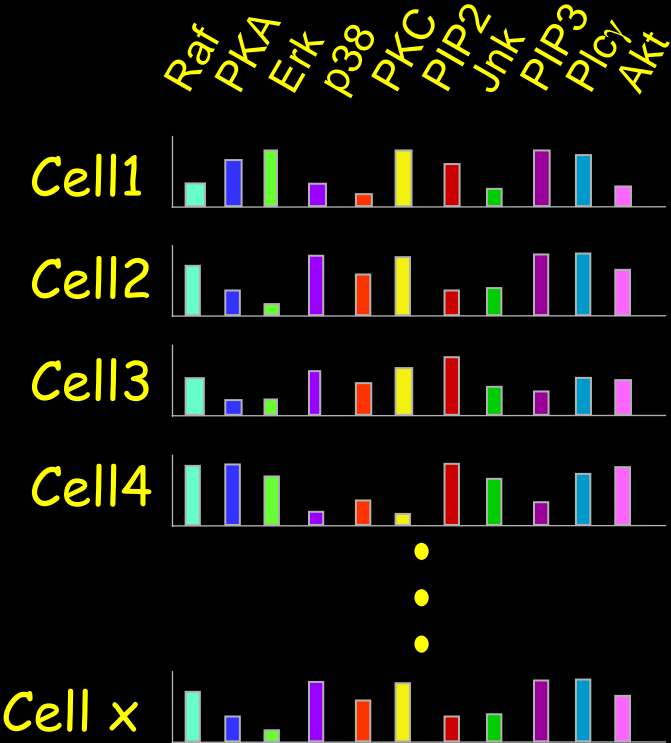
- Cells are computational devices
- Their inner workings (algorithms) cannot be determined by reading their steady-state values
- Cells must be interrogated to determine relationships of signaling components

Stem Cell Subpopulations

Determining How Stem Cells Differentiate

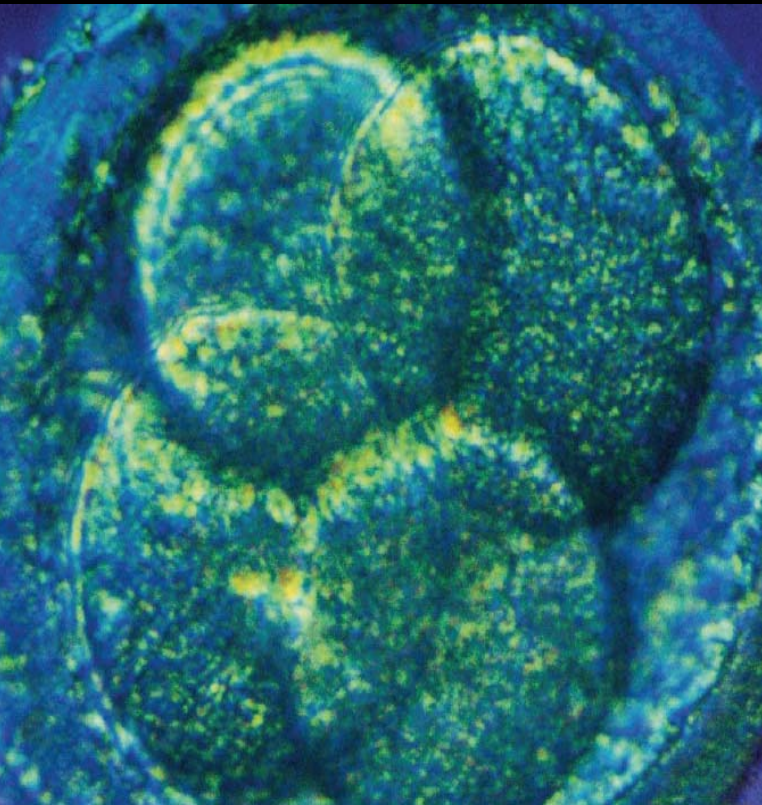


Flow
cytometry



Stem Cell Subpopulations

Determining How Stem Cells Differentiate




Computational
analysis

- Quantitative nature of flow data allows powerful statistical regimes to *derive* signaling maps.
- These inner relationships relate to
 - Differentiation state
 - Response to a given growth factor
 - Biosignatures and cell response profiles

Program Overview

- Gating strategies for analyzing stem cell populations
- Pluripotent cell analysis
- Neuronal progenitor cells
- Phosphorylation analysis





Gating Strategies for Stem Cell Populations Using Flow Cytometry

*For Research Use Only. Not for use in diagnostic or therapeutic procedures.
Class I (1) Laser Products*

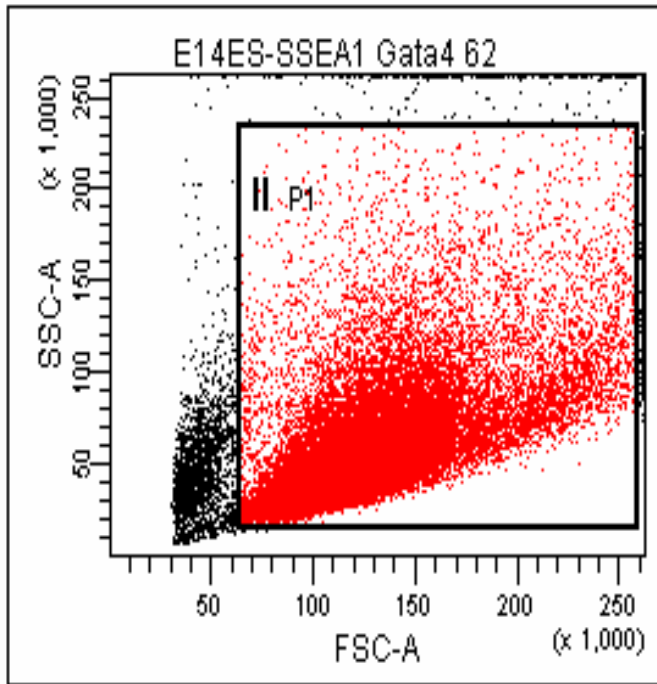


Gating Strategies for Analyzing Stem Cell Flow Data

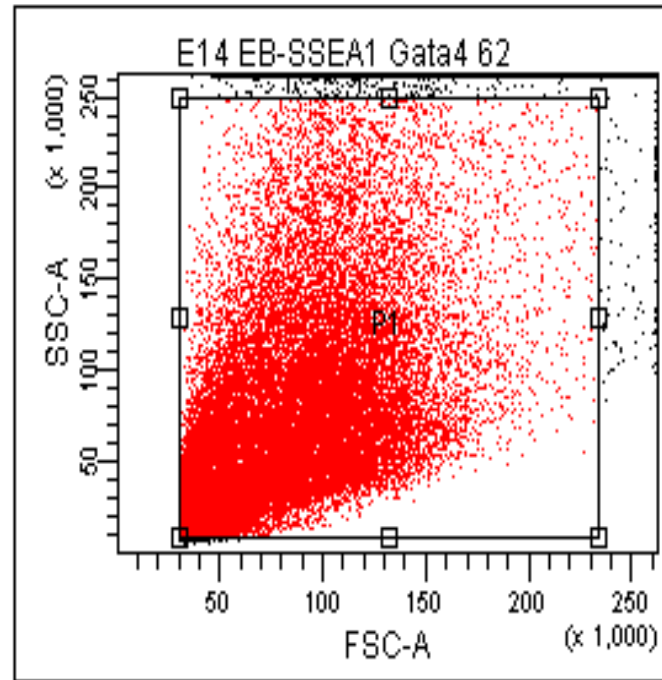
- mES and mEB E14 cell populations were stained and analyzed with:
 - FITC-conjugated SSEA-1
 - PE GATA4
- Populations were gated in different areas of the scatter plot to determine if size and granularity could be used to provide additional information
- Gating strategies can be created for sorting of homogeneous subpopulations



Gating Strategies for Analyzing Stem Cell Flow Data



ES

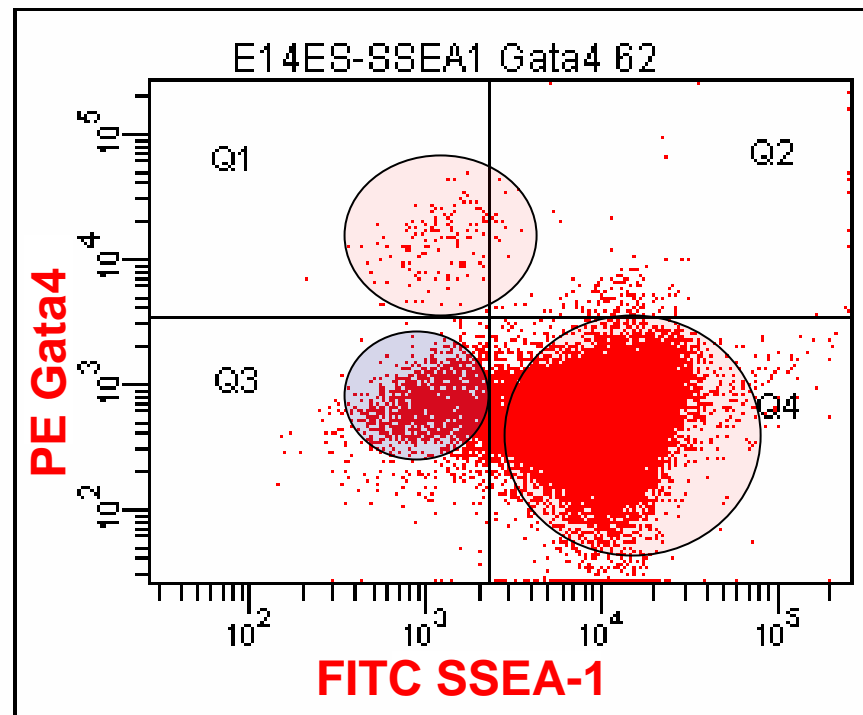
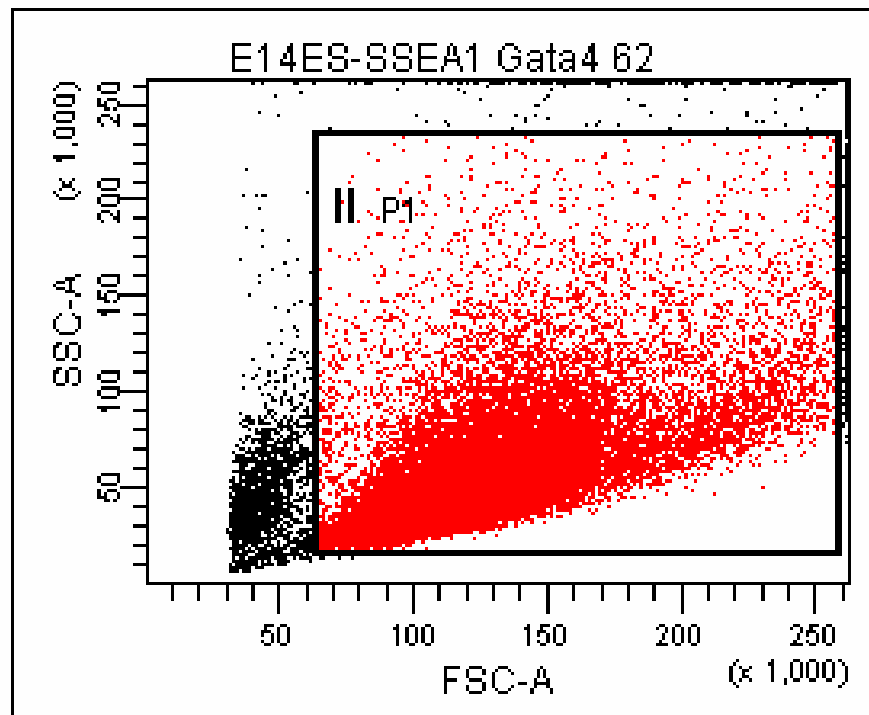


