

Generation of Monocytes from Human Pluripotent Stem Cells Using STEMdiff™ Medium and Supplements

Background

Monocytes are essential components of the innate immune system that provide defense against pathogens or tumors due to their ability to differentiate to macrophages and dendritic cells. While monocytes can be isolated from peripheral blood, human pluripotent stem cells (hPSCs) offer another, potentially unlimited, source of monocytes. These hPSC-derived monocytes can be used in disease modeling, development of cell therapy applications, and research into basic biology.

STEMdiff™ Monocyte Kit facilitates the differentiation of hPSCs to monocytes under feeder-free and serum-free culture conditions.

Differentiate hPSCs to Monocytes

STEMdiff™ Monocyte Kit is used to differentiate hPSCs to monocytes in a three-stage protocol as shown in Figure 1. In stage 1, Medium A differentiates hPSCs to the mesoderm lineage after 3 days of culture. In stage 2, the cells are cultured in Medium B for 4 days, promoting their specification to hematopoietic progenitor cells. During stage 3, Monocyte Differentiation Medium is used to promote the differentiation of progenitor cells to monocytes. These can be identified by the expression of CD14, as shown in Figure 2, and can be repeatedly harvested. STEMdiff™ Monocyte Kit is optimized for differentiation of multiple embryonic stem (ES) and induced pluripotent stem (iPS) cell lines maintained in TeSR™ PSC maintenance medium. After differentiation, hPSC-derived monocytes may be used in additional downstream assays.

Why Use STEMdiff™ for Generating Monocytes?

EFFICIENT. Generate up to 7 million CD14⁺ monocytes per plate in just 14 - 23 days.

CONSISTENT. Eliminate variation introduced by serum and feeder-cells by using serum- and feeder-free conditions.

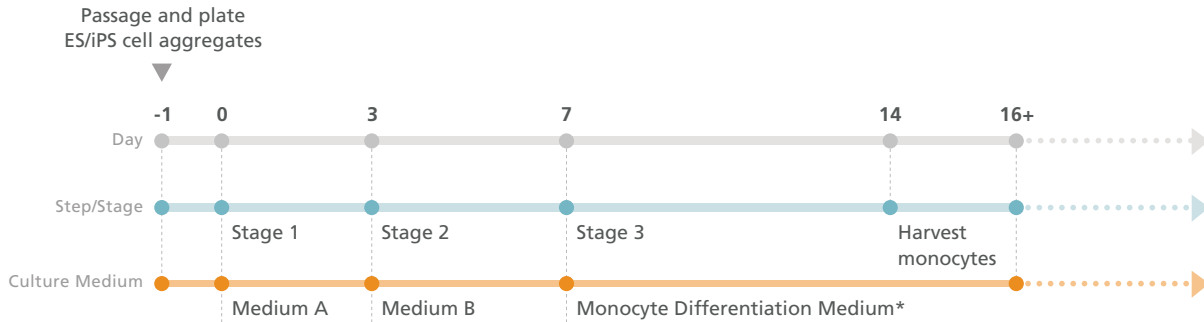
EASY-TO-USE. Produce monocytes in a simple monolayer culture for easier harvest of suspended cells.

FLEXIBLE. Achieve robust generation of monocytes across multiple ES and iPS cell lines.

Product Name	Catalog # (Size)
STEMdiff™ Monocyte Kit	05320 (1 Kit)
Kit Components	Component # (Size)
STEMdiff™ Hematopoietic Basal Medium	05311 (120 mL)
STEMdiff™ Hematopoietic Supplement A (200X)	05312 (225 µL)
STEMdiff™ Hematopoietic Supplement B (200X)	05323 (225 µL)
STEMdiff™ Monocyte Differentiation Supplement (100X)	05324 (3 x 1 mL)
StemSpan™ SFEM II	09605 (3 x 100 mL)

Protocol for Differentiation of hPSCs to Monocytes

This protocol is designed to promote the differentiation of hPSCs to monocytes over 17 - 23 days of culture in a 2-dimensional (2D) culture system. The three stages — mesoderm formation (stage 1), hematopoietic specification (stage 2), and monocyte differentiation (stage 3) — are shown in Figure 1.



*Perform a full medium change every 2 - 3 days as needed.

Figure 1. Monocyte Differentiation Protocol

One day prior to differentiation, hPSC colonies are harvested and seeded as small aggregates (100 - 200 µm in diameter) at 10 - 20 aggregates/cm² in mTeSR™1, TeSR™-E8™, or mTeSR™ Plus. After one day, the medium is replaced with Medium A (STEMdiff™ Hematopoietic Basal Medium + Supplement A) to induce mesodermal specification (stage 1). On day 3, the medium is changed to Medium B (STEMdiff™ Hematopoietic Basal Medium + Supplement B) to promote hematopoietic specification (stage 2). On day 7, the medium is replaced with Monocyte Differentiation Medium (StemSpan™ SFEM II + STEMdiff™ Monocyte Differentiation Supplement) to promote the production of CD14⁺ monocytes (stage 3). Monocyte Differentiation Medium is used for all medium changes for the remaining culture period. CD14⁺ cells can be detected in suspension starting after day 14, and their frequency gradually increases until day 17 - 23. CD14⁺ cells can be harvested directly from the culture supernatant during medium changes.

