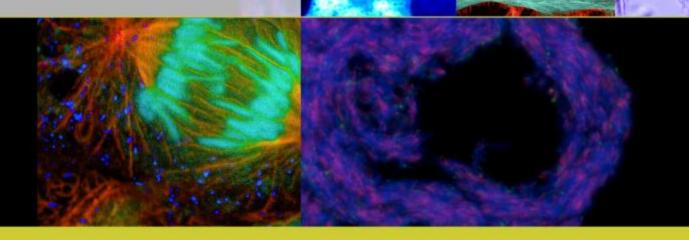


James Clinton, Ph.D. Scientist, ATCC February 19, 2015





### **About ATCC**

- Founded in 1925, ATCC is a non-profit organization with headquarters in Manassas, VA
- World's premiere biological materials resource and standards development organization
- ATCC collaborates with and supports the scientific community with industry-standard products and innovative solutions
- Broad range of biomaterials
  - Continuous cell lines, iPSCs, primary cells, and hTERT immortalized cells
  - Bacteria, fungi, yeasts, protists, and viruses
  - Microbial and tumor cell panels
  - Genomic and synthetic nucleic acids
  - Media, sera, and reagents





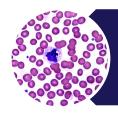




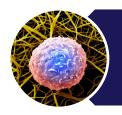
### **Outline**



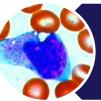
## Background



CD34+ hematopoietic stem & progenitor cells



Mononuclear cells

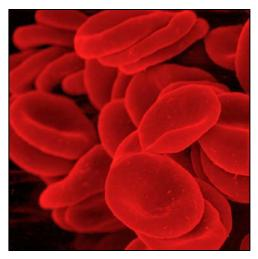


CD14+ monocytes



### **Blood cells**

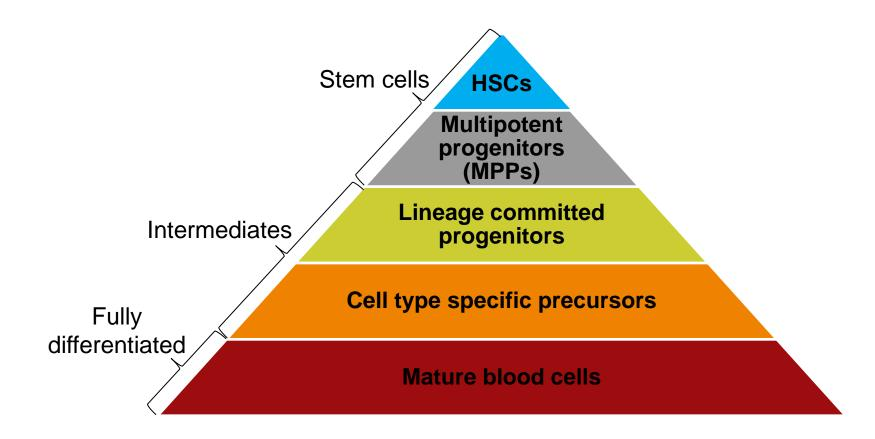
- Blood is comprised of a heterogeneous population of specialized cells
- Three major types
  - Leukocytes: Acquired and innate immunity
  - Erythrocytes: Gas transport
  - Thrombocytes: Wound healing
- Millions of blood cells are generated every second, approximately 1 trillion every day
- Hematopoiesis:
  - Dynamic process
  - Varies in response to injury or infection
  - Individual cells may live for hours to years
- Blood cells arise from hematopoietic stem cells (HSCs)



Erythrocytes

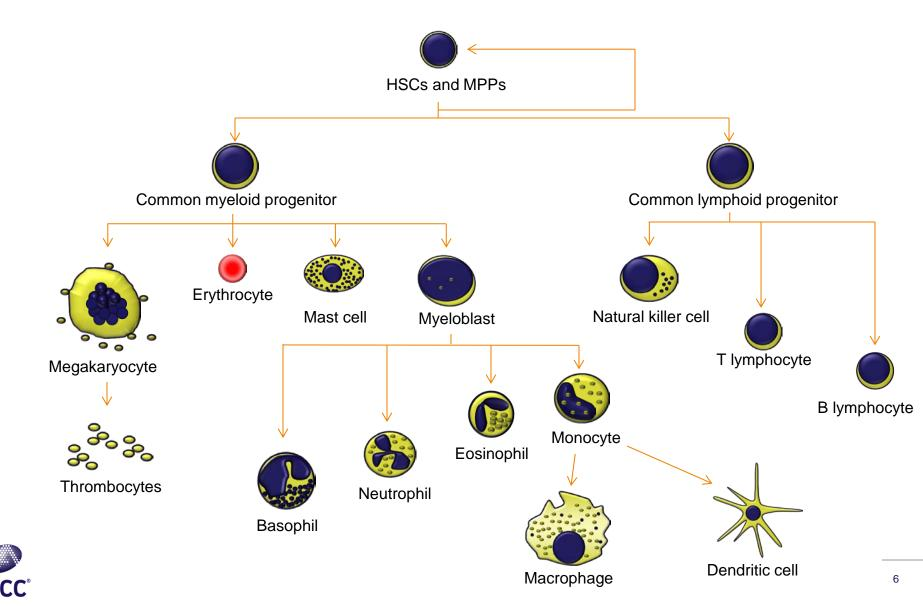


# Hematopoiesis: A hierarchal system





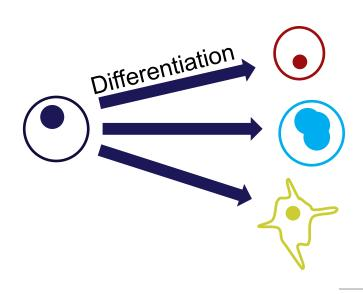
# Hematopoietic cell fate and lineage



# Hematopoietic stem cells: Characteristics

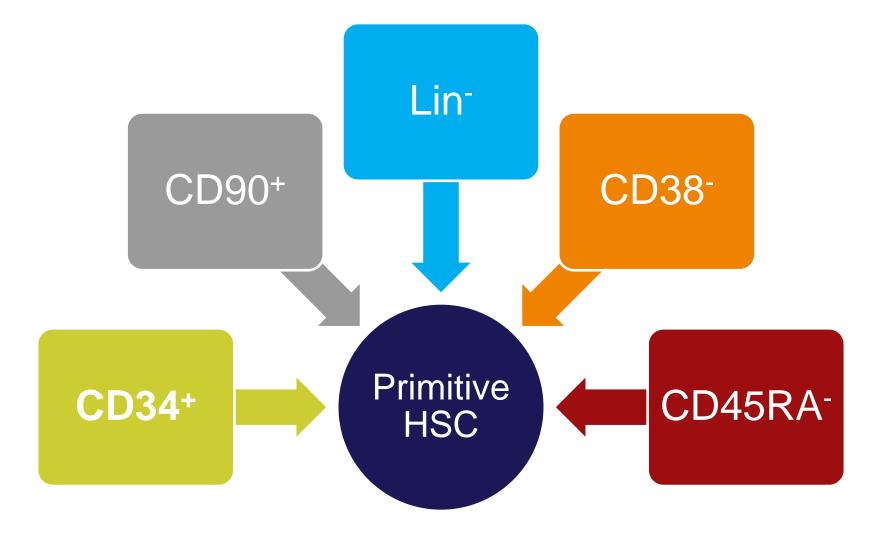
- Hematopoietic stem cells (HSCs) are multipotent cells that give rise to all other blood cells
- HSCs reside primary in bone marrow (major site of hematopoiesis in adults)
- True hematopoietic stem cells are rare
- True hematopoietic stem cells can only be confirmed via in vivo functional assays