EB

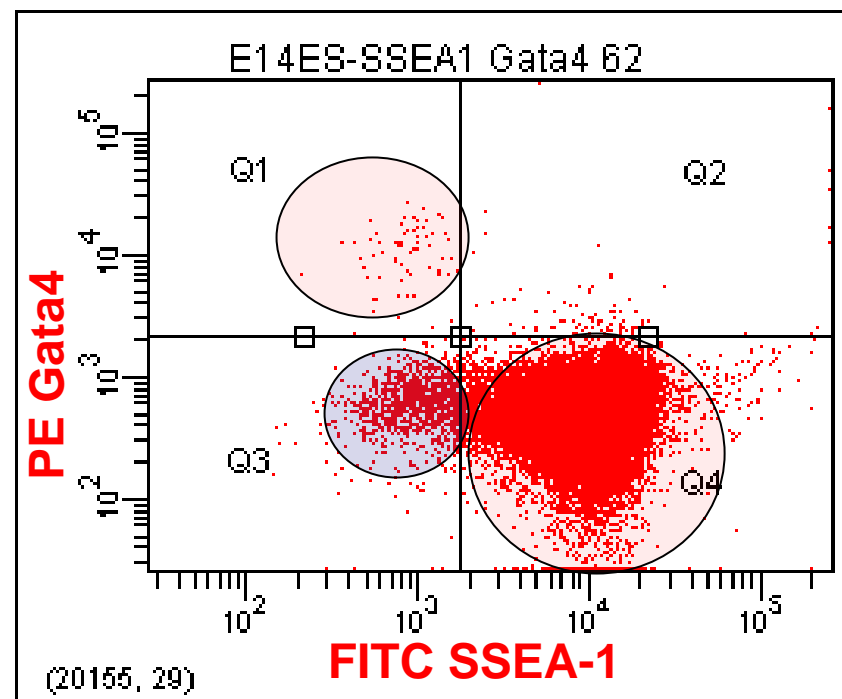
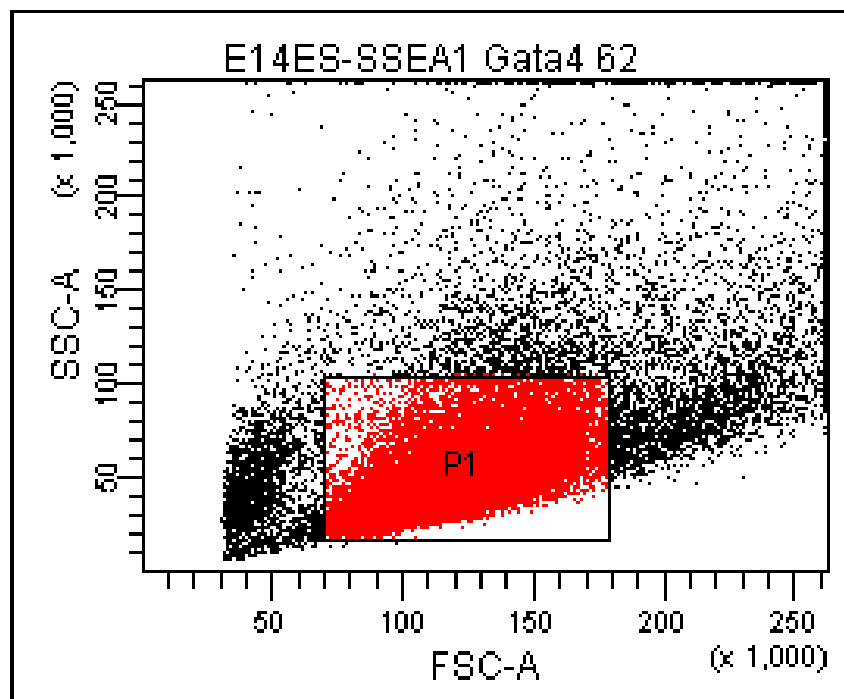
(Forward and side scatter patterns are different)

Can we find sub-populations of cells within each cell profile?

Analysis of SSEA-1 and GATA4 on mES E14

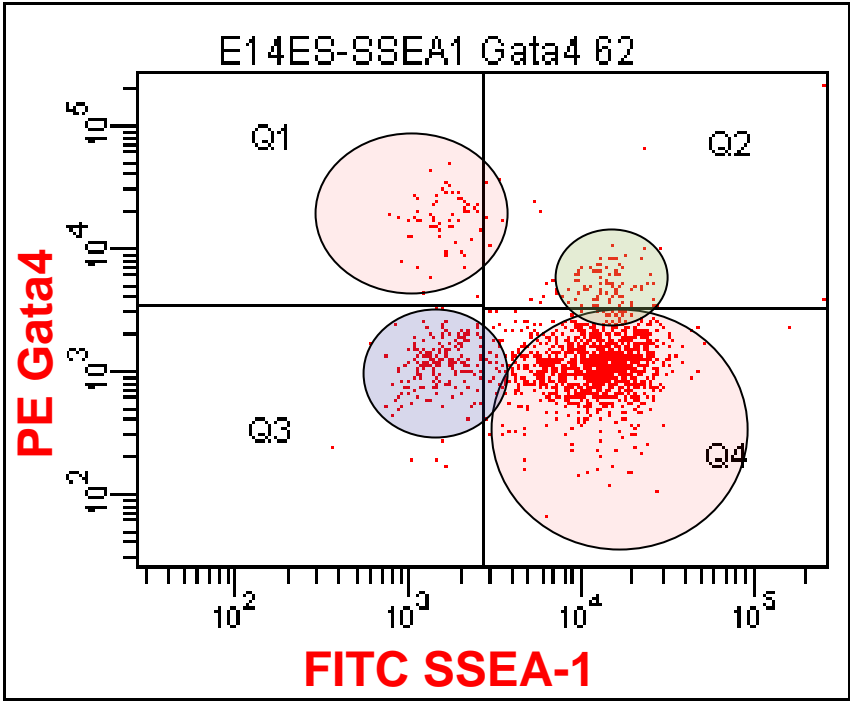
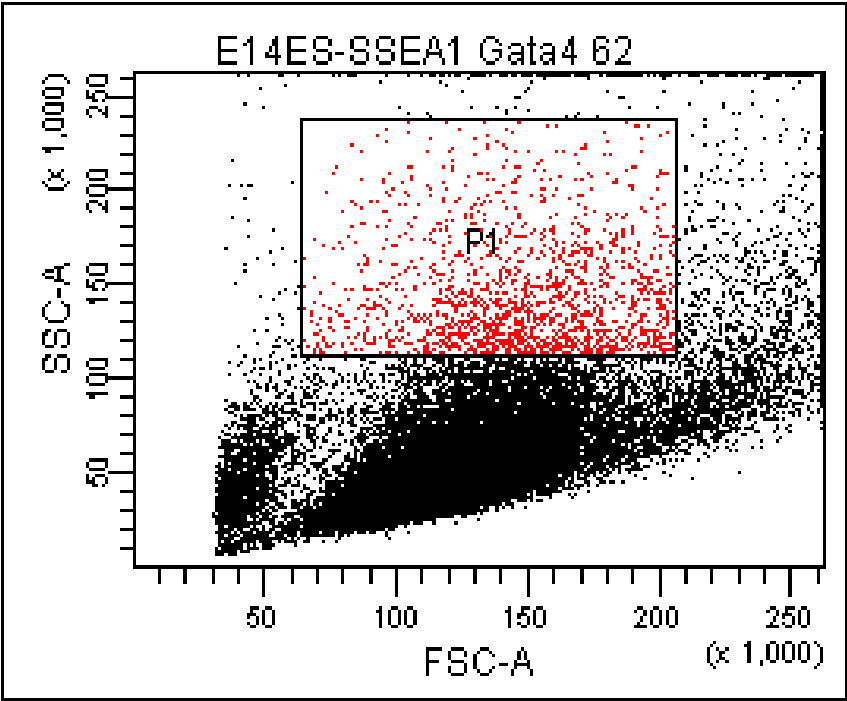


Analysis of SSEA-1 and GATA4 on mES E14

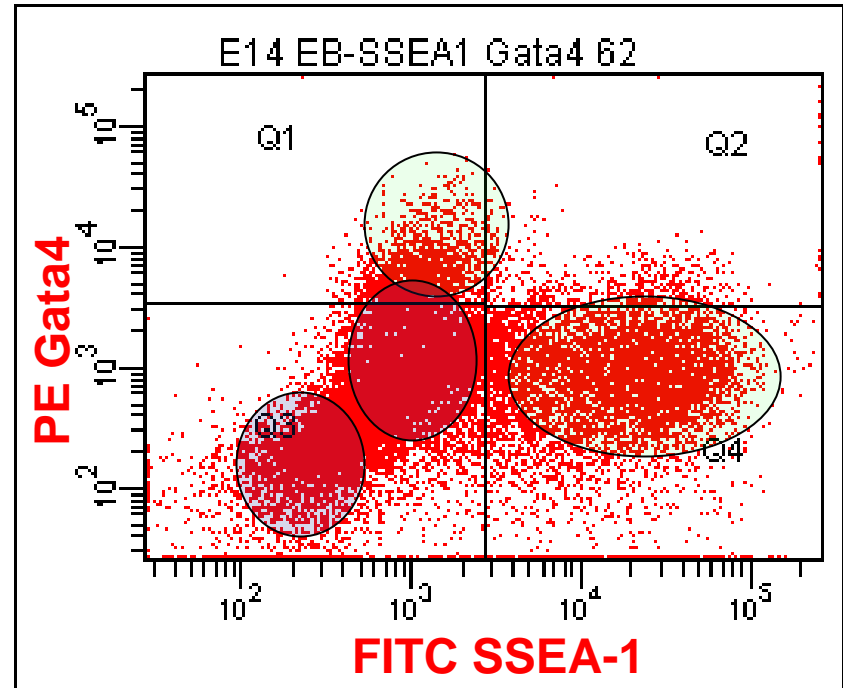
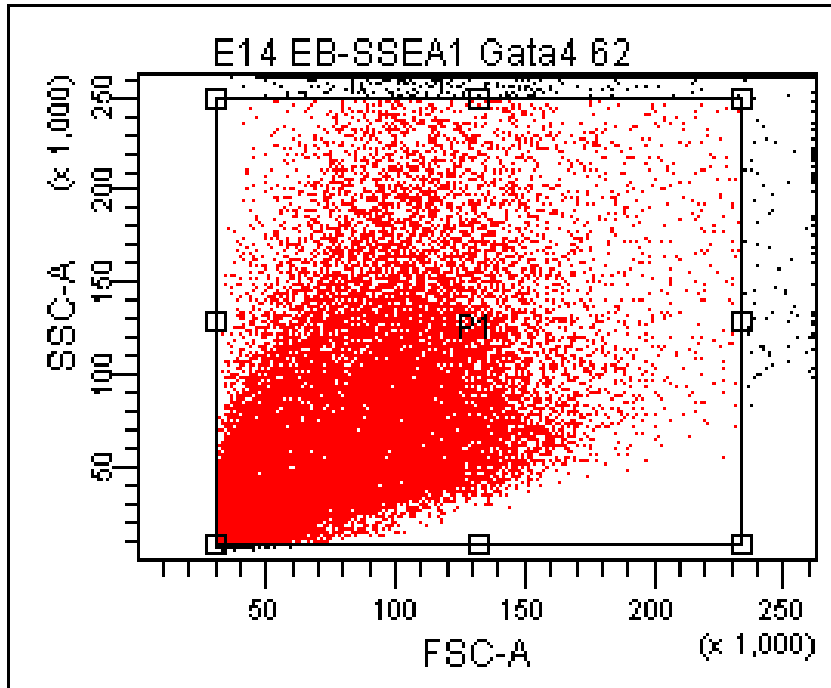