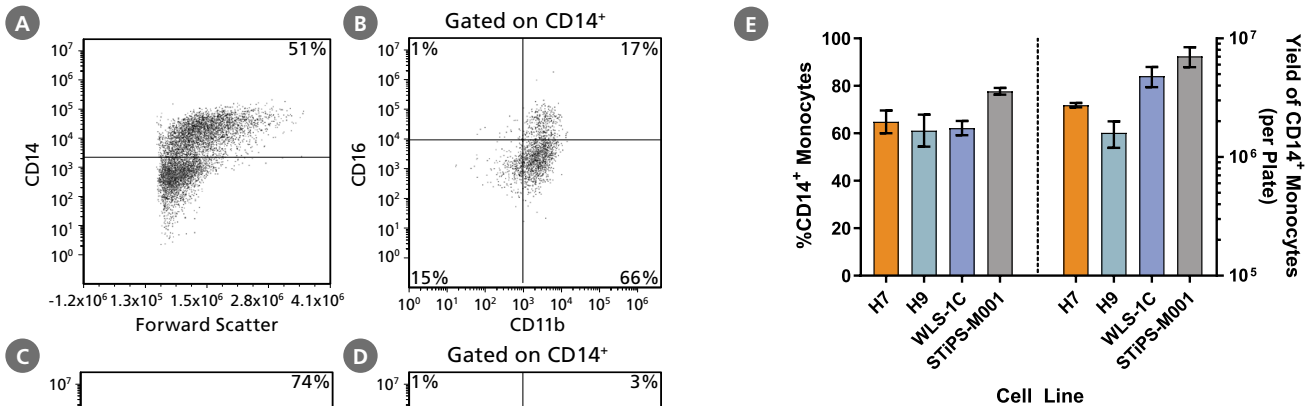


Figure 2. Robust and Efficient Generation of CD14⁺ Monocytes Using STEMdiff™ Monocyte Kit

hPSCs were differentiated to monocytes using the 2D culture system described in Figure 1. Between days 17 and 23, cells were harvested every 2 - 3 days and analyzed by flow cytometry for CD14 expression. Representative flow cytometry plots are shown for (A,B) iPS (WLS-1C)-derived cells and (C,D) ES (H9)-derived cells. (E) The average frequency of viable CD14⁺ monocytes at the peak harvest was 61 - 78%. The average yield of CD14⁺ monocytes produced per 6-well plate at the peak harvest was between 1.6 x 10⁶ and 7.1 x 10⁶ cells. Data are shown as mean ± SEM (n = 3 - 14).

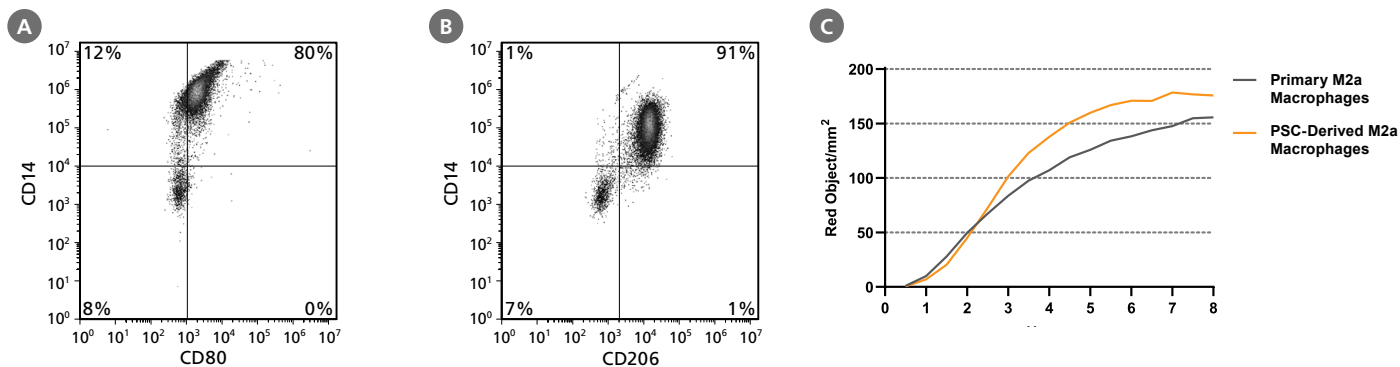


Figure 3. STEMdiff™ Monocyte Kit Generates Monocytes That Are Capable of Differentiation to Macrophages

hPSC-derived monocytes were harvested after 21 days of culture. These were then differentiated to macrophages using ImmunoCult™-SF Macrophage Medium with 100 ng/mL M-CSF for 4 days. Macrophages were then incubated for an additional 2 days with either 10 ng/mL of LPS and 50 ng/mL of IFN- γ , or 10 ng/mL IL-4, to become polarized to M1 or M2a macrophages, respectively. Representative flow cytometry plots of (A) M1 and (B) M2a macrophages produced from the WLS-1C iPS cell line are shown. (C) To measure phagocytosis, hPSC-derived M2a macrophages and peripheral blood (PB) monocyte-derived M2a macrophages (primary M2a macrophages), were incubated with pHrodo™ Red Zymosan A BioParticles® Conjugate and incubated at 37°C for 8 hours. Images were acquired using the IncuCyte® ZOOM every 30 minutes and analyzed for internalization of pHrodo™ Red Zymosan A BioParticles® (measured as red object/mm²). hPSC-derived and primary M2a macrophages show similar phagocytic activity.

