

Human Primary Cells

For Cell-Based Assays



Table of Contents

- 3 Introduction
- 4 Human Bone Marrow (BM) Products
- 5 Human Peripheral Blood (PB) Products
- 6 Human Umbilical Cord Blood (CB) Products
- 7 Other Primary Cell Types
- 8 Cell Separation
- 9 Contract Assay Services
- 10 Cell Culture Media and Reagents

STEMCELL Technologies Inc. is a biotechnology company that develops specialty cell culture media, cell separation products and ancillary reagents for life science research. Driven by science and a passion for quality, STEMCELL Technologies delivers over 1000 products to more than 70 countries worldwide. To learn more about how STEMCELL Technologies helps to make research work, visit www.stemcell.com.

Introduction

Human primary cells are invaluable components of cell-based assays. They are preferred over transformed or immortalized cell lines because they are more representative of cells in vivo.

STEMCELL Technologies offers a wide range of frozen (unprocessed and purified) human primary cells.* Primary hematopoietic cells are available from bone marrow (BM), peripheral blood (PB) and umbilical cord blood (CB). Marrow stromal cells are also available.

Benefits

REPRODUCIBILITY. Multiple vials of a single lot of cells can be requested, allowing the same cells to be used over a prolonged period of time.

FLEXIBILITY. Frozen cells allow experiments to begin at any time compounds are available for testing.

REDUCED COSTS. Save costs associated with collecting and culturing primary cells, including time, reagents and consumables.

RESOURCE ALLOCATION. The purchase of frozen cells allows reallocation of resources to other challenging projects.

NOTE: Treat all cell products as potentially contaminated biological specimens, even if available serological reports are negative. All human cell products should be handled at Biosafety Level 2 or higher.

STEMCELL Technologies is the exclusive worldwide distributor for AllCells™, LLC (excluding China & Japan), with co-distribution rights in North America.

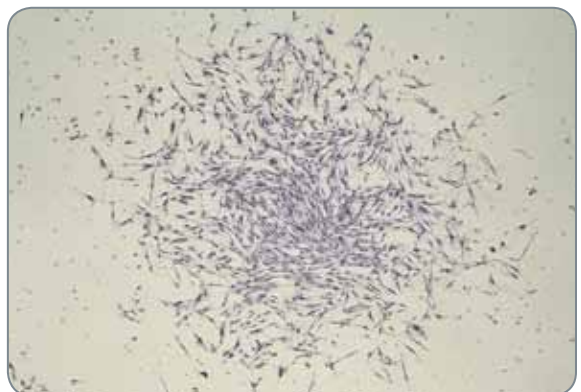
*Primary cell products are for in vitro research use only. Not approved for diagnostic, therapeutic, or clinical applications. Not approved for human or veterinary use in vivo.



Hematopoietic cells cultured in StemSpan™ serum-free expansion medium.



Colony derived from a human hematopoietic progenitor cell using MethoCult™ methylcellulose-based medium.



CFU-F derived from human bone marrow cells using MesenCult™.

Human Bone Marrow (BM) Products

Primary human cells are isolated from adult bone marrow (BM) drawn from the posterior iliac crest, 25 mL/site, from a maximum of four sites per donor. BM is diluted with PBS containing heparin (150 U/mL of BM). Mononuclear cells (MNC) are obtained by light density separation (1:2 dilution in PBS and centrifugation over Ficoll™). Lineage-specific cells are then positively or negatively selected using immunomagnetic cell separation procedures. All cell preparations are cryopreserved.

BM Mononuclear Cells (MNC)

DESCRIPTION	QUANTITY	CATALOG #
Frozen BM MNC	5 million	ABM007F
	15 million	ABM008F
	25 million	ABM009F
	50 million	ABM010F
	100 million	ABM011F

BM Stem/Progenitor Cells

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
CD34 ⁺ CELLS			
Frozen BM CD34 ⁺ Cells	Positive	0.1 million	ABM015F
		0.3 million	ABM021F
		0.5 million	ABM022F
		1 million	ABM017F
		2 million	ABM012F
		5 million	ABM018F
		10 million	ABM019F
CD133 ⁺ CELLS			
Frozen BM CD133 ⁺ Cells	Positive	0.5 million	ABM026F

Positive: cells isolated by positive selection using immunomagnetic isolation procedures.