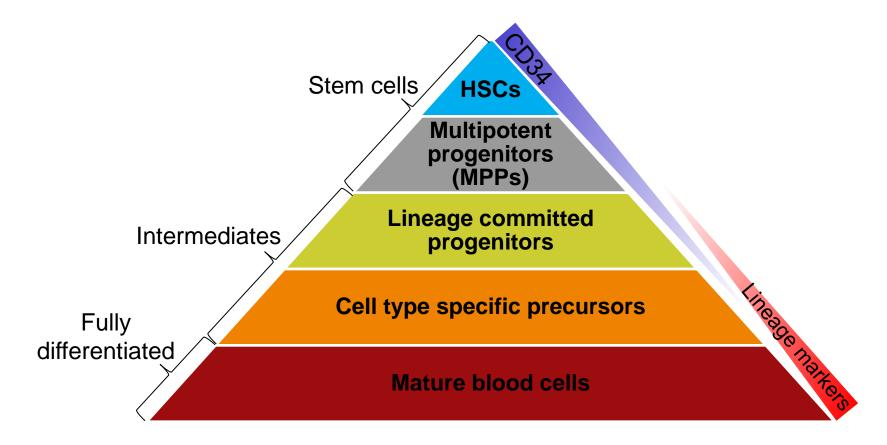


# Hematopoietic stem cells: Markers



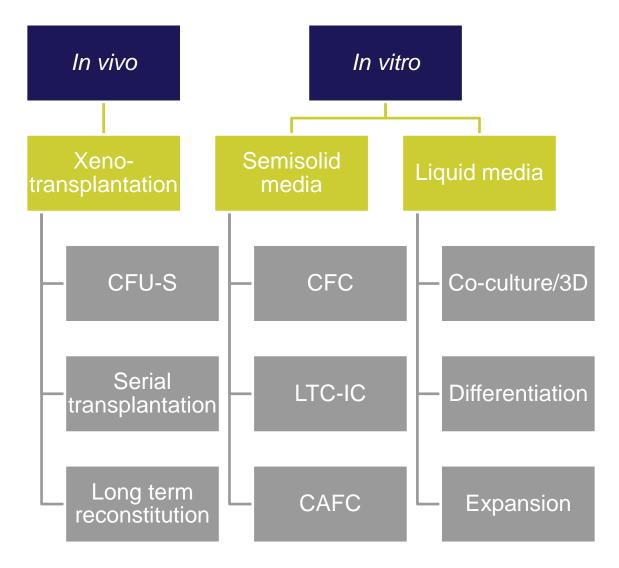


# CD34+ cells are a mixed population of stem and progenitor cells



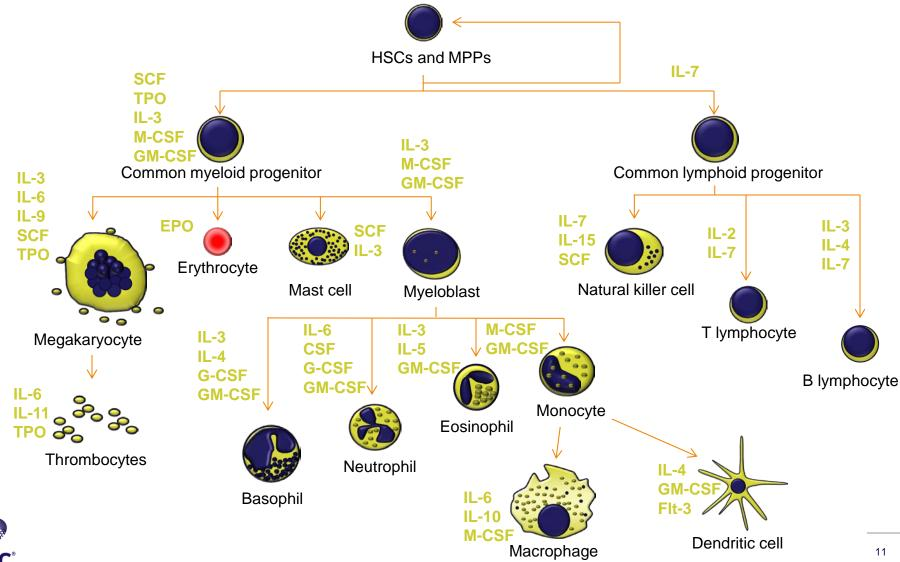


# Assays to study hematopoietic stem and progenitor cells





# Cytokines influence cell fate and lineage in-vitro





# **Blood and hematopoiesis summary**

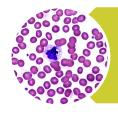
- Blood is a heterogeneous tissue its replacement in vivo is a complex process
- HSCs are responsible for the generation of all other blood cell types
- Recent advances allow for the identification and isolation of human HSCs as well as other blood cell types
- This process can be studied in vitro though the use of lineage directed differentiation of HSPCs



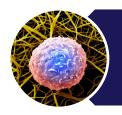
### **Outline**



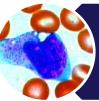
## Background



CD34+ hematopoietic stem & progenitor cells



Mononuclear cells



CD14+ monocytes



# **Primary CD34+ HSPCs**

#### **Applications**

- Ex vivo expansion and differentiation
- Stem cell markers
- Gene transfer
- Cytokine and chemokine expression and regulation
- Receptor expression

#### **Key research areas**

- Safely and efficiently expand HSCs in vitro or in vivo for transplantation
- Immune response
  - Graft-versus-host disease/transplant rejection
- Cancer
- Cell-fate determination



# Mature lymphocyte generation from CD34+ HSPCs on a 3D matrix

### A Simple Model System Enabling Human CD34\* Cells to Undertake Differentiation Towards T Cells

Antonio Lapenna<sup>1,2</sup>, Christopher B-Lynch<sup>1</sup>, Chrysa Kapeni<sup>1</sup>, Richard Aspinall<sup>1</sup>

1 Regenerative Medicine Group, Crunfield Health, Cranfield University, Cranfield, United Kingdom, 2 Department of Immunology and Cancer Research, Faculty of Medicine, the Hebrew University of Janussiem, Janussiem, Israel

#### Abstrac

Background: Channelling the development of haematopoietic progenitor cells into T lymphocytes is dependent upon a series of extrinsic prompts whose temporal and spatial sequence is critical for a productive outcome. Simple models of human progenitor cells development depend in the main on the use of xenogeneic systems which may provide some limitations to development.

Methods and Findings: Here we provide evidence that a simple model system which utilises both humen heratincoyte and fibroblast cell lines arrayed on a synthetic tartalum coated matrix provides a permissive environment for the development of human CD34\* haematopoietic cells into mature CD4\* or CD8\* T lymphocytes in the presence of Interleukin 7 (IL-7), Interleukin 15 (IL-15) and the Fms-like tyrosine kinase 3 sigand (Fil-SL). Interleukin 5 system was used to compare the ability of CD34\* cells to produce mature thymocytes and showed that whilst these cells derived from cord blood were able to producitively differentiate into thymocytes the system was not permissive for the development of CD34\* cells from adult peripheral blood.

Conclusions/Significance: Our study provides direct evidence for the capacity of human cord blood CD34\* cells to differentiate along the T lineage in a simple human model system. Productive commitment of the CD34\* cells to generate T cells was found to be dependent on a three-dimensional matrix which induced the up-regulation of the Notch delta-like ligand 4 (DII-4) by epithelial cells.

Citation: Lapenna A, B-Lynch C, Kapeni C, Aspinall R (2013) A Simple Model System Enabling Human CD34\* Celle to Undertake Differentiation Towards T Cells. PLoS ONE 5(7): e69572. doi:10.1371/journal.pone.0069572

Editor: Zoran Ivenovic, French Blood Institute, France

Received April 05, 2013; Accepted June 14, 2013; Published July 23, 2013

Copyright: © 2013 Lapenna et al. This is an open-access article distributed under the terms of the Creative Commons Attribution License, which permits unrestricted use, distribution, and reproduction in any medium, provided the original author and source are credited.