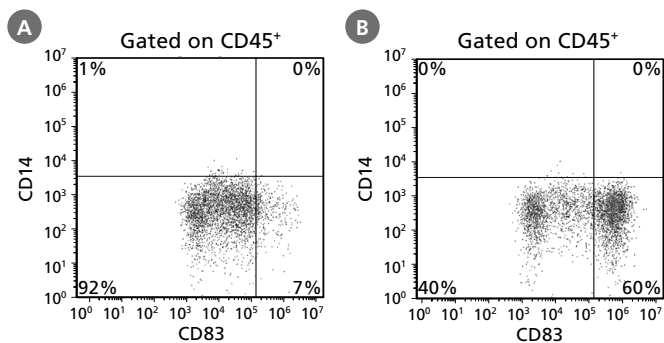


Figure 4. STEMdiff™ Monocyte Kit Generates Monocytes That Can Be Differentiated to Dendritic Cells

hPSCs were differentiated to monocytes, harvested after 21 days, and differentiated to dendritic cells using ImmunoCult™ Dendritic Cell Culture Kit. Half of the dendritic cells were harvested on day 7 and examined for CD14 and CD83 expression to identify CD14⁺CD83⁻ immature dendritic cells. The remaining dendritic cells were activated for 2 days and assessed for the presence of CD14⁺CD83⁺ mature dendritic cells at day 7. Representative cultures initiated with ES (H9) cells are shown for production of (A) immature dendritic cells and (B) mature dendritic cells.

Product Information

Materials Required But Not Included

Product Name	Size	Catalog #
mTeSR™1	500 mL Kit	85850
	1 L Kit	85857
	10 Kits	85870
	25 Kits	85875
OR		
TeSR™-E8™	1 Kit	05990
OR		
mTeSR™ Plus	1 Kit	05825

Product Name	Size	Catalog #
Gentle Cell Dissociation Reagent	100 mL	07174
OR		
ReLeSR™	100 mL	05872
OR		
Dispase (1 U/mL)	100 mL	07923

Product Name	Size	Catalog #
DMEM/F-12 with 15 mM HEPES	500 mL	36254
6-Well Flat-Bottom Plate, Tissue Culture-Treated	50 Plates	38015
96-Well Flat-Bottom Plate, Non-Treated	50 Plates	38044
Human Recombinant IFN-γ	100 µg	78020
Human Recombinant IL-4	100 µg	78045

Materials Recommended for Further Differentiation

Product Name	Size	Catalog #
ImmunoCult™-SF Macrophage Medium	250 mL	10961
Human Recombinant M-CSF	100 µg	78057

Product Name	Size	Catalog #
ImmunoCult™ Dendritic Cell Culture Kit	1 Kit	10985
Kit Components		
ImmunoCult™-ACF Dendritic Cell Medium	100 mL	10987
	500 mL	10986
ImmunoCult™-ACF Dendritic Cell Differentiation Supplement	1 mL	10988
ImmunoCult™ Dendritic Cell Maturation Supplement	0.5 mL	10989

Recommended Antibodies for Analysis

Product Name	Catalog #
Anti-Human CD14 Antibody, Clone M5E2	60004PE
Anti-Human CD16 Antibody, Clone 3G8	60041FI
Anti-Human CD11b Antibody, Clone ICRF44	60040AZ
Anti-Human CD83 Antibody, Clone HB15e	60107PE

Copyright © 2020 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies & Design, STEMCELL Shield Design, Scientists Helping Scientists, STEMdiff, ImmunoCult, ReLeSR, and StemSpan are trademarks of STEMCELL Technologies Canada Inc. IncuCyte is a registered trademark of Sartorius. pHrodo™ and BioParticles® are trademarks of Thermo Fisher Scientific. mTeSR, TeSR, and E8 are trademarks of WARF. All other trademarks are the property of their respective holders. While STEMCELL has made all reasonable efforts to ensure that the information provided by STEMCELL and its suppliers is correct, it makes no warranties or representations as to the accuracy or completeness of such information.

PRODUCTS ARE FOR RESEARCH USE ONLY AND NOT INTENDED FOR HUMAN OR ANIMAL DIAGNOSTIC OR THERAPEUTIC USES UNLESS OTHERWISE STATED. FOR ADDITIONAL INFORMATION ON QUALITY AT STEMCELL, REFER TO WWW.STEMCELL.COM/COMPLIANCE.