Lineage-Selected BM Cells

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
EARLY ERYTHROID PROGENITOR CELLS			
Frozen BM CD36 ⁺ Cells	Positive	1 million	ABM020F
ENDOTHELIAL CELLS			
Frozen BM CD105 ⁺ Cells	Positive	0.3 million	ABM028F
MYELOID CELLS			
Frozen BM CD33 ⁺ Cells	Positive	5 million	ABM032F
PLASMA CELLS			
Frozen BM CD138 ⁺ Plasma Cell Pellet	Positive	1 million	ABM029P
CD34 ⁺ DEPLETED CELL PELLETT*			
Frozen BM CD34 Depleted MNC	Negative	100 million	ABM023F

Positive: cells isolated by positive selection using immunomagnetic isolation procedures.
 Negative: cells isolated by negative selection using immunomagnetic isolation procedures.
 *Isolated cells are pelleted and snap frozen. Pelleted cells are not viable and cannot be cultured.

Marrow Stromal Cells

First passage marrow stromal cells are obtained by culturing BM MNCs for 10 - 14 days in a modified medium containing prescreened FBS. These marrow stromal cells have a fibroblast-like morphology and express differentiation markers CD105 (>90%), CD166 (>90%), CD29 (>90%) and CD44 (>90%). They lack expression of CD14 (<1%), CD34 (<1%) and CD45 (<1%). While these marrow stromal cells may exhibit a mesenchymal stem cell phenotype, they have not been functionally tested for multipotent differentiation capabilities.

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
Frozen Marrow Stromal Cells	Culture	0.75 million	MSC-001F

Human Peripheral Blood (PB) Products

Primary human cells are isolated from adult peripheral blood (PB) using a COBE Spectra™ apheresis machine. The resulting cell preparation, commonly called a “leukapheresis,” contains primarily mononuclear cells (MNC). The cells are treated with ammonium chloride to lyse red blood cells (RBCs), and lineage-specific cells are then positively or negatively selected using immunomagnetic cell separation procedures. All cell preparations are cryopreserved.

PB Mononuclear Cells (MNC)

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
Frozen PB MNC	n/a	15 million	PB003F
		25 million	PB006F
		50 million	PB004F
		100 million	PB005F

PB Stem/Progenitor Cells

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
CD34 ⁺			
Frozen PB CD34 ⁺ Cells	Positive	0.2 million	PB033F-1
		0.5 million	PB033F-2
		1 million	PB033F-3

Positive: cells isolated by positive selection using immunomagnetic isolation procedures.

PB Plasma

DESCRIPTION	QUANTITY	CATALOG #
Frozen PB Plasma	10 mL	PB017F
	20 mL	PB018F
	30 mL	PB019F
	40 mL	PB020F
	50 mL	PB021F
	100 mL	PB022F
	150 mL	PB023F

Positive: cells isolated by positive selection using immunomagnetic isolation procedures
 Negative: cells isolated by negative selection using immunomagnetic isolation procedures
 *Monocyte-derived CD14⁺ dendritic cells (DC) are cultured with 10% FBS in the presence of GM-CSF and IL-4 for 4 - 5 days. The cultured cells are assessed by flow cytometric analysis and found to express immature DC antigens including CD1a (>90%), CD11c (>90%), CD40 (>90%), CD86 (>90%), and lack the expression of CD14 (<5%).
 **Mononuclear cells from the same DC donor that have not been isolated and purified into dendritic cells.