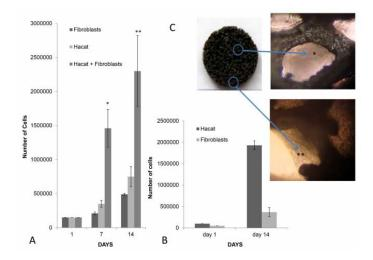
\* E-mail: antoniolapenna@hotmail.com

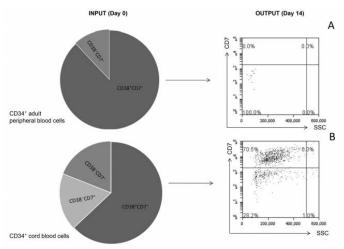
#### Introduction

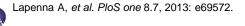
The generation of T cells from haematopoietic progenitor cells requires the positioning of progenitors within the thymus where a unique environment induces supports and directs their differentiation [1]. Production of new thymocytes continues throughout the and because the progenitions cannot be stored and maintained indefinitely within the thymus, continuation of production requires seeding of the thymus with these cells. Analysis of thymic output reveal that the rate of production of new T cells declines with age [2] and that as thymocyte production decreases so there is attorpty of the thymus.

In broad terms thymic atrophy has been linked to deficits in

(FTOC) systems or allogenetic cell lines such as mouse bore marrow-derived OP9 cells expressing the Notch delta-like ligand 1 (OP9-Dil1) [3-9]. But the experiments in human systems have proved more intractable. Analysis of the capacity of haematopoietic progenitor cell populations to produce Tels have proceeded but has been hampered, mainly through the use of xenogenetic model systems which by their very nature are limited and associated with incomplete or inefficient differentiation of the progenitors [5]. Some studies of thymic stromal cells have indicated changes with age in the thymic environment cell type composition and expression profile but these data were limited by the lack of culture methods which could effectively model the thymic architecture in vitro [6].



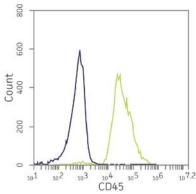


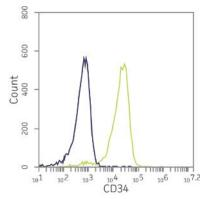


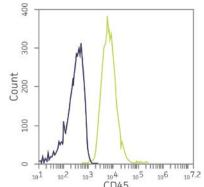
# **ATCC primary CD34+ HSPCs**

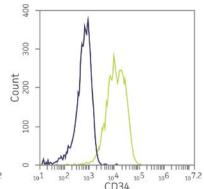
- Healthy human volunteer donors; IRB-approved informed consent
- Adult, non-pregnant (excluding cord blood)
- Cryopreserved at P0; Purity: ≥ 90% CD34+
- Age, gender, ethnicity, and blood type on CoA

ATCC® No.	Tissue	Туре	Size
PCS-800-012	Bone marrow	Hematopoietic stem/progenitor cells (CD34+); Normal	≥ 0.5 x 10 <sup>6</sup>
PCS-800-014	Cord blood	Hematopoietic stem/progenitor cells (CD34+); Normal	≥ 0.5 x 10 <sup>6</sup>











Bone marrow

Cord blood

# CD34+ HSPC lineage directed expansion and differentiation

#### Goals

- Demonstrate primary CD34+ HSPC capacity for lineage-directed expansion and differentiation in vitro
- Confirm multilineage differentiation (erythrocyte, megakaryocyte, and pan-myeloid)
- Compare CD34+ differentiation efficiency from multiple tissues
- Utilize a method amenable to high throughput assays

#### **Methods**

- Cryopreserved CD34+ primary cells from bone marrow and cord blood
- Serum-free liquid culture
- Analysis of phenotype by surface marker expression
- Commercially available cytokine cocktails



# General differentiation protocol workflow

Thaw & seed

**Maintenance** 

**Analyze** 

Thaw vial of CD34+ cells

Add fresh media every ~2 days

Collect cells for analysis

Seed directly into differentiation media

Re-suspend in fresh media every 4-7 days

Basal media: SFEM II

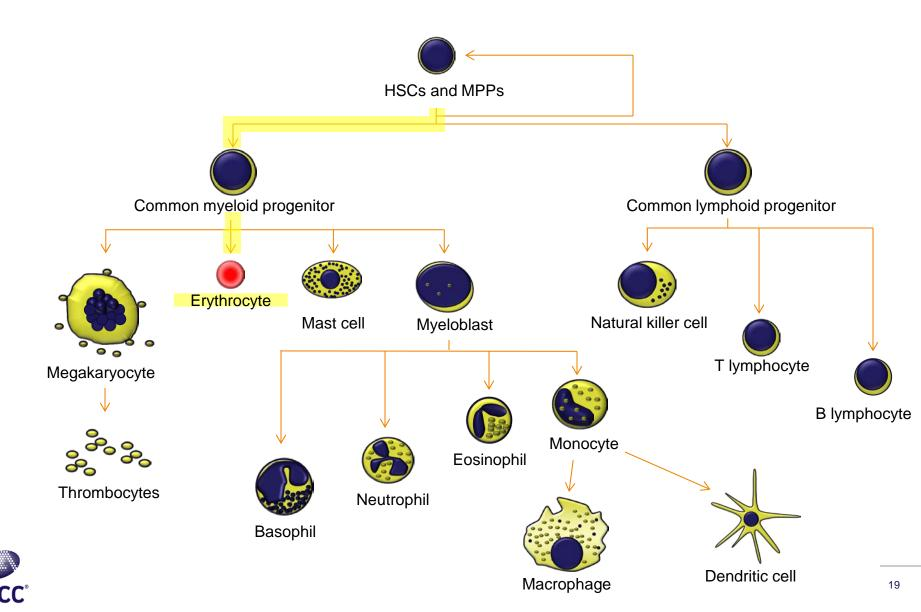
Maintain concentration below 1x10<sup>6</sup>/mL

Supplement with cytokines

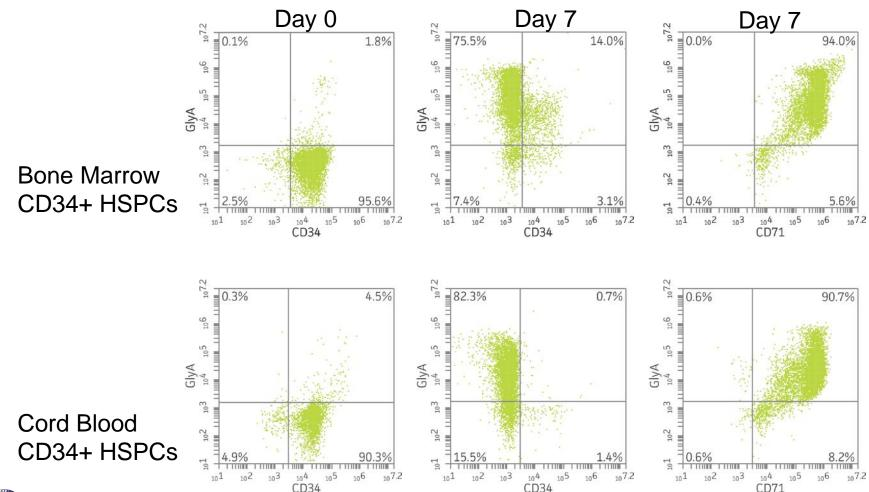
For detailed differentiation protocols see the ATCC website



