

The World Leader in Tools for Hematopoietic Stem Cell Research

A Complete Set of Tools for Your Hematopoietic Colony-Forming Cell (CFC) Assays

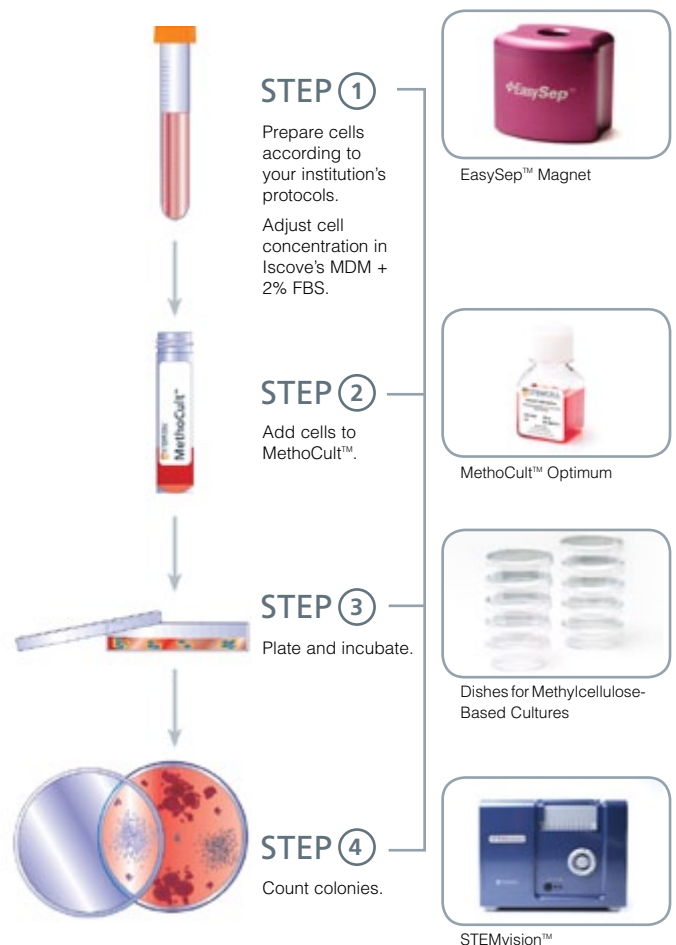
STEMCELL Technologies Inc. offers a complete set of tools for each step of the hematopoietic CFC assay. MethoCult™ methylcellulose-based medium is a key component of standardized hematopoietic CFC assays. Other tools, including proficiency testing programs, quality control kits and training courses, are available to help researchers and clinicians increase and maintain knowledge and proficiency in hematopoietic assays.

Products Associated with Procedure For Hematopoietic CFC Assays

STEP 1	
Ammonium Chloride Solution*	07800/07850
Ficoll-Paque™ PLUS*	07907/07917 07957/07967
EasySep™*	Immunomagnetic positive selection or enrichment
RosetteSep™*	Immunodensity enrichment
Iscove's MDM + 2% FBS	07700
MethoCult™ Cell Wash Medium**	87700
STEP 2	
MethoCult™ Express	04437/04447
MethoCult™ Optimum	04034/04044
MethoCult™ Optimum without EPO	04035/04045
MethoCult™ Enriched	04435/04445
MethoCult™ Enriched without EPO	04535/04545
MethoCult™ GF H84434**	84434/84444
MethoCult™ GF H84534**	84534/84544
MethoCult™ GF H84435**	84435/84445
MethoCult™ GF H84535**	84535/84545
STEP 3	
Syringes	28230/28240
Blunt-End Needles	28110/28120
Dishes for Methylcellulose-Based Cultures	27100/27150
Outer Dishes	27140/27141
STEP 4	
STEMvision™	22001/22002
Gridded Scoring Dishes	27500
Cord Blood Colony Atlas	29940
Mouse Hematopoietic Progenitor Wallchart Request a free copy at www.stemcell.com/methocultwallchart	

* Optional. Use of these reagents is dependent on institution's protocol for processing samples.

** For in vitro diagnostic use in the European Union (EU).



THE CELL EXPERTS™ | WWW.STEMCELL.COM

CATALOG #28790 | VERSION 1.0.0 | APRIL 2011

TOLL-FREE T. 1 800 667 0322 • T. +1 604 877 0713 • TOLL-FREE F. 1 800 567 2899 • F. +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.

The World Leader in Tools for Hematopoietic Stem Cell Research



STEP ① Prepare cells according to your institution's protocols.

- Removal of red blood cells using ammonium chloride lysis or density gradient separation is recommended.
- Progenitors can be enriched with EasySep™ or RosetteSep™ (optional).

Adjust cell concentration in Iscove's MDM + 2% FBS.

- Prepare a cell concentration that is 10X the final plating concentration (e.g. 2×10^6 cells/mL for 2×10^5 cells/dish). Note: A 10X cell concentration should only be used if adding 300 μ L of cells to 3.0 mL of MethoCult™ for duplicate dishes or 400 μ L of cells to 4.0 mL of MethoCult™ for triplicate dishes.

STEP ② Add cells to MethoCult™.

- Vortex tube to mix contents thoroughly.
- Let stand for at least five minutes to allow bubbles to rise to the top.

STEP ③ Plate and incubate.

- Dispense MethoCult™ and cell mixture into dishes for methylcellulose-based cultures.
- Place the culture dishes in an outer dish along with uncovered 35 mm dishes containing sterile water. This helps maintain humidity.
- Incubate for 14 - 16 days (7 days for MethoCult™ Express) at 37°C, in 5% CO₂, with \geq 95% humidity.

STEP ④ Count colonies.

- Use STEMvision™ or an inverted microscope and gridded scoring dish to enumerate colonies.

For detailed instructions, refer to the Technical Manual for Human Colony-Forming Cell (CFC) Assays Using MethoCult™ (Manual Catalog #28404).

Resources for the Hematopoietic CFC Assay

PRODUCT	CATALOG #
Human Bone Marrow Proficiency Testing Program	00602
Frozen Cord Blood Proficiency Testing Program	00608
Human Bone Marrow Quality Control Kit	00650
Human Cord Blood Quality Control Kit	00651
Hematopoietic Progenitor Assay Training Course	00215

Visit www.stemcell.com for more information.

VIDEO

Procedure for Setting Up the CFC Assay

www.stemcell.com/cfcmethodvideo

SCAN ME ▶