Lineage-Selected PB Cells

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
T CELLS			
Frozen PB Pan T Cells	Negative	20 million	PB009-1F
Frozen PB CD4 ⁺ Cells	Negative	15 million	PB009-2F
Frozen PB CD8 ⁺ Cells	Negative	10 million	PB009-3F
Frozen PB CD4 ⁺ /CD25 ⁺ T Cells	Negative/ Positive	2 million	PB009-4F
Frozen PB CD4 ⁺ /CD45RA ⁺ Naive T Cells	Negative/ Positive	5 million	PB009-5F
Frozen PB CD8 ⁺ /CD45RA ⁺ Naive T Cells	Negative	5 million	PB009-6F
Frozen PB CD4 ⁺ /CD45RO ⁺ Memory T Cells	Negative	5 million	PB009-7F
B CELLS			
Frozen PB B Cells	Negative	10 million	PB010F
Frozen PB CD19 ⁺ Cells	Positive	10 million	PB010-P-F
MONOCYTES			
Frozen PB Monocytes	Negative	10 million	PB011F
Frozen PB CD14 ⁺ Cells	Positive	10 million	PB011-P-F-1
		20 million	PB011-P-F-2
		40 million	PB011-P-F-3
NATURAL KILLER (NK) CELLS			
Frozen PB NK Cells	Negative	5 million	PB012F
Frozen PB CD56 ⁺ Cells	Positive	5 million	PB012-P-F
LYMPHOID DENDRITIC CELLS (DC)			
Frozen PB BDCA4 ⁺ Cells	Positive	0.5 million	PB013F
MONOCYTE-DERIVED IMMATURE DENDRITIC CELLS (DC)			
Frozen PB Immature Dendritic Cells	Positive*	1.5 million	PB-DC001F
Frozen PB Immature Dendritic Cells HLA-A2 Screened	Positive*	1.5 million	PB-DC001-HLA-A2F
Frozen PB MNC (from Same DC Donor**)	n/a	10 million	PB-DC002F
		50 million	PB-DC003F
		100 million	PB-DC004F
		200 million	PB-DC005F
MONOCYTE-DERIVED MACROPHAGES			
Frozen PB Macrophages	Negative	1.5 million	PB-MDM-001F

Human Umbilical Cord Blood (CB) Products

Primary human cells are isolated from umbilical cord blood (CB). CB from a full-term delivery is collected into a 300 mL standard blood bag containing acid-citrate-dextrose (ACD) anti-coagulant (60 mL/bag). Mononuclear cells (MNC) are obtained by light density separation (1:2 dilution in PBS and centrifugation over Ficoll™). Residual red blood cells (RBCs) are then lysed using ammonium chloride, and lineage-specific cells are positively or negatively selected using immunomagnetic cell separation procedures. Samples are processed and cryopreserved within 24 hours of collection.

CB Mononuclear Cells (MNC)

DESCRIPTION	QUANTITY	CATALOG #
Frozen CB MNC	15 million	CB002F
	50 million	CB004F
	150 million	CB003F

Cord blood products may be from pooled donors, please specify on the purchase order if you require the product to be from a single donor.

CB Stem/Progenitor Cells

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
CD34⁺ CELLS			
Frozen CB CD34 ⁺ Cells (Mixed Donor)	Positive	0.2 million	CB007F
		0.5 million	CB005F
		1 million	CB008F
		5 million	CB009F
Frozen CB CD34 ⁺ Cells (Single Donor)	Positive	0.2 million	CB007F-S
		1 million	CB008F-S
CD133⁺ CELLS			
Frozen CB CD133 ⁺ Cells (Mixed Donor)	Positive	0.5 million	CB013F
Frozen CB CD133 ⁺ Cells (Single Donor)	Positive	0.5 million	CB013F-S

Positive: cells isolated by positive selection using immunomagnetic isolation procedures

CB Plasma

DESCRIPTION	QUANTITY	CATALOG #
Frozen CB Plasma	10 mL	CB027F
	20 mL	CB028F
	30 mL	CB029F
	40 mL	CB030F
	50 mL	CB031F
	100 mL	CB032F

Cord blood products may be from pooled donors, please specify on the purchase order if you require the product to be from a single donor.

Lineage-Selected CB Cells

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
T CELLS			
Frozen CB Pan T Cells	Negative	15 million	CB020-1F
Frozen CB CD4 ⁺ Cells		15 million	CB020-2F
Frozen CB CD8 ⁺ Cells		5 million	CB020-3F
Frozen CB Naïve CD4 ⁺ /CD45RA ⁺ T Cells		15 million	CB020-4F
B CELLS			
Frozen CB CD19 ⁺ Cells	Positive	1 million	CB019F
		2.5 million	CB020F
		5 million	CB021F
EARLY ERYTHROID PROGENITOR CELLS			
Frozen CB CD36 ⁺ Cells	Positive	1 million	CB010M-F
MONOCYTES			
Frozen CB CD14 ⁺ Cells	Positive	5 million	CB022F
NATURAL KILLER (NK) CELLS			
Frozen CB CD56 ⁺ Cells	Positive	1 million	CB023F
CD34 DEPLETED CELLS			
Frozen CB CD34 Depleted MNC	Negative	100 million	CB011F
CD133 DEPLETED CELLS			
Frozen CB CD133 Depleted MNC	Negative	100 million	CB014F

Positive: cells isolated by positive selection using immunomagnetic isolation procedures
 Negative: cells isolated by negative selection using immunomagnetic isolation procedures
 Cord blood products may be from pooled donors, please specify on the purchase order if you require the product to be from a single donor.