# Erythroid differentiation and expansion

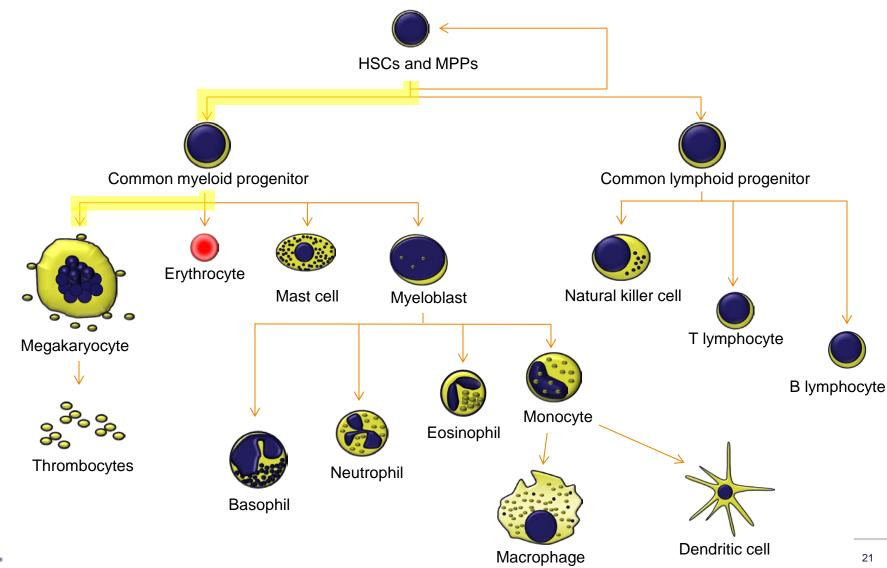


# Expression of erythroid lineage markers on differentiated BM and CB CD34+ cells



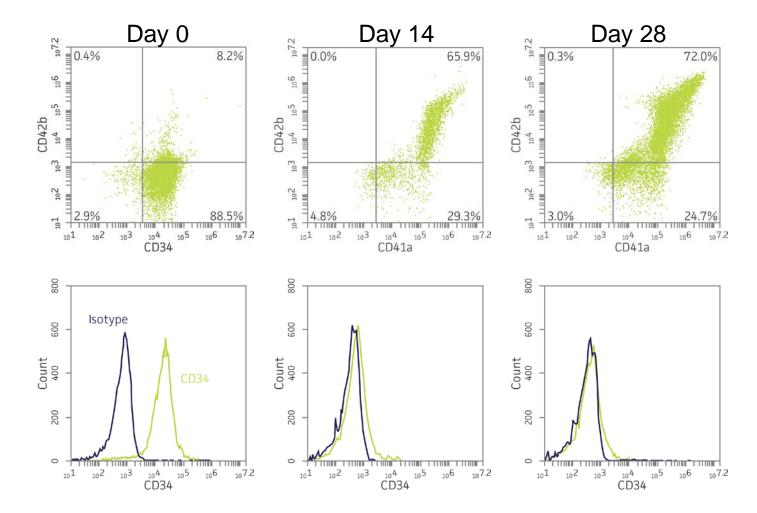


# Megakaryocyte differentiation and expansion



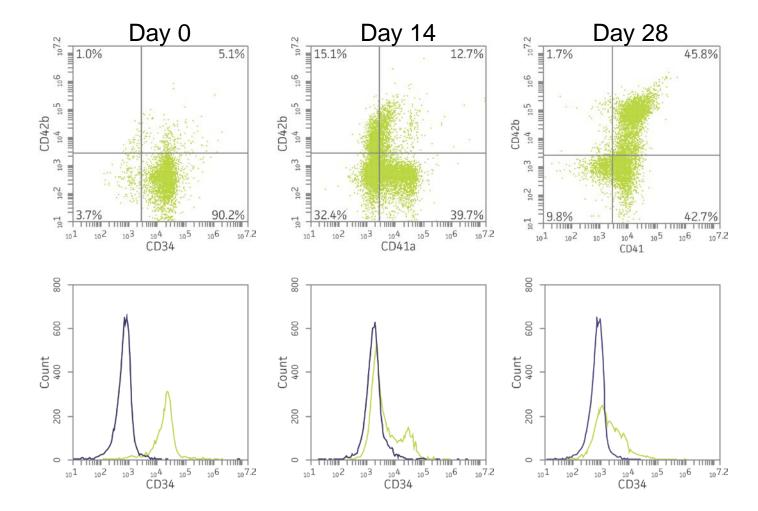


# Expression of megakaryocyte lineage markers on differentiated bone marrow CD34+ HPSCs



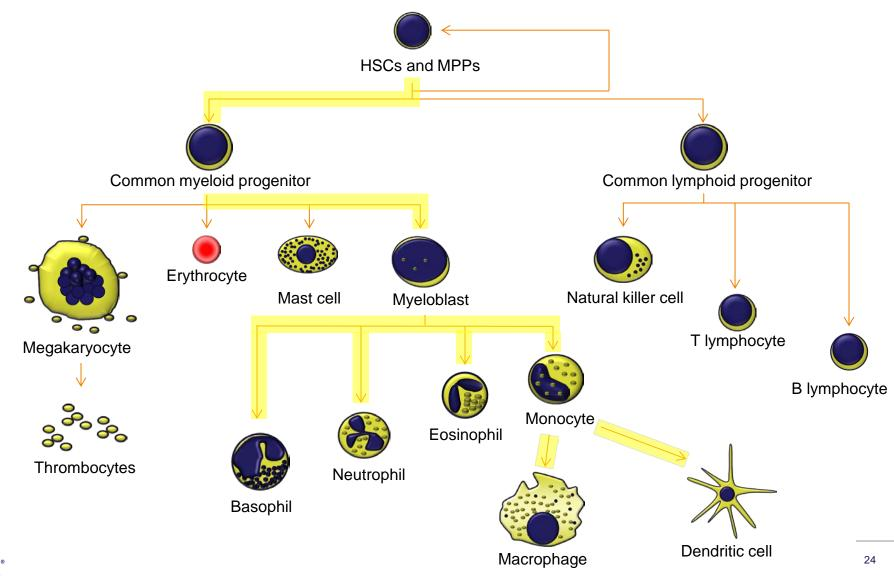


# Expression of megakaryocyte lineage markers on differentiated cord blood CD34+ HPSCs



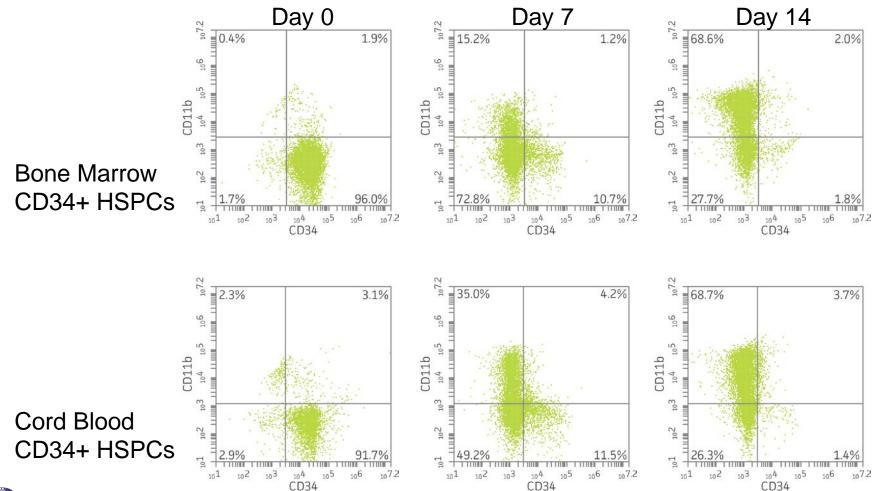


## Pan-myeloid differentiation and expansion





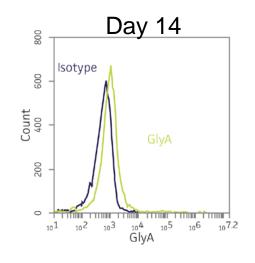
# Expression of pan-myeloid lineage markers on differentiated BM and CB CD34+ HPSCs