Analysis of SSEA-1 and GATA4 on mES E14

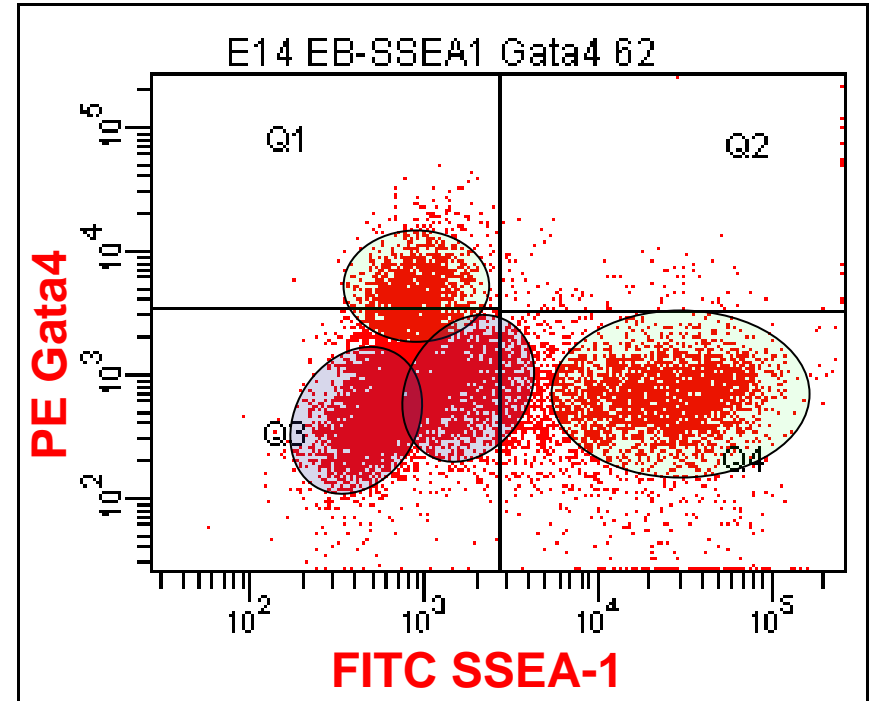
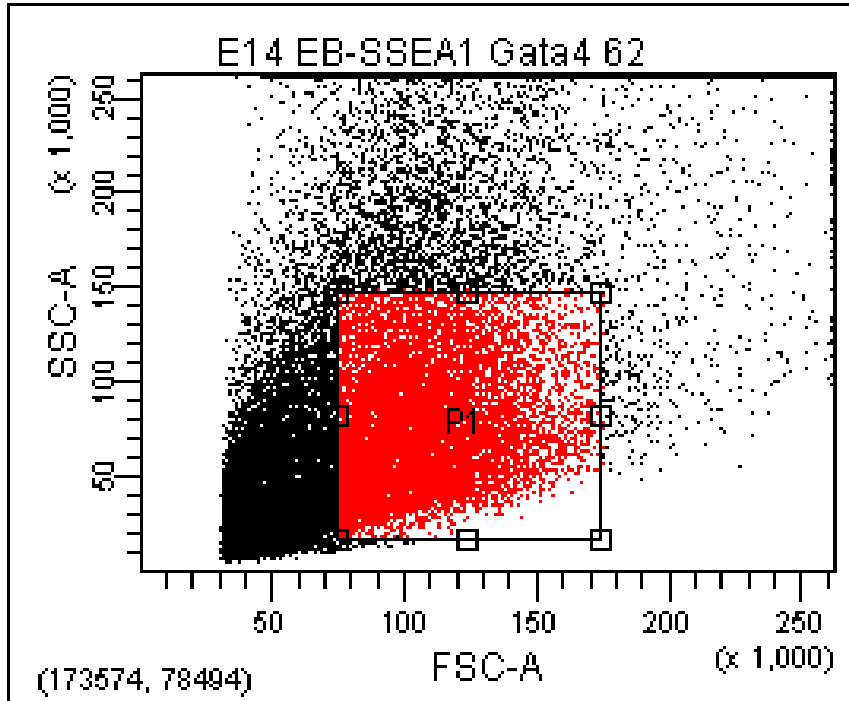
Gating subpopulations reveals SSEA-1-negative and GATA4-positive cells in naïve mES culture



Analysis of SSEA-1 and GATA4 on mES E14-derived Embryoid Bodies

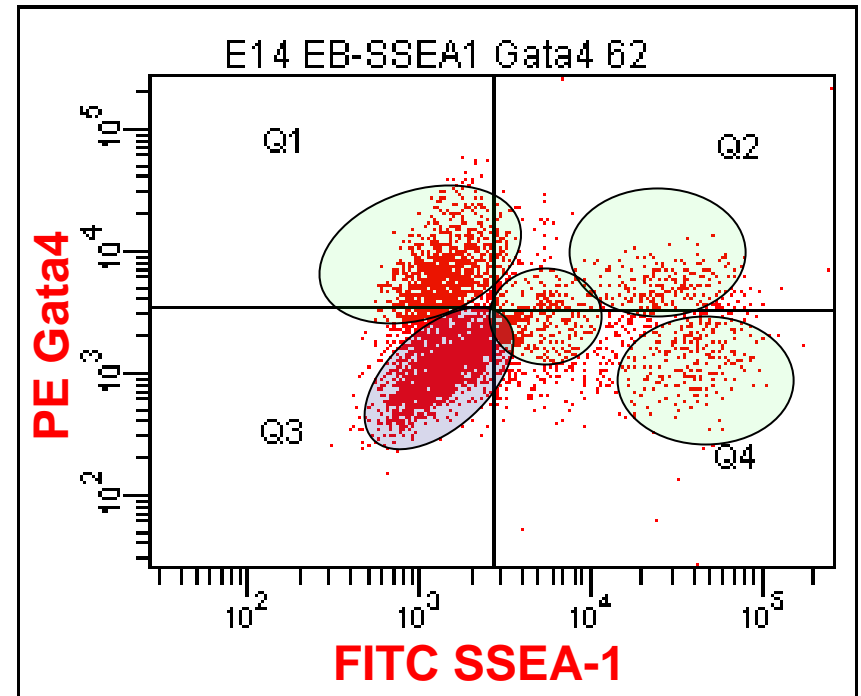
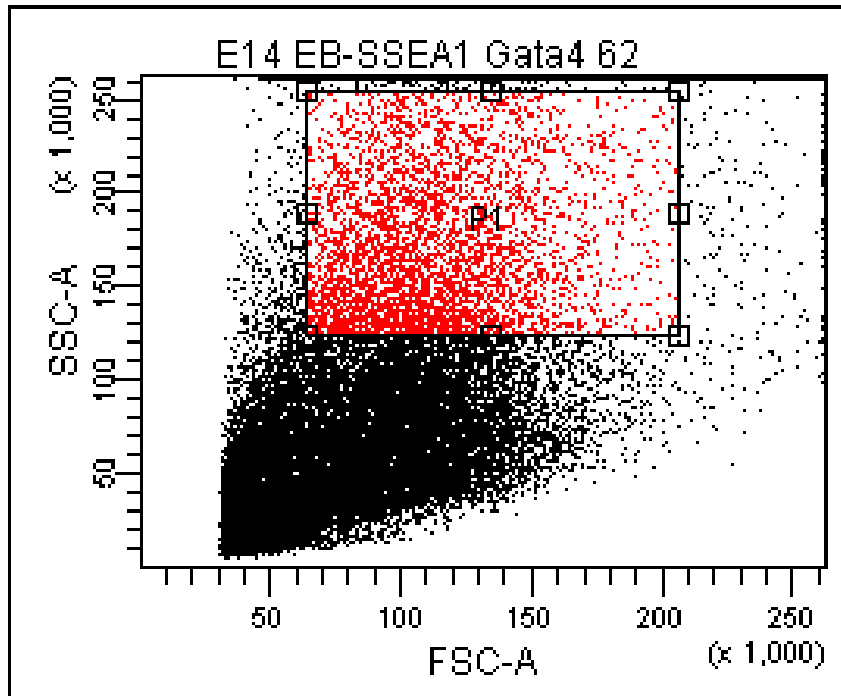


Analysis of SSEA-1 and GATA4 on mES E14-derived Embryoid Bodies



Analysis of SSEA-1 and GATA4 on mES E14-derived Embryoid Bodies

Gating reveals subpopulations of GATA4-positive cells in embryoid bodies





Pluripotent Cell Analysis

*For Research Use Only. Not for use in diagnostic or therapeutic procedures.
Class I (1) Laser Products*

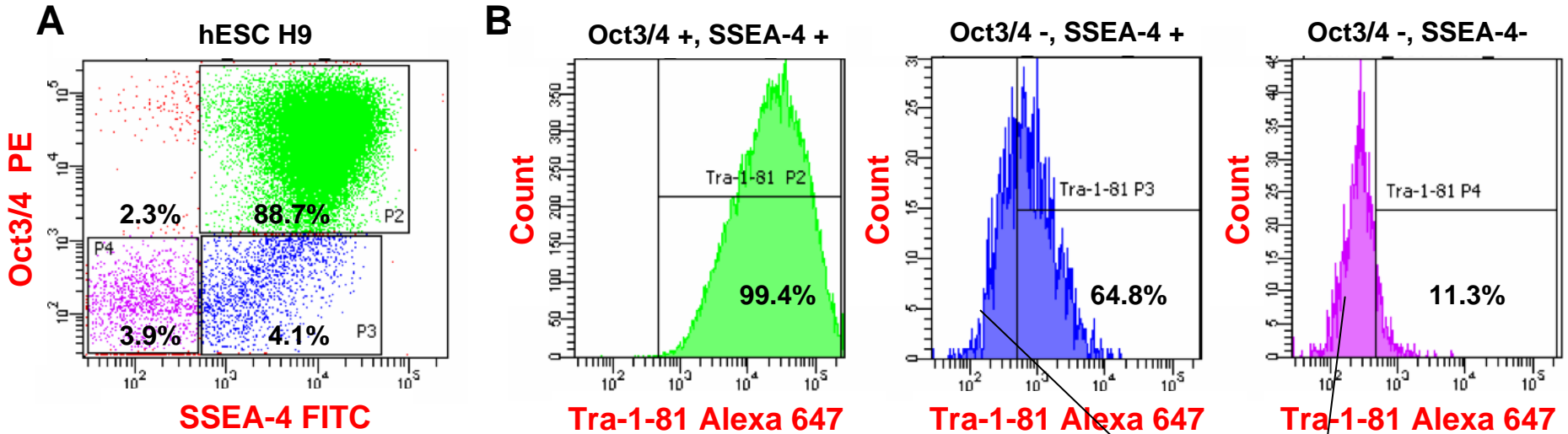


Markers of Pluripotency

Cell surface markers
SSEA-3
SSEA-4
Tra-1-60
Tra-1-81
CD9
CD81
SSEA-1/CD15 (neg)
CD30 (neg/low)