Other Primary Cell Types

Human Umbilical Vascular Endothelial Cells (HUVEC)

Human umbilical cord veins are processed to isolate endothelial cells via collagenase digestion and culture. HUVECs are cryopreserved at the end of a primary culture and can be cultured for up to 6 passages. HUVECs are characterized by morphological observation and flow cytometric analysis of differentiation markers, including Dil-Ac-LDL (LDL-uptake, >90%), CD31 (PECAM-1, >90%), CD54 (ICAM-1, >90%), CD62e (E-selectin, >90%) and von Willebrand factor (vWf, >90%).

DESCRIPTION	SELECTION	QUANTITY	CATALOG #
Frozen HUVEC	Culture	0.5 million	HUVEC-001F
Frozen HUVEC from mixed donors	Culture	0.5 million	HUVEC-001F-M

Customized Cell Types

In addition to the products listed, we offer other cell types from primary human hematopoietic sources, including products in custom volumes and additional donor testing. Please email us at techsupport@stemcell.com with your custom request.

DESCRIPTION	FORMAT	CATALOG #
Custom Request	Frozen Primary cells	ANYCELL-F

Product Warranty

STEMCELL Technologies warrants primary cell products meet the claimed product specifications, including viable cell number and purity, when the recommended protocols are followed. STEMCELL Technologies assures its cells to be viable and cell numbers recovered to be accurate, when handled exactly according to our instructions for thawing and counting on the product information sheets (available on request and on our website at www.stemcell.com). STEMCELL Technologies cannot guarantee biological function or any other properties associated with performance of cells in researchers' individual assay systems.

Donor Criteria And Screening

Primary cell products are obtained from volunteers participating in an Institutional Review Board (IRB) or Human Subject Committee approved donor program. All donors sign an informed consent form and are aware that these cells are used for research purposes. Donors must be between 18 and 55 years of age, and may not be pregnant (with the exception of CB products). Each volunteer is pre-screened for general health, HIV, hepatitis B and hepatitis C using FDA approved methods. When infectious serological tests for donors are not available, the cell products are tested for the presence of viral DNA by PCR.

WARNING: Treat cell products as potentially contaminated biological specimens even if available serological reports are negative. All human cell products should be handled at Biosafety Level 2 or higher. For more information, please contact your site Safety Officer.

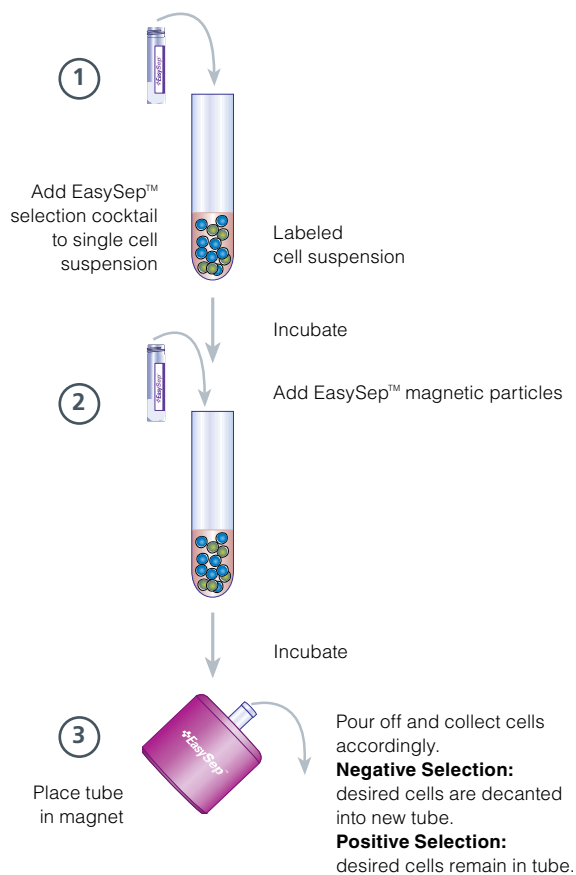
STEMCELL Technologies is the exclusive worldwide distributor for AllCells™, LLC (excluding China & Japan), with co-distribution rights in North America.