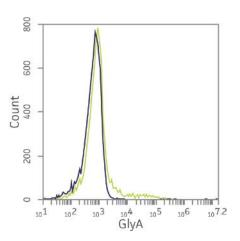


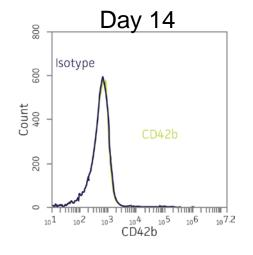


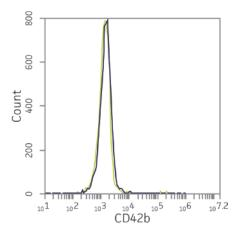
# Specificity of pan-myeloid-directed differentiation of CD34+ HPSCs

Bone Marrow CD34+ HSPCs





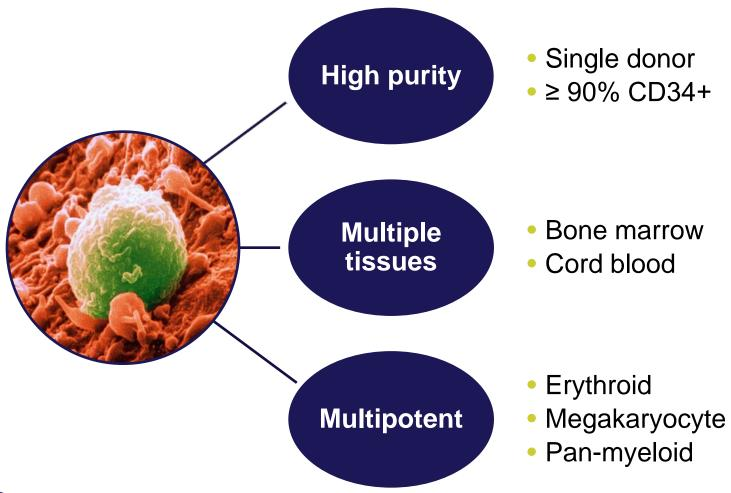




Cord Blood CD34+ HSPCs



# **ATCC primary CD34+ HSPCs: Summary**

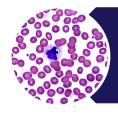




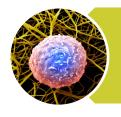
### **Outline**



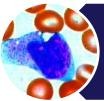
## Background



CD34+ hematopoietic stem & progenitor cells



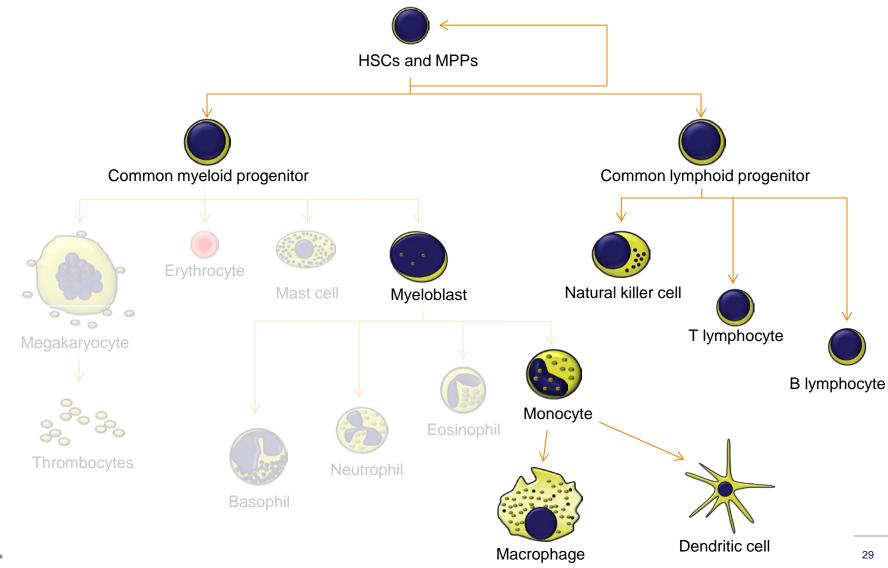
Mononuclear cells



CD14+ monocytes



### **Mononuclear Cells**





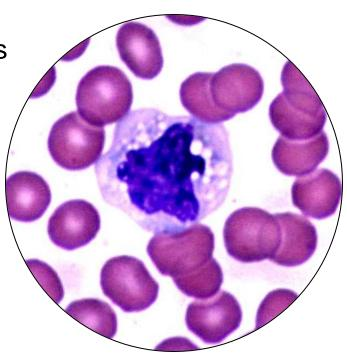
# Primary mononuclear cells

### **Applications**

- Isolation and study of cell subpopulations
- Molecular expression profiling

### Key research areas

- Infectious disease
- Blood pathologies
- Immunology
- Vaccine development
- Toxicology
- Regenerative medicine



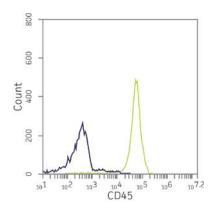
Monocyte and erythrocytes

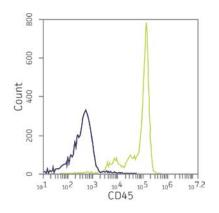


# Primary mononuclear cells

- Healthy human volunteer donors; IRB-approved informed consent
- Adult, non-pregnant
- Cryopreserved at P0; Purity: ≥ 90% CD45+
- Age, gender, ethnicity, and blood type on CoA

ATCC® No.	Tissue	Туре	Size
PCS-800-011	Peripheral blood	Mononuclear cells (PBMCs); Normal, Human	≥ 25 x 10 <sup>6</sup>
PCS-800-013	Bone marrow	Mononuclear cells (BMMCs); Normal, Human	≥ 25 x 10 <sup>6</sup>







Bone marrow mononuclear cells

# Primary mononuclear cells

#### Reported on CoA (lot specific)

Specific marker expression

#### **Peripheral Blood**

- CD45+ (%)
- CD3+ (%)
- CD8+ (%)
- CD14+ (%)
- CD19+ (%)
- CD56+ (%)

#### **Bone Marrow**

- CD45+ (%)
- CD3+ (%)
- CD8+ (%)
- CD14+ (%)
- CD19+ (%)
- CD34+ (%)
- CD56+ (%)



Human pathogenic virus testing

(by flow cytometry)

for HIV (VII), HepB, HepC, and HTLV (VII)

Characterization / call apecific staining

1	
	***************************************
ATCC (American Type Culture Collection)	800-838-8597 or 703-365-2700
P.O. Box 1549	Fax: 703-365-2750
Managaga, VA 20108 USA	E-mail: tech@atoc.org
www.aicc.org	or contact your local distributor
Pane 1 of	

Negative

CD45+: Positive (2 70%)

CO3+: Report results

CD4+: Report results

COS+: Report results

CO14+: Report results

CO19+: Report results

CO34+: Report results

CO56+: Report results

HIV (VII) - Negative HepS - Negative

HepC - Negative HTLV (UE) - Negati

CD45+: 98.25%

CD3+: 53.80%

CD4+: 44.77%

CD6+: 22.27%

CD14+: 9.26%

CD19+: 10.38%

CD34+: 6.15%

CD58\*: 7.40%



# Non-viral method increases natural killer cells' anti-cancer cell cytotoxicity

#### Enhanced Cytotoxicity of Natural Killer Cells following the Acquisition of Chimeric Antigen Receptors through Trogocytosis