Intracellular markers
Sox2
Oct3/4
Nanog

Three-color flow-cytometric analysis of human embryonic stem cells using fluorochrome-conjugated antibodies to pluripotent stem cell markers

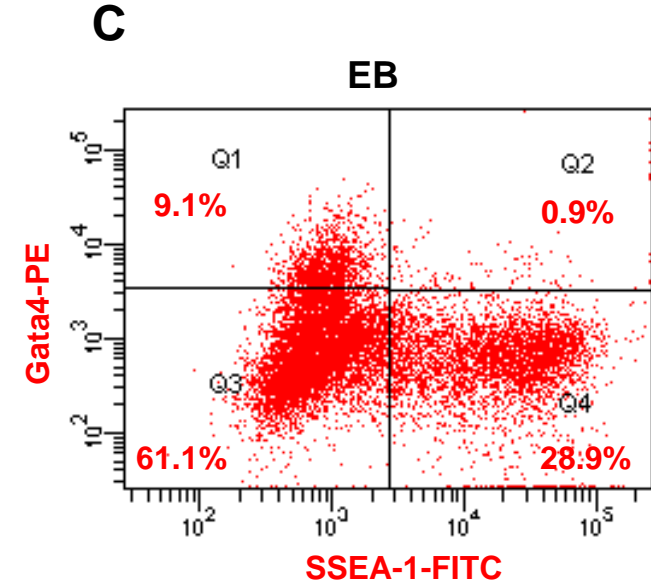
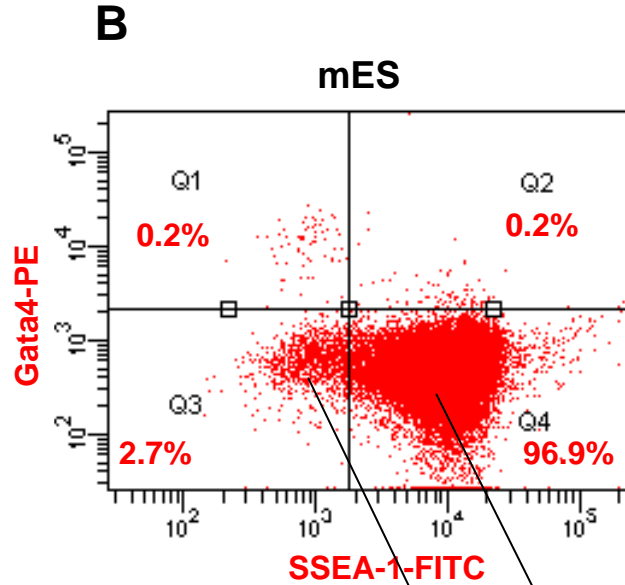
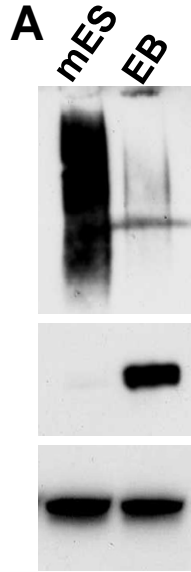


-Human ESC H9 grown on mouse embryonic fibroblast feeder layer

-Cells dissociated into a single cell suspension

Differentiated Cells

Analysis of GATA4 expression during differentiation of mouse embryonic stem cells (mESs) into embryoid bodies (EBs)



SSEA-1: mES pluripotent marker

GATA4: marker of cardiac mesoderm

Majority of cells are undifferentiated

Differentiated cells



Neuronal Progenitor Cells

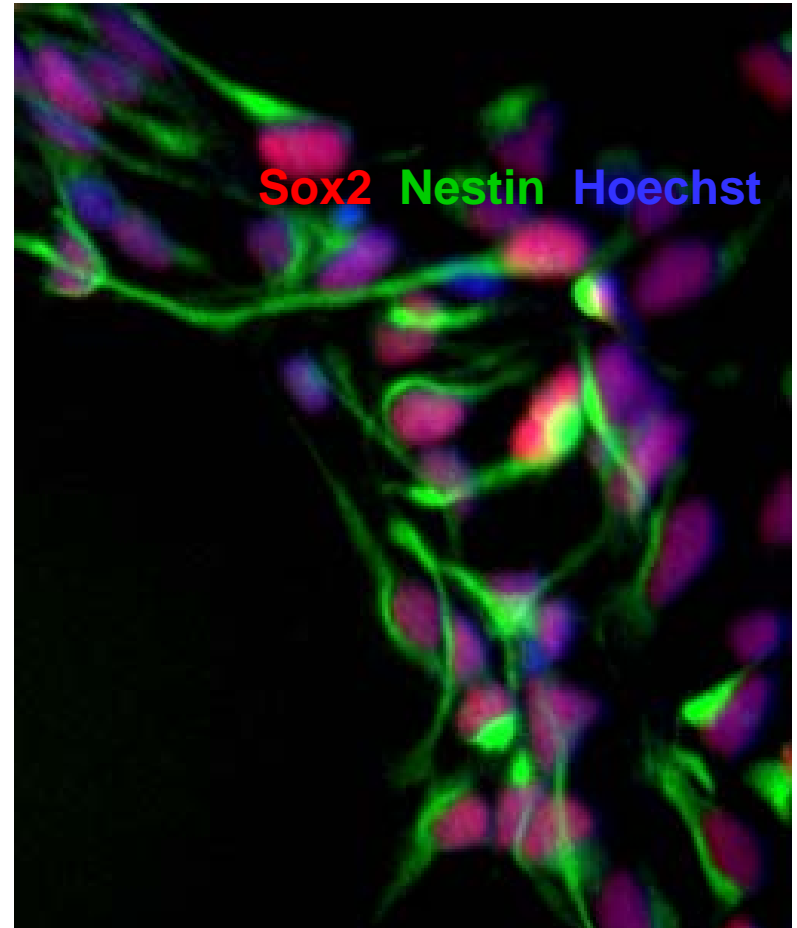
*For Research Use Only. Not for use in diagnostic or therapeutic procedures.
Class I (1) Laser Products*