*Primary Cell products are for in vitro research use only. Not approved for diagnostic, therapeutic, or clinical applications. Not approved for human or veterinary use in vivo.

Cell Separation

Additional cell types may be positively or negatively selected from the MNC fraction of human BM, PB or CB using immunomagnetic cell separation. Cells are targeted for selection or depletion by crosslinking them to EasySep™ magnetic particles using STEMCELL Technologies' Tetrameric Antibody Complex (TAC) technology. A tube containing the the magnetically labeled cells is then placed in an EasySep™ magnet. After a short incubation, the unlabeled cells are poured off, leaving the magnetically labeled cells behind. EasySep™ magnetic particles do not interfere with flow cytometry and isolated cells are available for use immediately.

FIGURE 1. Typical EasySep™ Procedure for Human Cells



EasySep™ Equipment

EasySep™ Magnet • Catalog #18000



The EasySep™ magnet is designed for cell separation procedures using EasySep™ reagents. The EasySep™ magnet generates a high-gradient magnetic field in the interior cavity that is strong enough to separate cells labeled with EasySep™ Magnetic Particles. This magnet is designed to hold a standard FACS tube (12 x 75 mm polystyrene tube).

“The Big Easy” EasySep™ Magnet • Catalog #18001



“The Big Easy” EasySep™ Magnet is designed for cell separation procedures using EasySep™ reagents when processing a large number of cells, or when using RoboSep™. Like the EasySep™ Magnet, “The Big Easy” EasySep™ Magnet generates a high-gradient magnetic field in the interior cavity that is strong enough to separate cells labeled with EasySep™ Magnetic Particles. The “Big Easy” EasySep™ Magnet is designed to hold a standard 17 x 100 mm (14 mL) polystyrene tube.

“Easy 50” EasySep™ Magnet • Catalog #18002



“Easy 50” EasySep™ Magnet is the largest of the EasySep™ magnets. Like the EasySep™ Magnet, the “Easy 50” EasySep™ Magnet generates a high-gradient magnetic field in the interior cavity that is strong enough to separate cells labeled with EasySep™ Magnetic Particles. “Easy 50” is designed to hold a standard 50 mL conical tube.

Note: This magnet is for use with certain negative selection kits only. For a complete list of compatible kits, please refer to the “Easy 50” EasySep™ Magnet Product Information Sheet available on our website.

RoboSep™ • Catalog #20000



RoboSep™ is the first instrument to offer true walk-away automation of magnetic cell separation. Using column-free EasySep™ technology to isolate cells by either positive or negative selection, RoboSep™ can label and separate up to four samples at one time.

Advantages

FAST. Isolate untouched cells in as little as 25 minutes.

HIGH CELL PURITIES AND RECOVERIES. Achieve up to 99% cell purity.

EXTREMELY VERSATILE. Select or deplete virtually any cell type from any species in a standard polystyrene tube.

FLOW CYTOMETRY COMPATIBLE. Use isolated cells immediately for culture, flow cytometry and other downstream applications.

Contract Assay Services

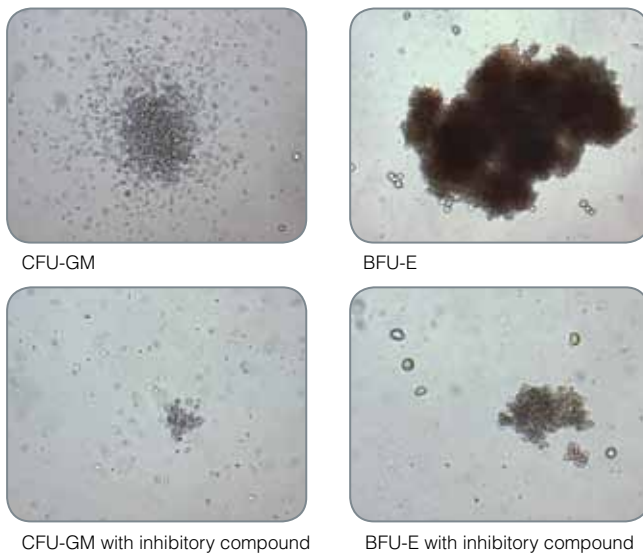
Contract Assay Services combines the power of quality STEMCELL Technologies reagents with the practical knowledge of more than 50 highly qualified scientists with expertise in hematopoietic, neural and mesenchymal stem cell biology to provide customized assay services to meet your needs. We have performed more than 500 studies for more than 100 organizations around the world.