Fu-Nan Cho<sup>1</sup>, Tsung-Hsien Chang<sup>2</sup>, Chih-Wen Shu<sup>2</sup>, Ming-Chin Ko<sup>3</sup>, Shuen-Kuei Liao<sup>4</sup>, Kang-Hsi Wu<sup>3</sup>, Ming-Sun Yu<sup>6</sup>, Shyh-Jer Lin<sup>6</sup>, Ying-Chung Hong<sup>6</sup>, Chien-Hsun Chen<sup>7</sup>, Chien-Hui Hung<sup>3</sup>, Yu-Hsiang Chang<sup>3,6,6</sup>

1 Department of Christinics and Gymecology, Calchisung Visterans General Hospick, Sachshaing, Taisow, 2 Department of Medical Education and Research, Kachshaing, Taisow, 2 Department of Pedicities, Cachshaing, Taisow, 2 Department of Medicities, Cachshaing, Taisow, 2 De

#### Abstract

Natural killer (NO) cells have the capacity to target tumors and are ideal candidates for immunotherapy. Viral vectors have been used to generically modify in vitroe expanded Nic cells to express chimeric entiripes mecapros (CARI), which competed the control of the control of

Citation: Cho F-N, Chang T-H, Shu C-W, Ko M-C, Liao S-K, et al. (2014) Enhanced Cytotoxicity of Natural Killer Cells following the Acquisition of Chimeric Antigen Receptors through Trogocytoxis. PLoS ONE 9(10): e109352. doi:10.1371/journal.pone.0109352

Editor: Jacques Zimmer, Centre de Recherche Public de la Santé (CRP-Santé), Luxembourg

Received April 25, 2014; Accepted September 4, 2014; Published October 14, 2014

Copyright: © 2014 Cho et al. This is an open-access article distributed under the terms of the Creative Commons Attribution License, which permits unrestricted use, distribution, and reproduction in any medium, provided the original author and source are credited.

Data Availability: The authors confirm that all data underlying the findings are fully available without restriction. All relevant data are wi

Fundings: This study was supported by grann from National Health Research Institutes (95680), http://english.nhi.org.tu/hHRJLV88/hhrise001Action.doi, and professional transfer of the Study Stu

Compation Interests: The authors have declared that on connection interests exist

\* Email: yhchang@vghks.gov.tw

These authors contributed equally to this work

#### Introduction

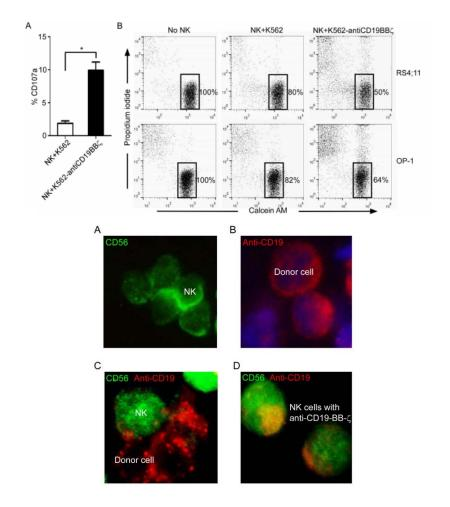
Natural killer (NK) cells have the ability to recognize and diminate numor cells, making them ideal candidates for tumor immunotherapy [1,2]. NK cell activity is regulated by the rumulative effects of multiple activating and inhibitory signals that are transmitted through the receptors on the NK cell surface. We have previously genetically modified in wire expanded NK cells to express DAP10 and the chimeric NKG2D receptor rotataining the CD35, signal domain, which altered the balance between the activating and inhibitory signals of NK cells and ruhanced the cytotoxicity against NKC2D ligand-bearing tumors [3], Further, expression of anti-CD19 chimeric antigen receptors CARs) containing 41BB and CD35, signal domains on NK cells ruhanced the activating signals originating from CD19 antigen

engagement, leading to cytotoxicity specifically against B-cell leukemia [4].

Trogosytosis is a process in which membrane patches are exchanged between target and immune cells [3–7]. When an NK cell interacts with a target cell, an immune synapse, which is strong enough to allow the transfer of small membrane patches from one cell to its partner cell, is formed [8,9]. Therefore, target cell suffice molecules can be found on the surface of NK cells. The chemokine receptor CCR7 has been shown to be transferred from donor cells onto the surface of NK cells via trogocytosis, and this transfer stimulated NK cell inguistics, Icading to evaluate of lymph node homing [10,11]. Similarly, T cells captured NKG2D and NKpd6 [ligands on timore cell through trogosytosis and promoted NK cell.

PLOS ONE | www.plosone.org

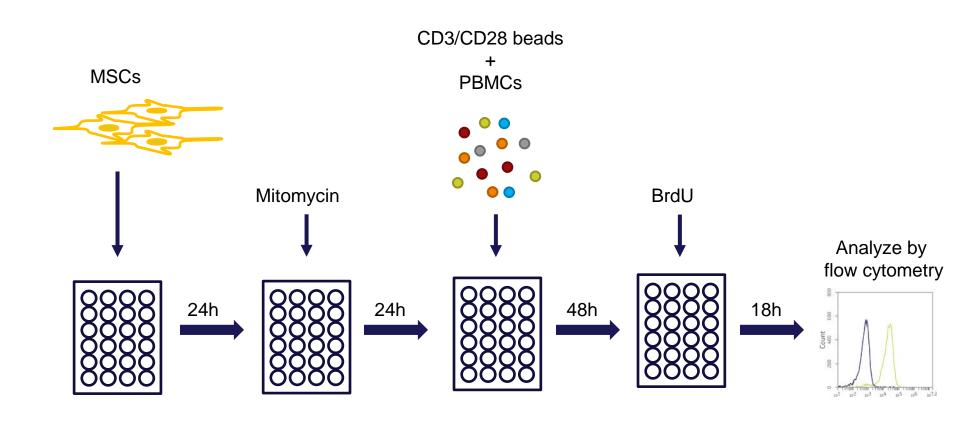
October 2014 | Volume 9 | Issue 10 | e109352





Cho, Fu-Nan, et al. PloS one 9.10 (2014): e109352.

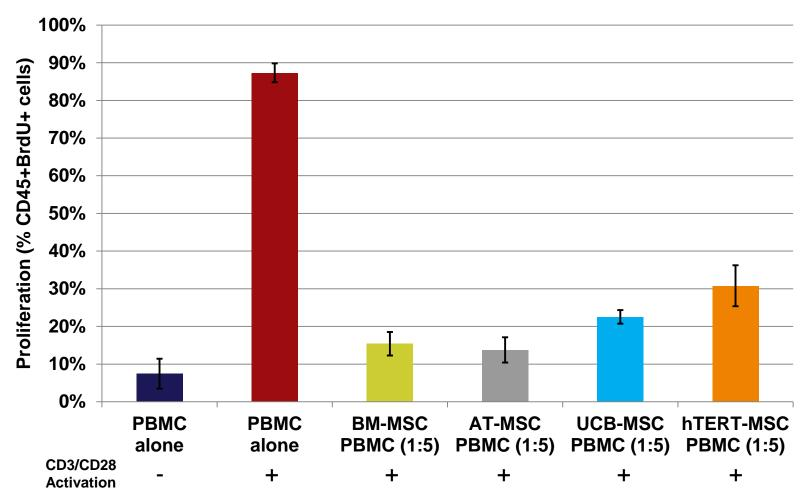
# Immunosuppressive assay using PBMCs



For detailed differentiation protocols see the ATCC website



# MSCs suppress activated T-cell proliferation





# **ATCC Mesenchymal Stem Cells**

The complete study, presented at ISSCR 2014, is available on the ATCC website:

Comparative analysis of cell proliferation, immunosuppressive action, and multi-lineage differentiation of immortalized MSC and MSC from bone marrow, adipose tissue, and umbilical cord blood



Dezhong Yin, Ph.D., Joy A. Wells, James Clinton, Ph.D. and Chaozhong Zou, Ph.D. *ATCC Cell Systems*, 22 Firstfield Rd, Suite 180, Gaithersburg, MD 20878, USA