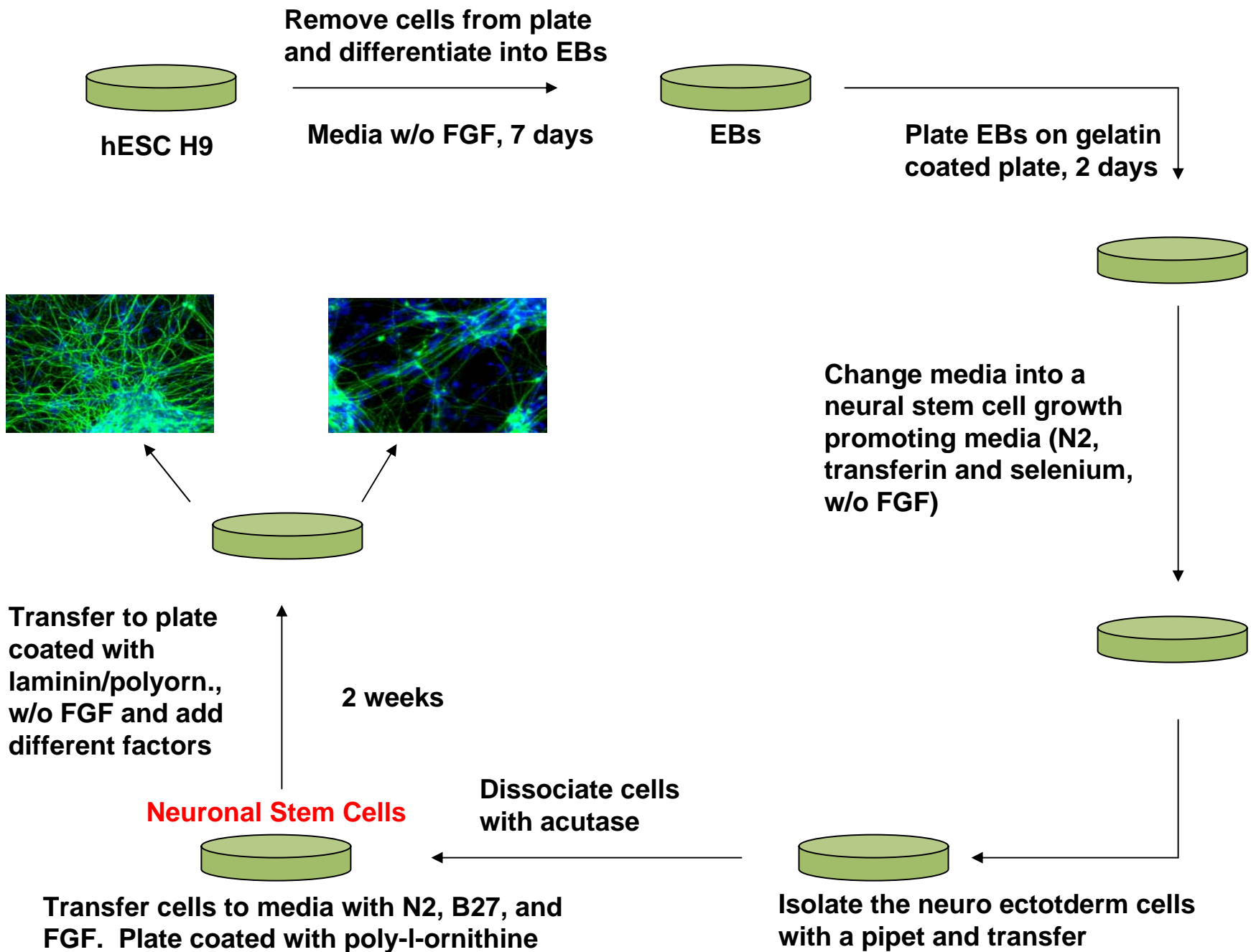


Analysis of neural differentiation of human embryonic stem cells

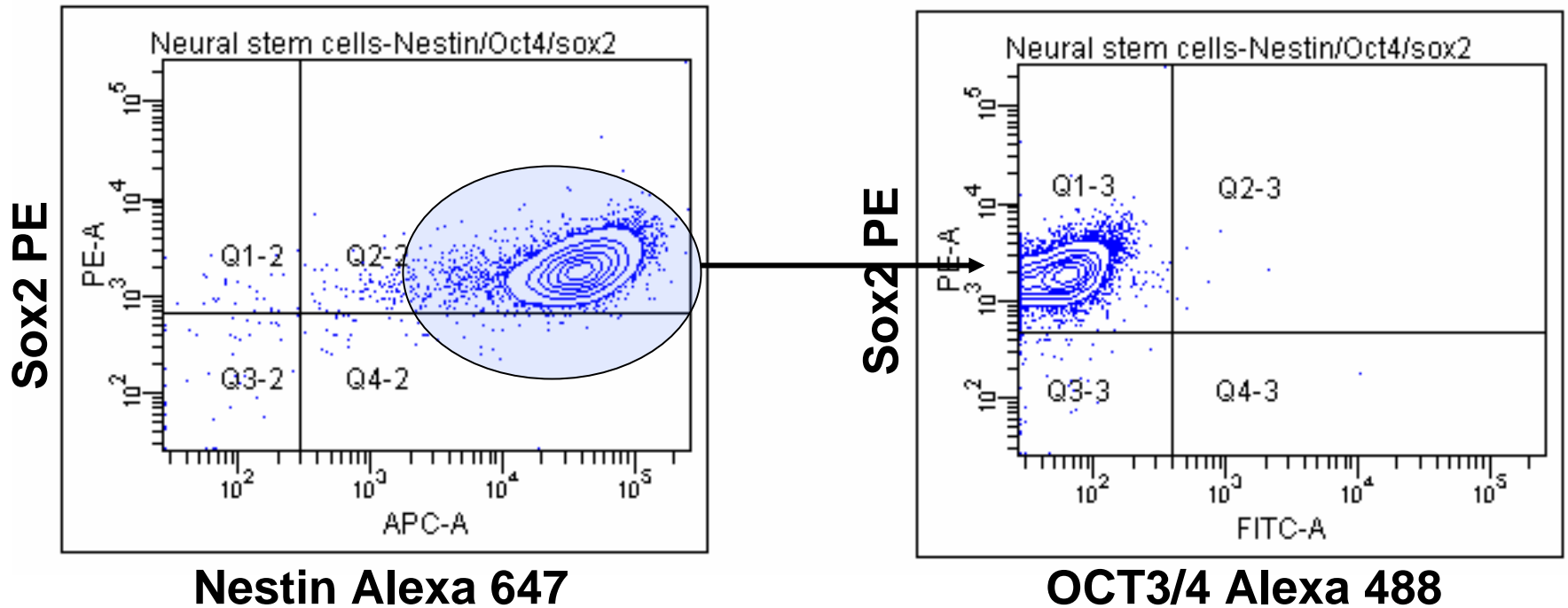
- Neural stem cells (NSCs) can be differentiated from hESCs.
- NSCs can self-renew for many passages and can be induced to differentiate into a variety of post-mitotic neural cell types.
- Assessing purity of NSCs is crucial to control for batch-to-batch variability.
- Identifying cell surface molecular signatures of NSCs and neurons will allow for rapid purification of specific neural subtypes from heterogeneous populations.

hESC (H9)-derived neural stem cells





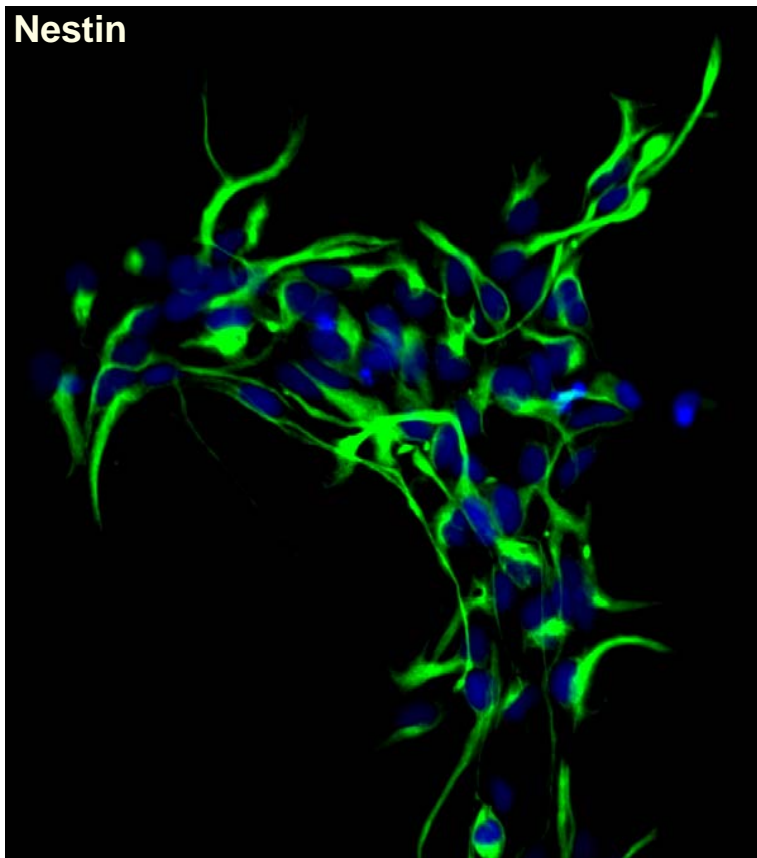
Flow cytometric quantification of hESC-derived neural stem cells



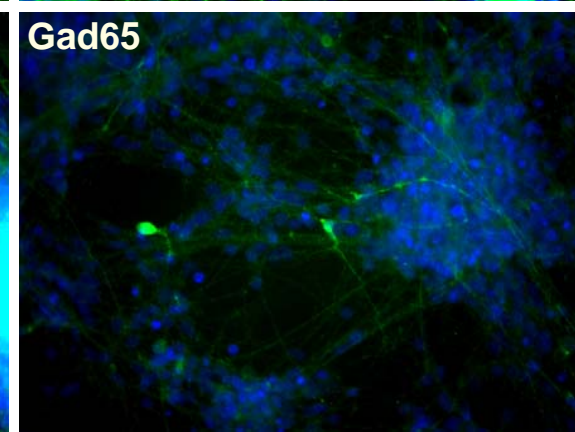
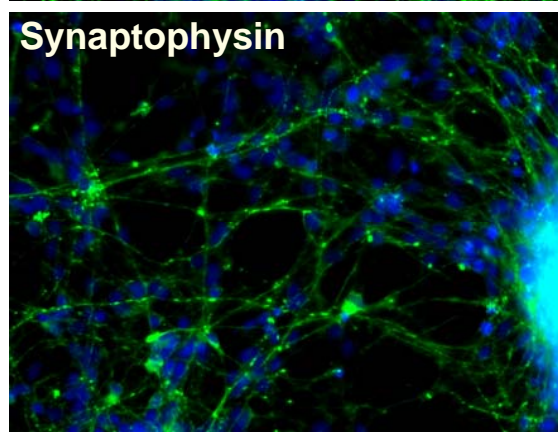
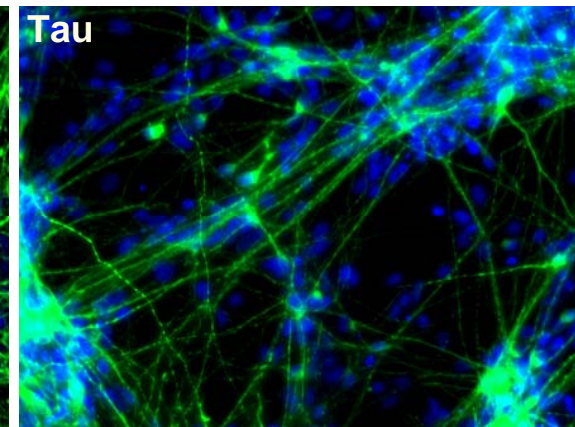
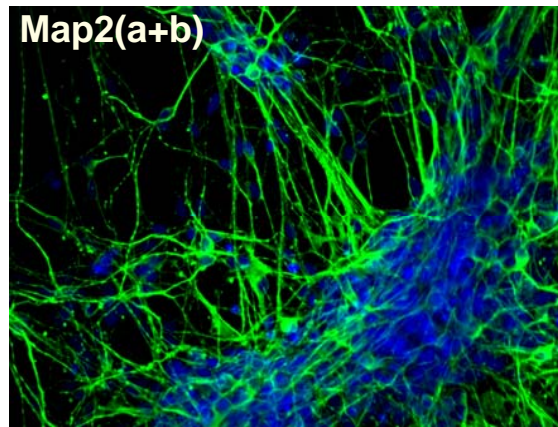
We are able to generate a near-pure population of neural stem cells expressing Sox2 and Nestin

Human ESC-derived neural stem cells can also be differentiated into neurons

A




B



A. hESCs (H9) were grown on hESC-qualified BD Matrigel™ and mTeSR1 media (Stem Cell Technologies). hESCs were differentiated into neural stem cells similarly as described by Yeo, et al. (2007)

B. Differentiation of H9-derived neural progenitors into neurons (2 weeks)



Finding New Stem Cell Markers using Flow Cytometry

*For Research Use Only. Not for use in diagnostic or therapeutic procedures.
Class I (1) Laser Products*



CD marker analysis of hESC (H9) and neural stem cells by flow cytometry

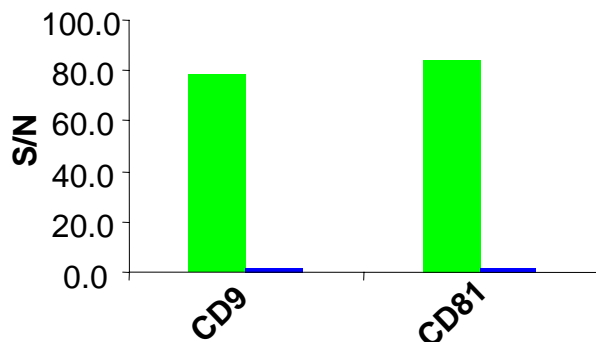
- Pilot screen of 21 CD markers was performed (partial analysis of >250 cell surface markers)
- Goal is to define cell surface molecular signatures of both cell types and potentially identify subpopulations
- Next step is to include differentiated neurons



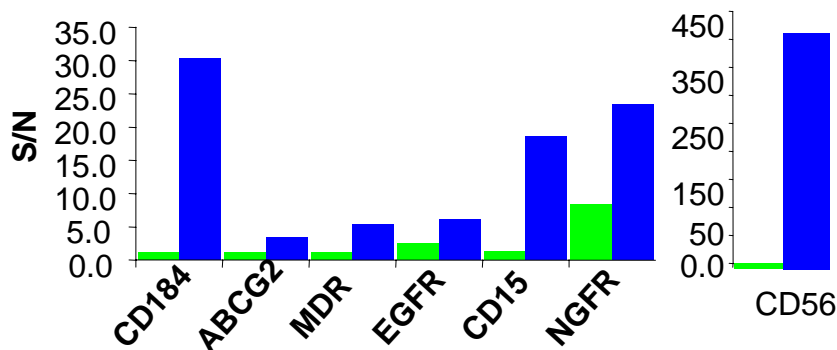
CD marker analysis of hESC (H9) and neural stem cells by flow cytometry

Signal to Noise over Isotype control

High in hESC, negative in NSC



High in NSC, negative to low in hESC

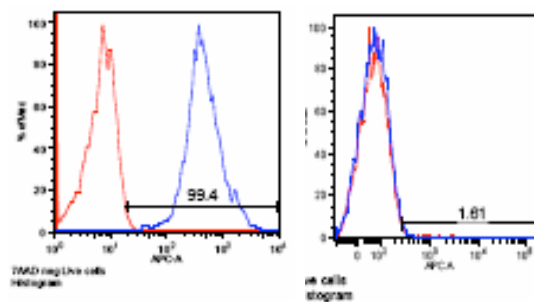


■ Hu ESC (H9)
■ NSC

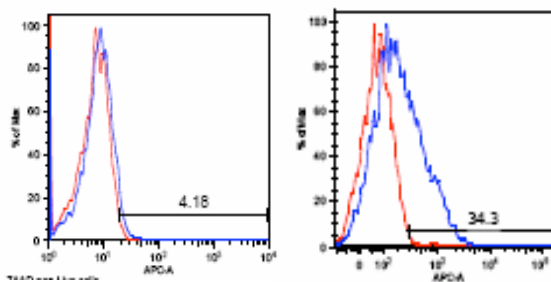
Representative data

H9

NSC



CD81



ABCG2

Significance

Potential markers of pluripotency

Potential markers of neuroectoderm, NSC or of adult SC but not of pluripotent stem cells