Contract Assay Services offers validated and customized cell-based assays to:

- Determine safety and efficacy of compounds
- Provide clinically relevant, cost-effective information through the drug discovery process
- Reduce animal testing



FIGURE 2. Colony size changes in the presence of an inhibitory compound



Hematopoietic Stem and Progenitor Cell Assays

- Screen for toxic effects of compounds in blood or bone marrow
- Determine the hematopoietic modulating activity of small molecule compounds
- Assess hematological parameters and progenitor content in clinical samples

Mesenchymal Stem and Progenitor Cell Assays

- Determine progenitor frequency and proliferative potential
- Examine overall expansion potential
- Identify differentiation potential of mesenchymal stem cells into adipose, osteogenic and chondrocyte lineages

Neural Stem and Progenitor Cell Assays

- Determine proliferation potential using the Neural Colony-Forming Cell (NCFC) assay, which enables identification and quantification of both neural stem and progenitor cells
- Assess differentiation potential of neural stem and progenitor cells into neurons, astrocytes and oligodendrocytes

Cell Culture Media and Reagents

Hematopoietic Cell Expansion Media

STEMCELL Technologies provides two serum-free media formulations to support the maintenance and expansion of hematopoietic stem and progenitor cells.

Serum-Free Medium for Expansion of Hematopoietic Cells

PRODUCT: StemSpan™ SFEM
CATALOG #: 09600 100 mL/bottle
 09650 500 mL/bottle



RECOMMENDED FOR:

- The culture and expansion of human hematopoietic progenitors in samples of bone marrow, cord blood and mobilized peripheral blood.
- The expansion of mouse colony-forming cells (CFCs), long-term culture-initiating cells (LTC-IC) and long-term repopulating cells.
- The generation and culture of human and mouse dendritic cells.
- The culture of rat and non-human primate hematopoietic cells.

CONTAINS:

- Bovine Serum Albumin
- rh Insulin
- Human Transferrin (Iron-Saturated)
- 2-Mercaptoethanol
- Supplements
- Iscove's MDM

Xeno-Free Defined Medium for Expansion of Hematopoietic Cells

PRODUCT: StemSpan™ H3000
CATALOG #: 09800 100 mL/bottle
 09850 500 mL/bottle



RECOMMENDED FOR:

- The culture and expansion of human CD34⁺ cells and hematopoietic progenitors from bone marrow, mobilized peripheral blood and cord blood.

CONTAINS:

Contains only human-derived or recombinant human proteins.

Note: Recombinant cytokines for the optimal growth and expansion of hematopoietic cells must be added depending on researcher's requirement. STEMCELL Technologies offers a full range of recombinant cytokines. Visit www.stemcell.com for more information.

Hematopoietic Cell Colony Assay Media

The CFC assay allows for in vitro quantitative assessment of hematopoietic progenitors. In the CFC assay, progenitor cells, in response to cytokines and supplements in the culture medium, proliferate and differentiate into mature cell types that can be distinguished morphologically. The culture medium consists of a semi-solid matrix that allows for the clonal progeny of each progenitor cell to form distinct colonies that can be enumerated. The morphological nature of each colony represents the developmental potential of the original progenitor cell.

PRODUCT NAME	DESCRIPTION	CELL TYPE	APPLICATIONS
MethoCult™	Methylcellulose-based semi-solid medium for hematopoietic CFC assays	MNCs, CD34 ⁺ , CD133 ⁺	Quantitate and characterize human hematopoietic progenitors from BM, CB and PB. Screen effects of growth factors, inhibitors and drugs
CollagenCult™	Collagen-based semi-solid medium for hematopoietic CFC assays	MNCs, CD34 ⁺ , CD133 ⁺	Same applications as MethoCult™, however, samples can be dehydrated and fixed for immunocytochemical staining and long-term storage
MegaCult™-C	Collagen-based semi-solid medium for detection and quantitation of megakaryocytic progenitors	MNCs, CD34 ⁺ , CD133 ⁺	Detection and quantitation of megakaryocytic progenitors

MNC = mononuclear cells; BM = bone marrow; CB = cord blood; PB = peripheral blood

Mesenchymal Stem Cell (MSC) Media

CFU-F Assay and Expansion Kit

PRODUCT: MesenCult™ Proliferation Kit
(Human)

CATALOG #: 05411 500 mL



RECOMMENDED FOR:

- The in vitro culture and expansion of human mesenchymal stem cells, as well as their detection and enumeration using the CFU-F assay.