ISSCR Poster #: F-3115

#### For more information on our MSC products: <a href="www.atcc.org/stemcells">www.atcc.org/stemcells</a>

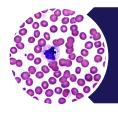
ATCC® No.	Туре	Description
PCS-500-010	Primary	Umbilical Cord-derived Mesenchymal Stem Cells
PCS-500-011	Primary	Adipose-derived Mesenchymal Stem Cells
PCS-500-012	Primary	Bone Marrow-derived Mesenchymal Stem Cells
SCRC-4000	Immortalized	hTERT Immortalized Adipose-derived Mesenchymal Stem Cells



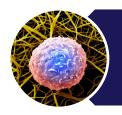
### **Outline**



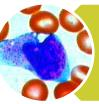
## Background



CD34+ hematopoietic stem & progenitor cells



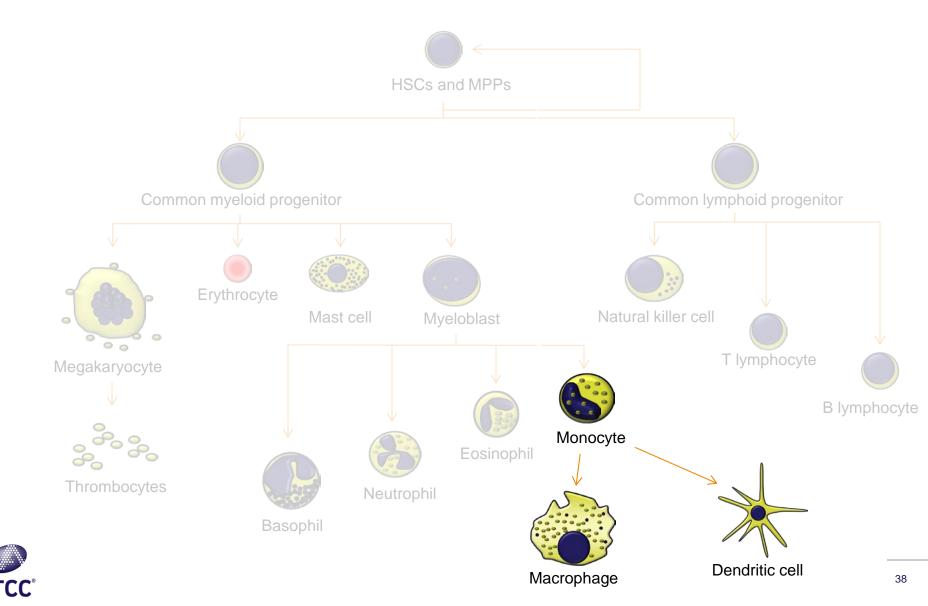
Mononuclear cells



CD14+ monocytes



# **Primary monocytes (CD14+)**



# **Primary CD14+ monocytes**

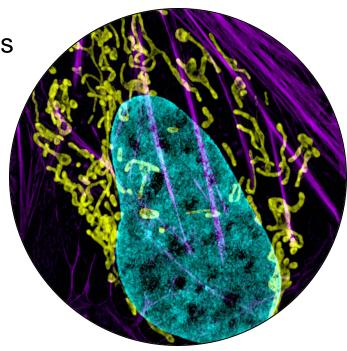
#### **Applications**

Isolation and study of monocyte subtypes

- Differentiation
- Phagocytosis
- Chemotaxis/migration assays

#### **Key research areas**

- Immunology
- Monocyte polarization
- Inflammation associated pathologies
- Infectious disease
- Cytokine release



Bone cancer cell, courtesy of Burnett Lippincott-Schwartz, National Cancer Institute



# Primary monocyte-derived macrophages are suitable for studying *Lm*, cell lines are not

OPEN & ACCESS Freely available online

PLOS ONE

#### CD14-Dependent Monocyte Isolation Enhances Phagocytosis of *Listeria monocytogenes* by Proinflammatory, GM-CSF-Derived Macrophages

Caroline Neu<sup>1</sup>, Anne Sedlag<sup>1</sup>, Carina Bayer<sup>2</sup>, Sabine Förster<sup>3</sup>, Peter Crauwels<sup>3</sup>, Jan-Hendrik Niess<sup>4</sup>, Ger van Zandbergen<sup>3</sup>, Giada Frascaroli<sup>2</sup>, Christian U. Riedel<sup>1</sup>\*

1 Institute of Microbiology and Biotechnology, University of Ulm, Ulm, Germany, 2 Institute of Virology, University Medical Center Ulm, Ulm, Germany, 3 Division of Immunology, Paul-Ehrlich-Institute, Fuderal Institute for Vaccines and Biomedicines, Langen, Germany, 4 Department of Visceral Medicine and Surgey, Insetspital, Bern, Seitzvatand

#### Abstract

Macrophages are an important line of defence against invading pathogens, Human macrophages derived by different methods were tested for theis visualishily as models to investigate interior amongospense (Inr.) infection and compared to macrophage-like THP-1 cells, Human primary monocytes were isolated by either positive or negative immunomagnetic selection and differentiated in the presence of granulocyte macrophage colony-stimulating factor (Ind.CSF) or macrophage colony-stimulating factor (Ind.CSF) or macrophage (Ind.Med) stained positive for C206s and McSF-derived macrophages (Ind.Med) for CD163. THP-1 cells did not express CD360 or CD163 following incubation with PMA, M- or GM-CSF alone or in combination (CD163. THP-1 cells did not express CD360 or CD163 following incubation with PMA, M- or GM-CSF alone or in combination of the PMF or CD163. THP-1 cells did not express CD360 or CD163 following incubation with PMA, M- or GM-CSF alone or in combination of the PMF or CD163. THP-1 was severely reduced even at lower bacterial numbers. M-Meg generally showed high phagocytosis of LM (SM-Mey was markedly influenced by treatmed used for isolation of monocytes, GM-Mey derived from negatively isolated monocytes showed pool support of the Medical monocytes of LM (Medical CM) and the Medical monocytes and the protocoan parasite Leishmania major by GM-Mey was not inducating that this effect is specific for LM. Based on these observations we propose macrophages derived by ex who differentiation of negatively selected human primary monocytes as the most suitable model to study In infection of macrophages.

Citation: Neu C, Sedlag A, Bayer C, Förster S, Crauwels P, et al. (2013) CD14-Dependent Monocyte Isolation Enhances Phagocytosis of Listeria monocytogenes by Proinflammatory, GM-CSF-Durived Macrophages. PLoS ONE 8(6): e66898. doi:10.1371/journal.pone.0066898

Editor: Jörn Coers, Duke University Medical Center, United States of America Received March 7, 2013; Accepted May 13, 2013; Published June 11, 2013

Copyright: © 2013 Neu et al. This is an open-access article distributed under the terms of the Creative Commons Attribution License, which permits unrestricted use, distribution, and reproduction in any medium, provided the original author and source are credited.

Funding: This study was partially funded by a grant of the Carl Zeiss Foundation in the "Programm zur 3 Stärkung von Forschungsstrukturen an Universitäten" to GF, GVZ and CUB. The funders had no role in study design, data collection and analysis, decision to publish, or preparation of the manuscript. No additional actual funding mass received for this crudy.

Competing Interests: The authors have declared that no competing interests exist.