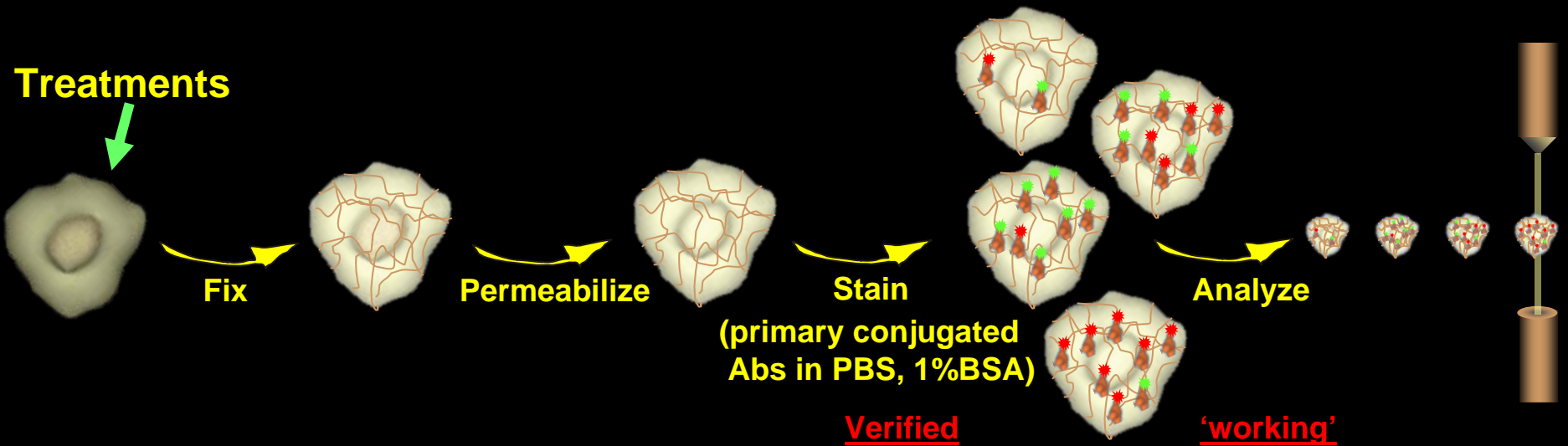


Phosphorylation Studies

*For Research Use Only. Not for use in diagnostic or therapeutic procedures.
Class I (1) Laser Products*



Cell Surface and Intracellular Staining for Flow Cytometry



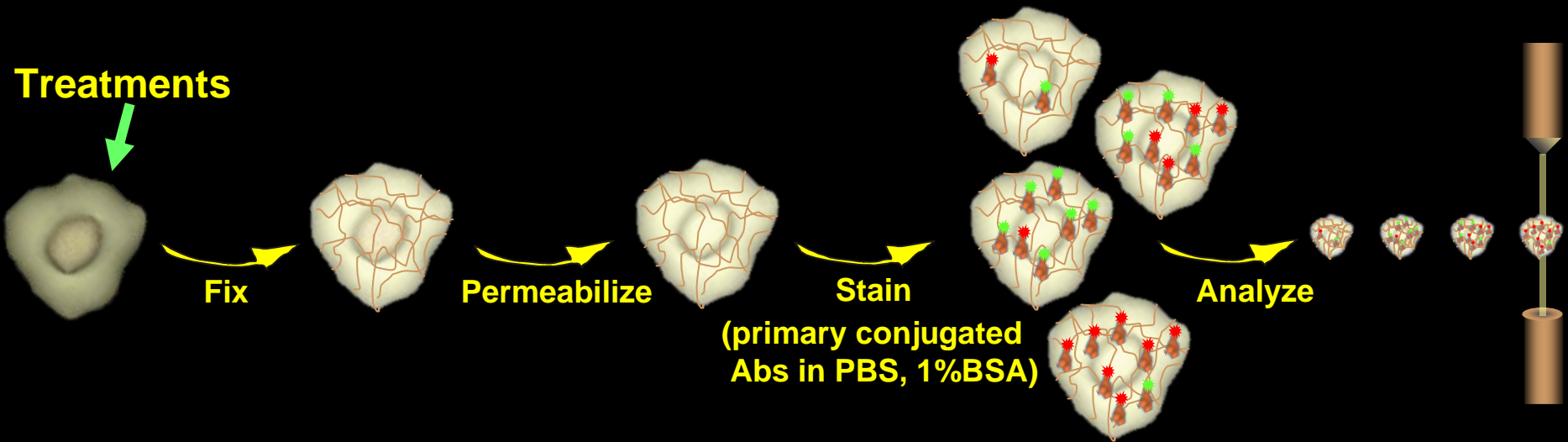
Thorough development of fixation protocols for cell lines and whole blood (immediately out of fresh samples).

p38 MAPK
 JNK, cJun
 AKT, PIP2, PIP3,
 PKC $\alpha/\beta/\theta/\delta$, Rsk
 Raf, Mek, ERK, ELK
 Rsk, Creb,
 STATs, SRC
 CREB, cJUN, IKK α
 p53 s15, s20 s37, s392
 Pyk2, Shc, Fak, Src
 Slp76, Zap70, Syk, Lat, Vav,
 Lck, PLC γ
 Beta-integrins

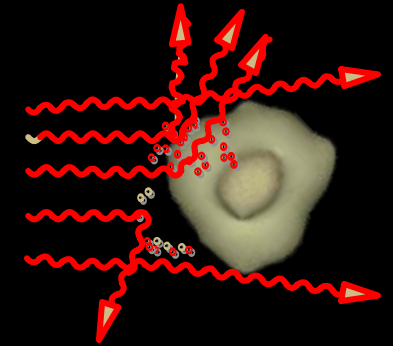
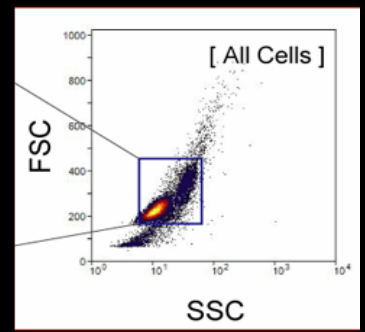
EGFR
 PDGFR
 cKit
 VEGFR
 PKA
 RB
 NFAT
 NF- κ B p65
 Caveolin
 Paxillin
 FLT3
 MEKS

>80 specificities

Cell Surface and Intracellular Staining for Flow Cytometry

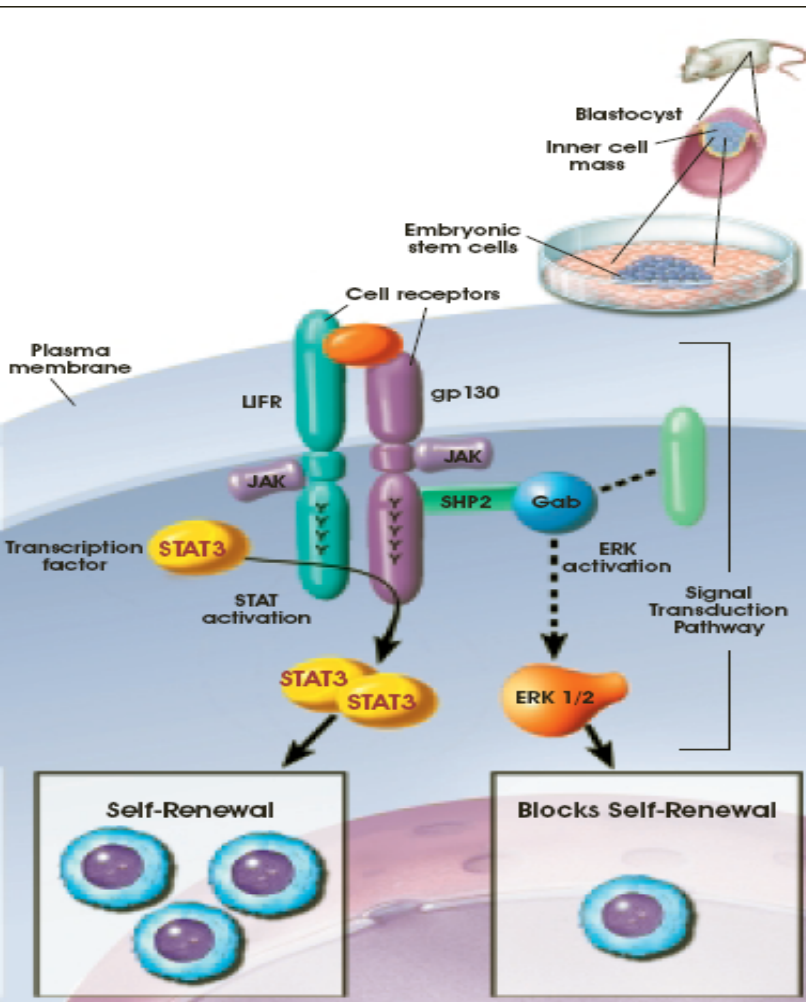


1. State-specific antibodies: phospho-specific antibodies and others
2. Adopting entirely new fluorophores
3. Generation of efficient conjugation, purification, and testing protocols



- Analysis Steps:**
1. Identify live cells
 2. Identify cell type
 3. Evaluate cell signal

Activation of STAT3 is Essential for Self-Renewal of Mouse Embryonic Stem Cells

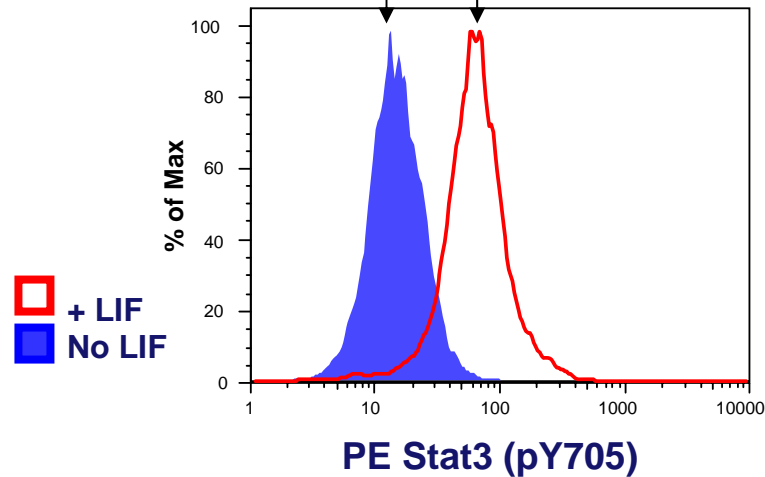
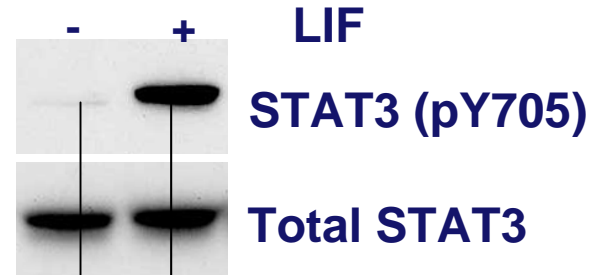
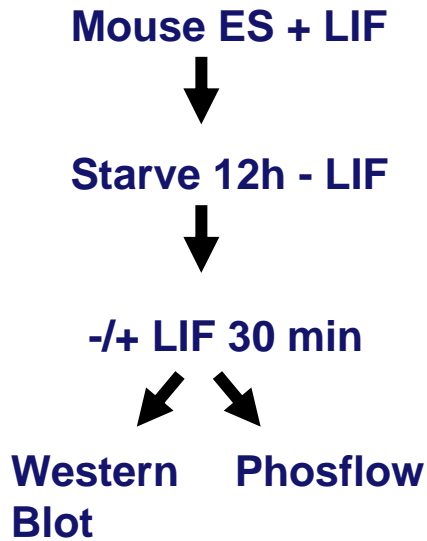