CONTAINS:

- MesenCult™ MSC Basal Medium (Human; Catalog #05401)
- Mesenchymal Stem Cell Stimulatory Supplement (Human; Catalog #05402)
- Kit components are also available separately.

Balanced Salt Solutions

PRODUCT	QUANTITY	CATALOG #
D-PBS	500 mL	37350
D-PBS, 10X Concentrate	500 mL	37354
D-PBS with 2% FBS	500 mL	07905
HBSS Modified (Ca ⁺⁺ & Mg ⁺⁺ free)	500 mL	37250
HBSS Modified (with HEPES, without Phenol Red)	500 mL	37150

Miscellaneous Culture Reagents

PRODUCT	QUANTITY	CATALOG #
L-glutamine, 200 mM	100 mL	07100
Sodium Pyruvate, 100 mM	100 mL	07000
DNase I, 1 mg/mL	1 mL	07900
Trypsin-EDTA (0.25%)	500 mL	07901
Trypsin-EDTA (0.05%)	500 mL	07910
Penicillin and Streptomycin Solution (100X)	100 mL	07500
3% Acetic Acid with Methylene Blue	100 mL	07060
Trypan Blue	100 mL	07050
Ficoll-Paque™ PLUS*	100 mL	07907
	500 mL	07957

*Ficoll™ and Ficoll-Paque™ PLUS are trademarks of GE Healthcare Ltd.

Copyright © 2011 by STEMCELL Technologies Inc. All rights reserved including graphics and images. STEMCELL Technologies and Design, STEMCELL shield, THE CELL EXPERTS and EasySep, RoboSep, StemSpan, MethoCult, CollagenCult, MegaCult-C, and MesenCult are trademarks of STEMCELL Technologies Inc. Ficoll is a trademark of GE Healthcare Ltd. All other trademarks are the property of their respective holders.

Xeno-Free Culture Kit

PRODUCT: MesenCult™-XF Culture Kit
(Human)

CATALOG #: 05429 1 Kit



RECOMMENDED FOR:

- MSC expansion from primary human bone marrow or culture-expanded cells.
- Long-term culture of MSCs (>8 passages).
- Detection and enumeration of CFU-F from primary human bone marrow and culture-expanded MSCs.

USAGE:

- The MesenCult™-ACF Dissociation Kit (Catalog #05426) must be used when passaging cells. Components are prescreened to ensure optimal cell attachment when MSCs are cultured with MesenCult™-XF Medium.
- MesenCult™-XF Medium must be supplemented with L-glutamine (Catalog #07100).

CONTAINS:

- MesenCult™-XF Medium (includes Basal Medium and Proliferation Supplement; Catalog #05420)
- MesenCult™-XF Attachment Substrate (Catalog #05424)
- Kit components are also available separately.

General Cell Culture Media

PRODUCT	QUANTITY	CATALOG #
DMEM with 4500 mg D-glucose/L	500 mL	36250
DMEM with 1000 mg D-glucose/L	500 mL	36253
DMEM/F-12	500 mL	36254
Iscove's MDM (IMDM)	500 mL	36150
IMDM with 2% FBS	100 mL	07700
McCoy's 5A Medium (Modified)	500 mL	36350
MEM Alpha Modification with Nucleosides	500 mL	36450
MEM Alpha Modification without Nucleosides	500 mL	36453
MEM	500 mL	36550
RPMI 1640	500 mL	36750

For more information on cell culture media and reagents, visit www.stemcell.com or contact us at info@stemcell.com.

THE CELL EXPERTS™ | WWW.STEMCELL.COM

TOLL-FREE PHONE 1 800 667 0322 • PHONE 1 604 877 0713

TOLL-FREE FAX 1 800 567 2899 • FAX 1 604 877 0704

ORDERS@STEMCELL.COM • INFO@STEMCELL.COM

FOR FULL CONTACT DETAILS WORLDWIDE VISIT OUR WEBSITE

FOR RESEARCH USE ONLY. NOT FOR THERAPEUTIC OR DIAGNOSTIC USE. CATALOG #28929 VERSION 3.0.0 APRIL 2011