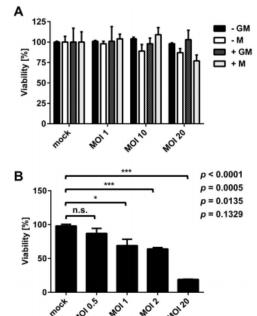
#### \* E-mail: christian.riedel@uni-ulm.de

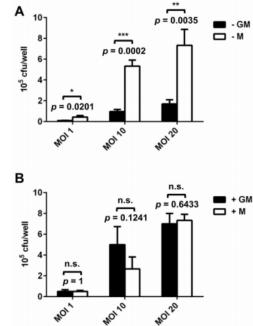
#### Introduction

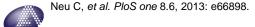
Littrie sossoptopue (La) is a food-borne Gram-positive obligate intracellular pathogen that is able to cope with a wide range of environmental conditions and is thus found in various of labilists [1]. In humans, the disease caused by L in is termed listeriosis and manifests primarily in immunocompromised individuals, pregnant women, new-borns, and elderly patients with a mortality of 20–30% in these art risk groups [2]. Infections with L are usually acquired upon consumption of contaminated food products and thus the first habitat inside the host is the gastroinestical tract [3]. L m is able to cross the intestinal barrier, subsequently enters the blood and lymph stream, and finally colonizes liver and splicen where it is primarily phagocytosed by resident unacrophages [4].

the secretion of two phospholipases,  $\Re C \lambda$  and  $\Re C R_i$ , and the poreforming toxin listeriolysin O (LLO) [5]. This results in the release of L m into the cytoplasm where it starts to replicate and spread from one cell to another by hijacking the host cell actin cytoskeleton [6].

Macrophages play a central role in activating and finely balancing the proc and anti-fullammatory pathways of the host immune system to mount effective host responses against invading pathogens. In vise, macrophage differentiation is driven by GMand McSF [7,8]. High levels of GMcSF induce a proinflammatory phenotype resulting in high II.-12 secretion. These pro-inflammatory cells are also termed MI macrophages. By contrast, McSF polarizes macrophages to an anti-inflammatory phenotype characterized by III-10 secretion, which is referred to



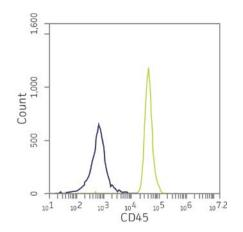


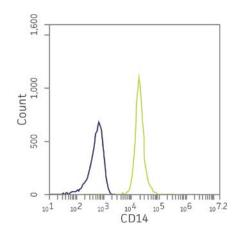


# **ATCC** primary CD14+ monocytes

- Healthy human volunteer donors; IRB-approved informed consent
- Adult, non-pregnant
- Cryopreserved at P0; Purity: ≥ 90% CD14+
- Age, gender, ethnicity, and blood type on CoA

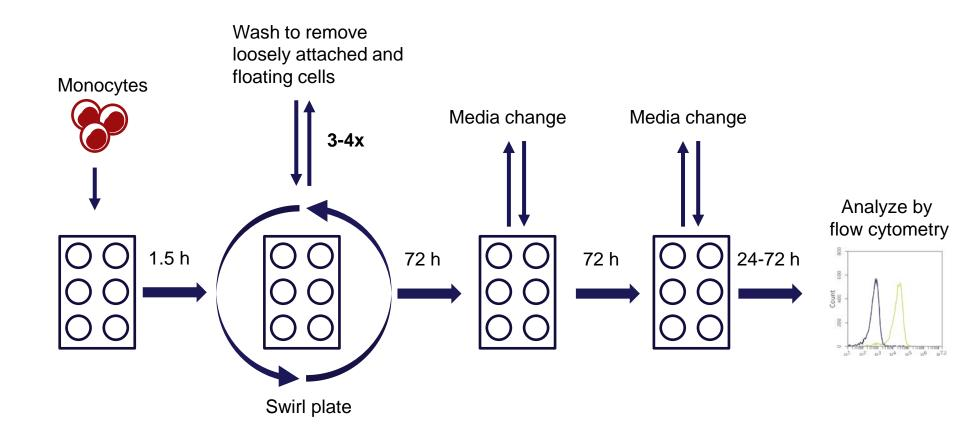
ATCC® No.	Tissue	Туре	Size
PCS-800-010	Peripheral blood	Monocytes (CD14+)	≥ 50 x 10 <sup>6</sup>







# Macrophage differentiation protocol



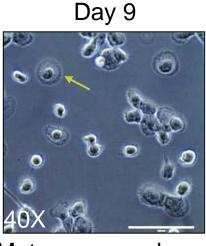


For detailed differentiation protocols see the ATCC website

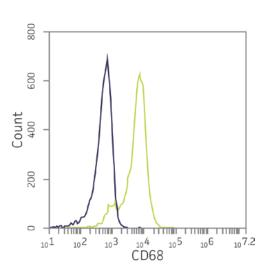
# Generation of CD68+ macrophages from monocytes

Day 1

Monocytes



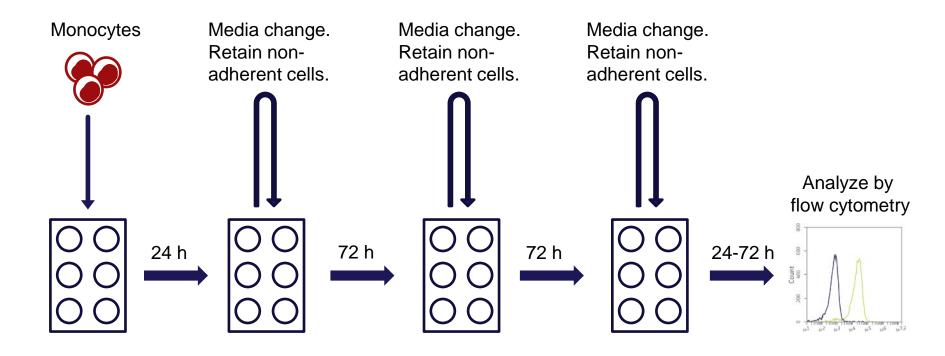
Mature macrophage



- Macrophages differentiated from CD14+ monocytes for 9 days.
- Morphology characteristic of type M1-polarized macrophages.
- > 80% of cells were CD68+



# Dendritic cell differentiation protocol



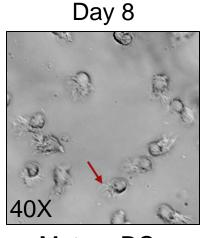


For detailed differentiation protocols see the ATCC website

# **Generation of CD83+ DCs from monocytes**

Day 1

Day 6

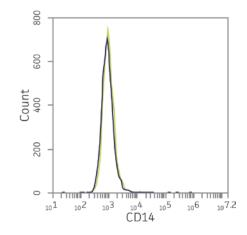


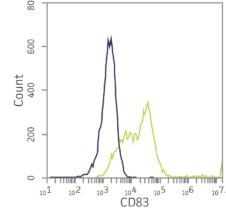
Monocytes

Immature DCs

Mature DCs

- After 8 days differentiation 95% of non-adherent cells were CD14-
- >70% of cells were CD83+







# **Summary**

ATCC offers a variety of well-characterized and functionally validated primary hematopoietic cell types

- CD34+ HSPCs
- BMMCs and PBMCs
- CD14+ monocytes

ATCC provides hematopoietic lineage-specific differentiation protocols

- Erythroid
- Megakaryocyte
- Non-specific myeloid
- Dendritic
- Macrophage

ATCC hematopoietic cells are useful in numerous areas of research



# Thank you!

Register for more webinars in the ATCC "Excellence in Research" webinar series at www.atcc.org/webinars.



March 5, 2015
10:00 AM, 3:00 PM ET
David Clawson, Lead Biologist, ATCC
Enhancing vector-borne research with biological and molecular standards



March 19, 2015
10:00 AM, 3:00 PM ET
Scott Sutton, Ph.D., *Principal*, Microbiology Network and
Tracy Vandenbroek, *Product Line Business Manager*, ATCC
Microbiology quality control as described in the Compendia

Thank you for joining today!

Please send additional questions to <a href="mailto:tech@atcc.org">tech@atcc.org</a>