© 2001 Terese Winslow, Lydia Kibiuk

- Niwa H. et al. *Gene & Dev.* 12, 2048, 1998
- Raz R. et al. *PNAS*, 96, 2846, 1999
- Matsuda T. et al. *EMBO J.* 18, 4261, 1999

- Mouse ES cells can be maintained in a proliferative undifferentiated state in vitro with the addition of LIF into the culture medium
- LIF binds to a two-part receptor complex composed of the LIF receptor and the gp130 receptor
- The binding of LIF triggers the activation of STAT3

LIF-Dependent STAT3 Signaling Can Be Measured Using BD™ Phosflow Technology





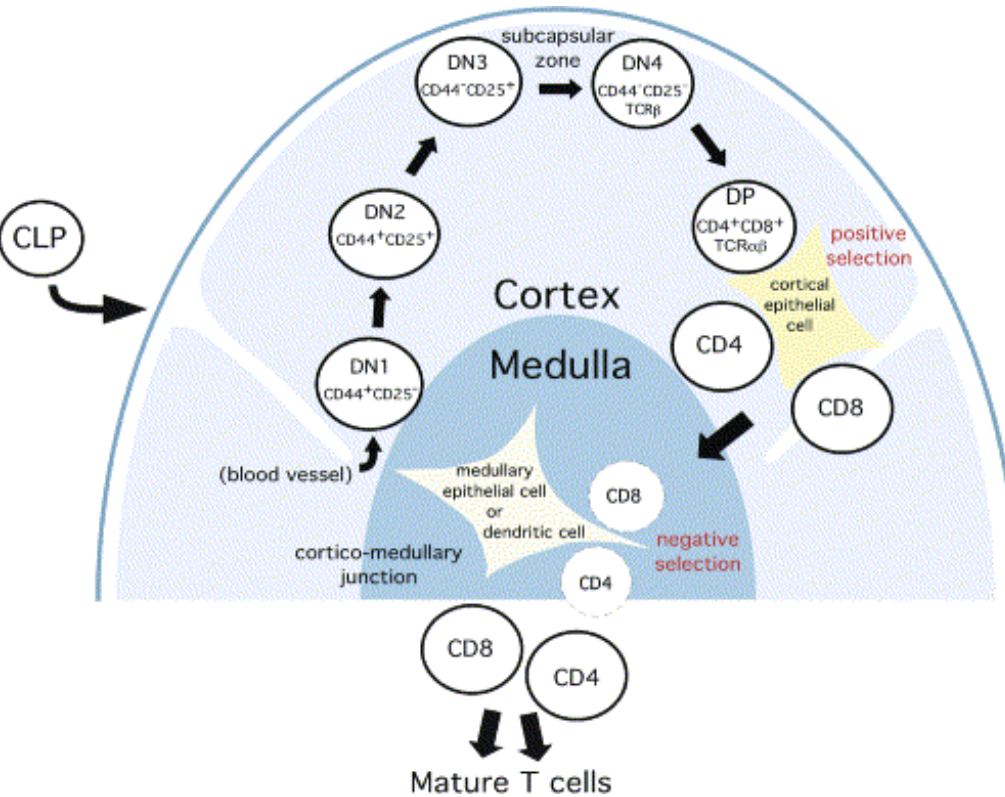
Intrathymic T-cell Development

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

Class I (1) Laser Products



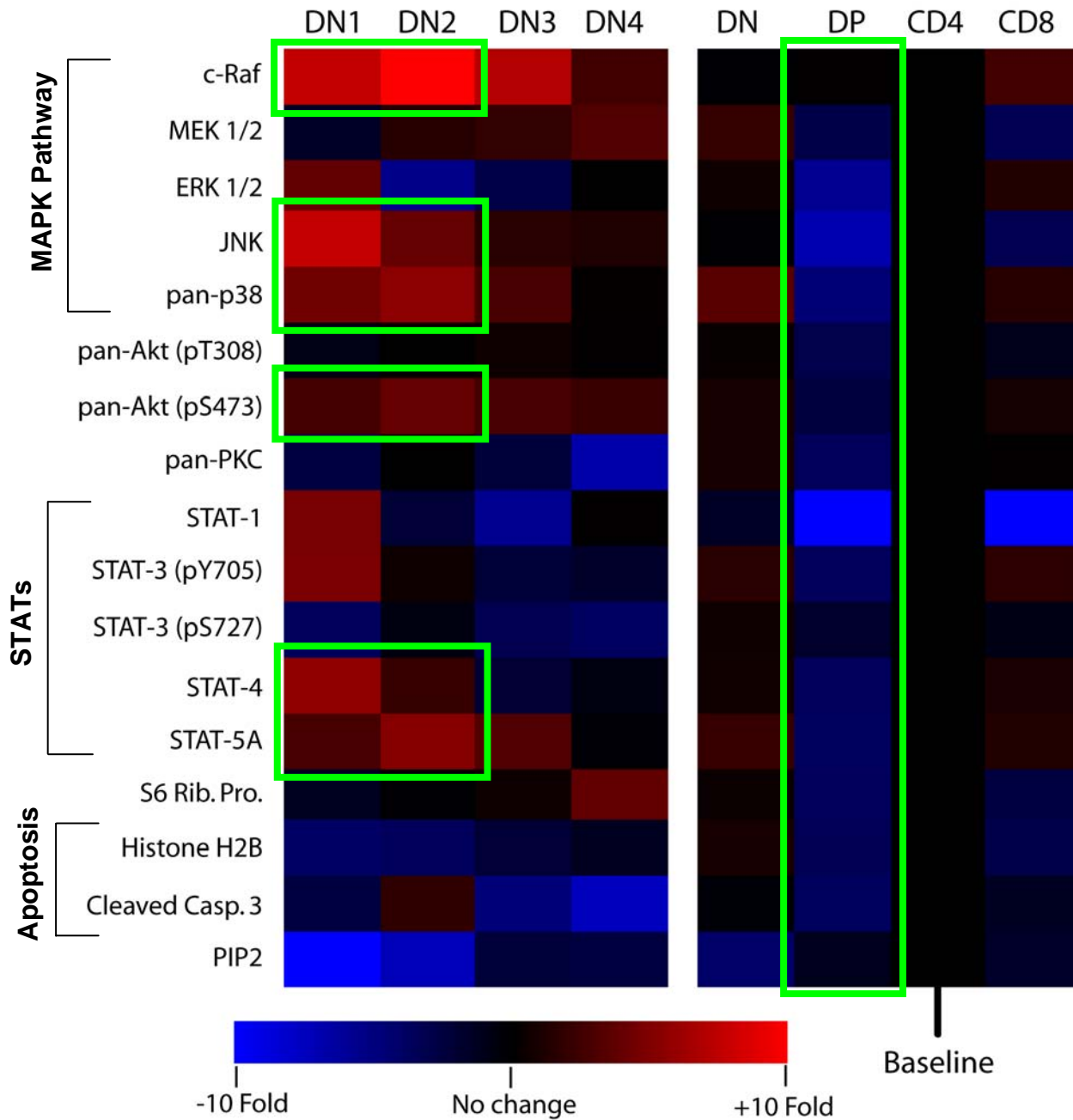
Intrathymic T-Cell Development



- **Commitment to T lineage: Earliest stage is ETP** ($Sca-1^+ Lin^- Thy1.1-c-kit^{hi} CD62L^+$)
- **Migration of maturing T cells in the thymic cortex: DN1-4, DP, positive selection of DP cells**
- **Medulla: Negative selection of SP cells**
- **Migration of mature SP cells from the thymus**

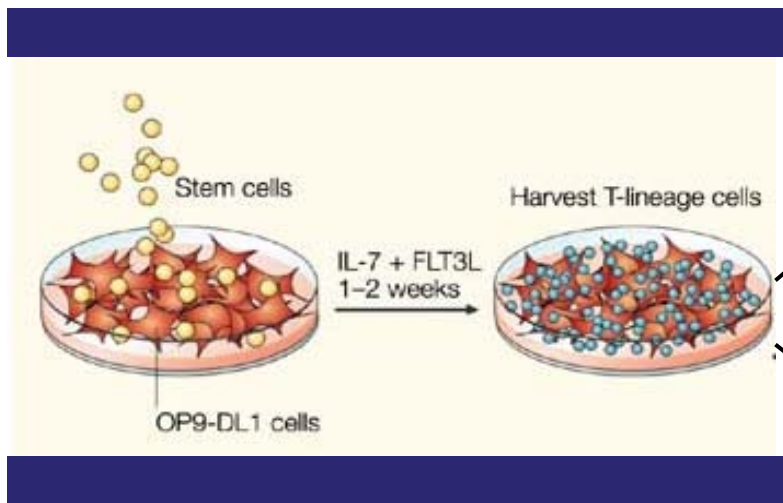


Normal B6 Thymus



T-Cell Development on OP9-DL1 Cells

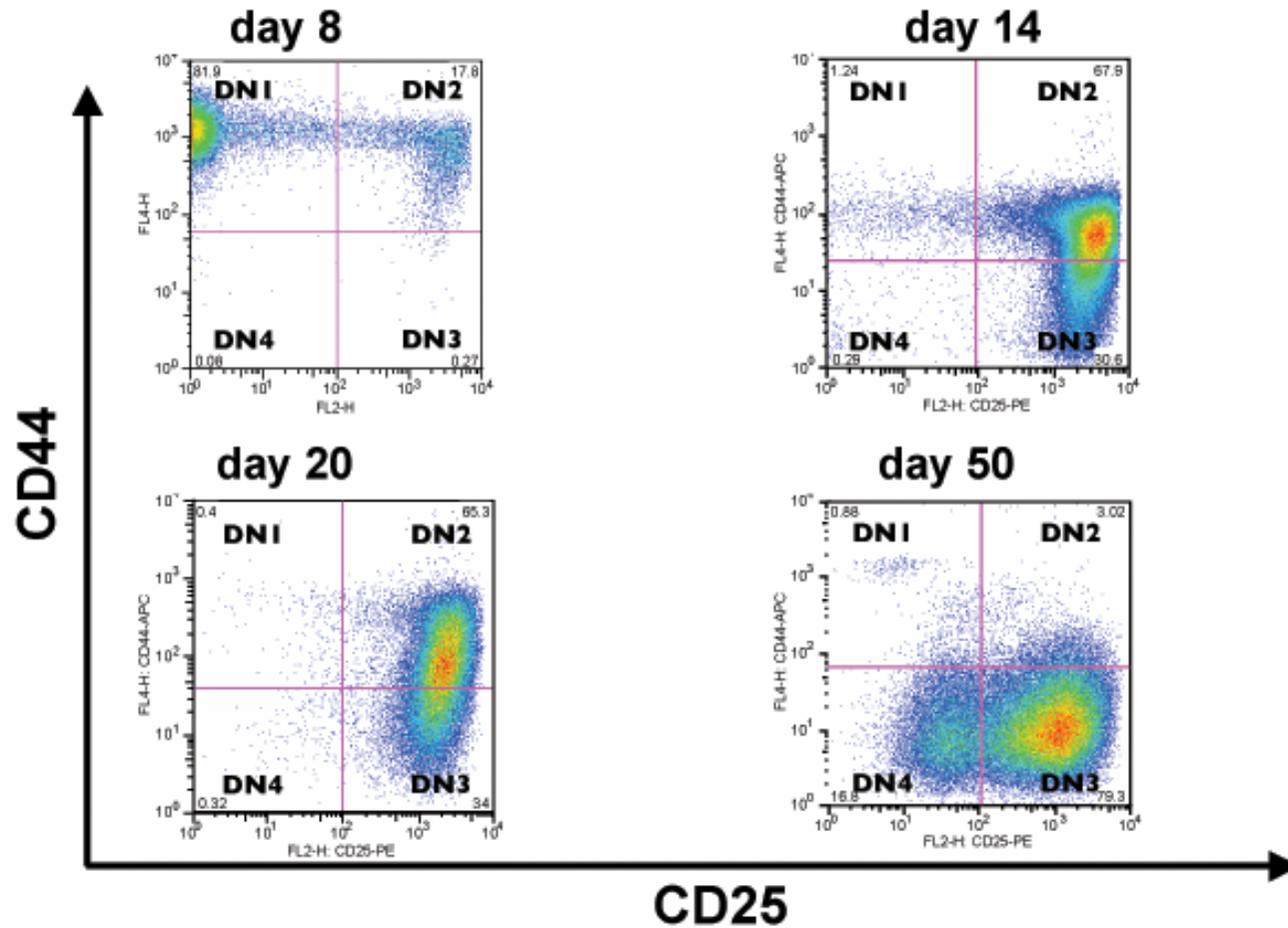
- OP9-DL1: Murine BM stromal cell line that has been retrovirally transduced to express the Notch 1 receptor ligand Delta-like 1 (DL1)
- Culture of HSC on OP9-DL1 cells in the presence of IL-7 and FLT3L induces T-cell differentiation



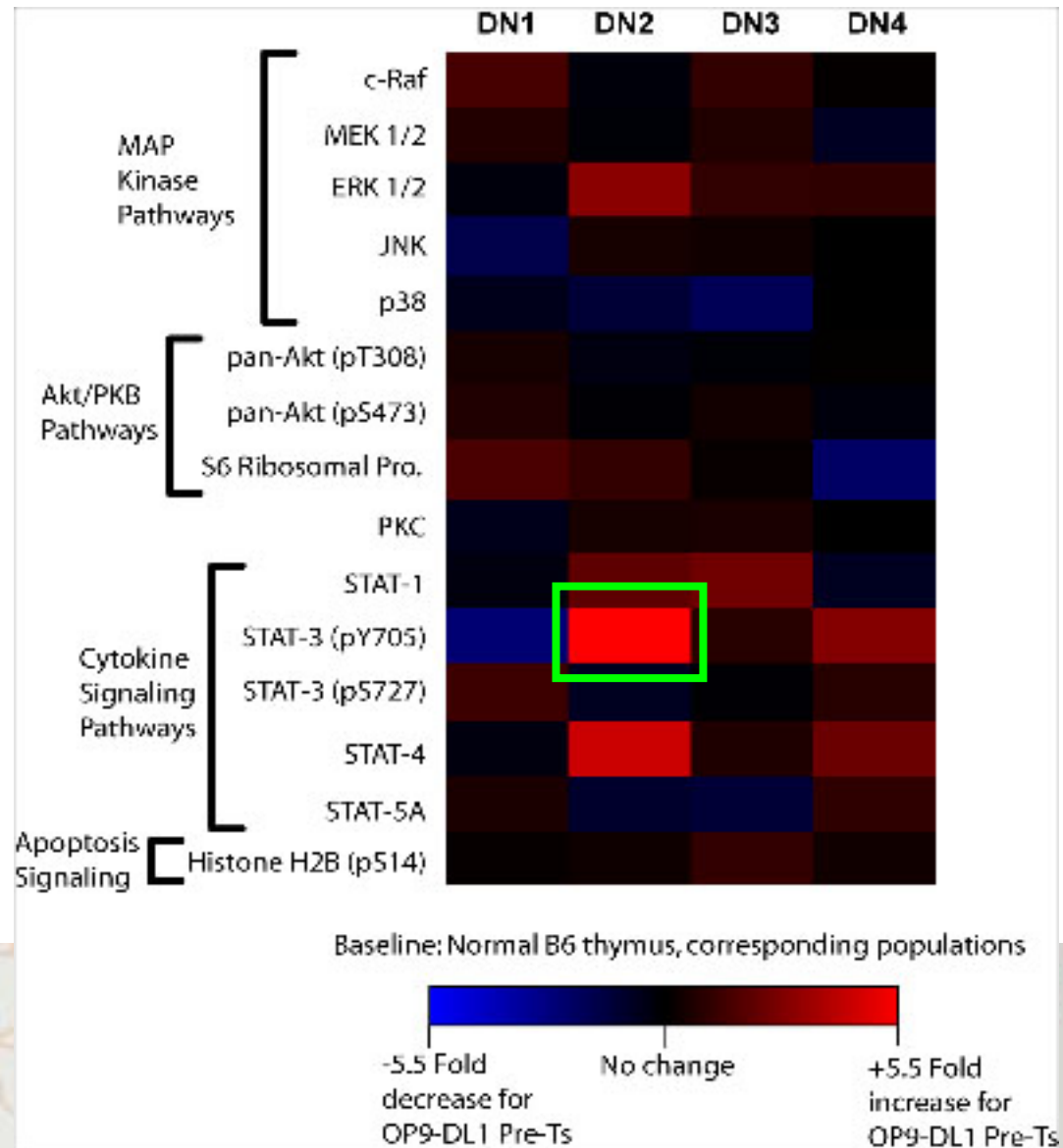
In vitro analysis

Adoptive transfer into allogeneic HSCT recipients

T-Cell Development from Adult Murine BM Derived HSC In Vitro Using OP9-DL1 Cells Generates Primarily DN2-DN3 Precursors



OP9-DL1 Derived T-Cell Precursors



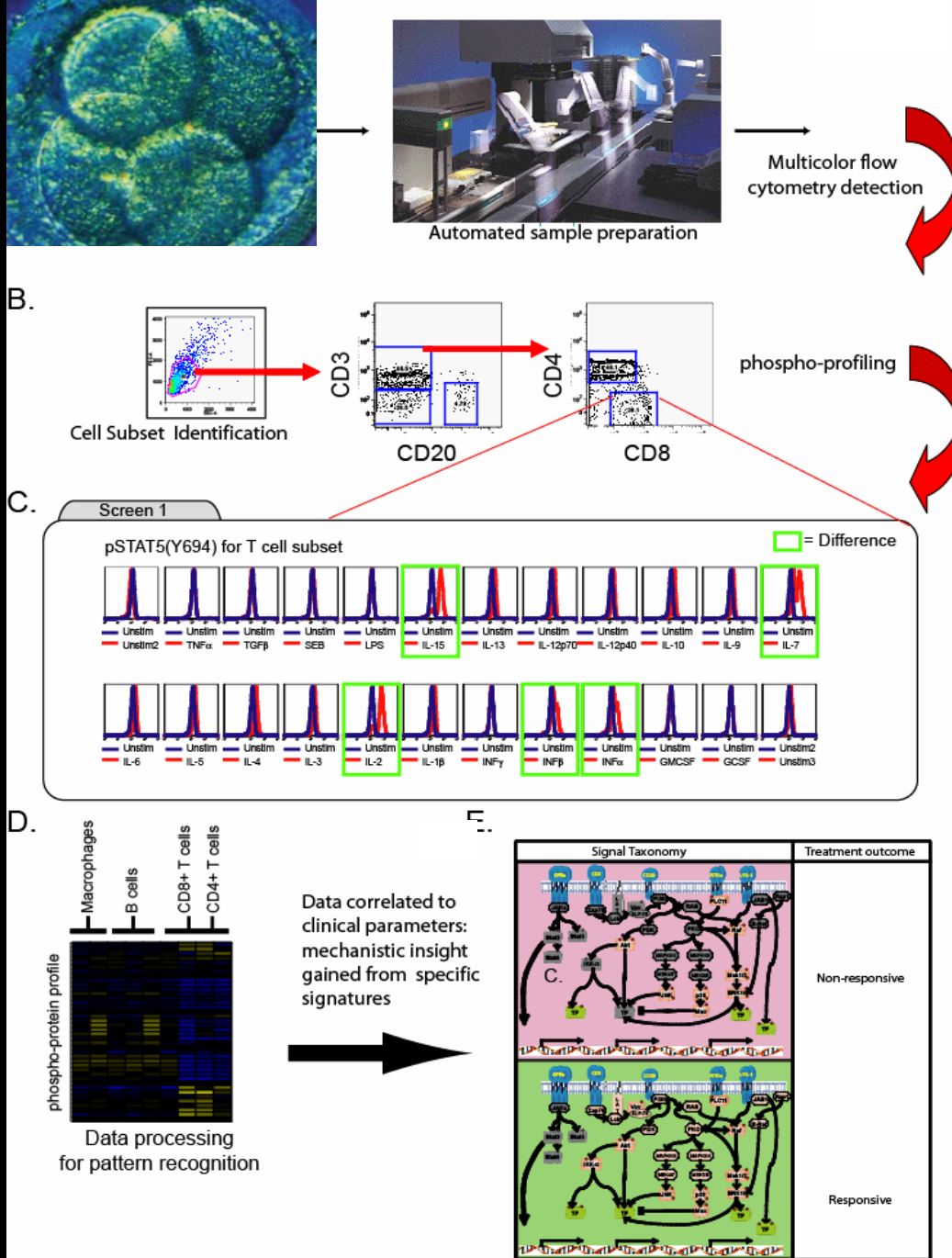
Summary of Findings for the Thymus

- DP thymocytes: Low signaling profile
- DN1/DN2 thymocytes: Increased Raf, Jnk, p38, Akt, and STAT-4 and -5 signaling
- DN and SP thymocytes phosphorylate STAT-5A in response to challenge with IL-7

1

2

3



Acknowledgments

BD Biosciences

- Christian Carson
- Frederick Princen
- Guo-Jian Gao
- Li Li
- Chad Sisouvanthong
- Jerome Zawadski
- Anissa Agadir

Memorial Sloan Kettering Cancer Center

- Sydney Lu
- Onder Alpdogan
- Marcel van den Brink

Stanford University

- Peter Krutzik
- Garry Nolan





Technical Support (US)

e-mail: techserv@bd.com

Phone: 877-232-8995

prompts #3, then #2

Please visit our Stem Cell Source web page:

www.bdbiosciences.com/stemcellsource/

