# Dynabeads® CD34 Positive Isolation Kit

#### Catalog no. 11301D

#### Store at 2 to 8°C

Rev. Date: November 2011 (Rev. 004)

## **Kit Contents**

Kit contents	Volume		
Dynabeads® CD34	5 mL		
DETACHaBEAD® CD34	5 mL		
Kit capacity			

MNC: ~5 × 10<sup>9</sup>

Dynabeads<sup>®</sup> CD34 contains  $4 \times 10^8$  beads/mL in phosphate buffered saline (PBS), with 0.1% bovine serum albumin (BSA) and 0.02% sodium azide as a preservative. DETACHaBEAD<sup>®</sup> CD34 contains a polyclonal anti-Fab antibody in 0.15 M PBS.

## **Product Description**

This product is intended for positive magnetic isolation of CD34<sup>+</sup> cells from mononuclear cells (MNCs) from human bone marrow, peripheral blood or cord blood. In the first step CD34<sup>+</sup> cells are captured by the Dynabeads<sup>®</sup> and isolated using a magnet. In the second step beads are removed from the cells.

Isolated cells are bead- and antibody-free, phenotypically unaltered, and suitable for any downstream application including flow cytometry, functional studies and cell culture.

## Required Materials

- Magnet (DynaMag<sup>™</sup>): See www. lifetechnologies.com/magnets for recommendations.
- Mixer allowing tilting and rotation of tubes (e.g. HulaMixer<sup>®</sup> Samp4 x le Mixer).
- Isolation Buffer: PBS

   (Ca<sup>2+</sup> and Mg<sup>2+</sup> free with
   0.1% BSA 2 mM EDTA, pH 7.4.

   Note: BSA can be replaced by

   human serum albumin (HSA)
   or fetal calf serum (FCS). EDTA
   can be replaced by 0.6% sodium
   citrate. PBS containing
   Ca<sup>2+</sup> or Mg<sup>2+</sup> is not

   recommended, except when
   required for DNase treatment.
- Culture Medium: e.g. RPMI 1640 with 2% FCS.
- DNase I (when required).

## **General Guidelines**

- Visit www.lifetechnologies. com/samplepreparation for recommended sample preparation procedures.
- Use a mixer that provides tilting and rotation of the tubes to ensure that Dynabeads<sup>®</sup> do not settle in the tube.
- This product should not be used with MPC<sup>™</sup>-1 (Cat. no. 12001D).
- Avoid air bubbles (foaming) during pipetting.
- Carefully follow the recommended pipetting volumes and incubation times.
- Keep all buffers cold.

## Protocol

#### Wash Dynabeads®

See Table 1 for volume recommendations.

- 1. Resuspend the Dynabeads<sup>®</sup> in the vial (vortex >30 sec or tilt and rotate for 5 min).
- 2. Transfer the desired volume of Dynabeads<sup>®</sup> to a tube.
- 3. Add the same volume of Isolation Buffer, or at least 1 mL, and mix.
- 4. Place the tube in a magnet for 1 min and discard the supernatant.
- 5. Remove the tube from the magnet and resuspend the washed Dynabeads<sup>®</sup> in the same volume of Isolation Buffer as the initial volume of Dynabeads<sup>®</sup> (step 2).

#### Isolate CD34<sup>+</sup> Stem Cells

The isolation and release protocol is based on 1 mL MNC  $(4 \times 10^7 - 1 \times 10^8 \text{ cells/mL})$  as starting sample, but can be scaled up according to Table 1.

- 1. Resuspend the prepared cell sample thoroughly with a narrow pipette.
- 2. Add 100 μL washed and resuspended Dynabeads<sup>®</sup> to the prepared sample and vortex 2–3 sec.
- 3. Incubate at 2°C to 8°C for 30 min with gentle tilting and rotation.
- 4. Fill the tube with cold buffer to the height of the magnet (or at least 1 mL) and resuspend the cell-bead complexes.
- 5. Place the tube in a magnet for 2 min and discard the supernatant.
- 6. Resuspend the bead-bound cells in 2 mL buffer by vortexing or pipetting. Separate on a magnet for 1 min.
- 7. Repeat step 6 twice to increase the purity.
- 8. Resuspend the bead-bound cells in  $100 \,\mu\text{L}$  buffer.

#### Prepare MNC

- Prepare MNC according to "General Guidelines".
- Prepare MNC to  $4 \times 10^7 1 \times 10^8$  MNC/mL. Allways use fresh samples.

#### Release CD34<sup>+</sup> Stem Cells

Tubes and buffers can be kept at room temperature for the release procedure.

- Add 100 μL DETACHaBEAD<sup>®</sup> (never use less than 100 μL DETACHaBEAD<sup>®</sup>).
- 10. Incubate at room temperature for 45 min with tilting and rotation. Keep the sample in the bottom of the tube.
- 11. Add 2 mL Isolation Buffer and vortex 2–3 sec to enhance detachment of beads from the cells.
- 12. Place the tube in a magnet for 2 min.
- 13. Transfer the supernatant containing released cells to a fresh tube. To obtain residual cells, wash the beads 3 times in 500  $\mu$ L Isolation Buffer and pool the supernatants.
- 14. Wash the cells thoroughly by resuspending in 10 mL Isolation Buffer. Centrifuge for 10 min at  $400 \times g$  to remove excess DETACHaBEAD<sup>®</sup>.
- 15. Resuspend the cells in Isolation Buffer or Culture Medium.

The isolated cells are pure, viable, and free from antibody bound to the surface, and may be used in any downstream application.

Table 1: Volumes for human CD34<sup>+</sup> stem cell isolation from MNC. Use  $4 \times 10^7 - 1 \times 10^8$  cells/mL. For lower volumes than 1 mL, use the same volumes as the 1 mL column. For higher cell numbers, scale up the volumes accordingly.

Step	Step description	Cell volume: 1 mL	Cell volume: 5 mL
	Recommended tube size	5 mL tube	15 mL tube
	Recommended magnet	DynaMag <sup>™</sup> -5	DynaMag <sup>™</sup> -15
2	Dynabeads® CD34	100 µL	500 μL
4	Resuspend cells (Isolation Buffer)	~3 mL	~8 mL
6-7	Wash cells (Isolation Buffer)	2 mL × 2	10 mL × 2
8	Resuspend cells (Isolation Buffer)	100 μL	500 μL
9	Release cells (DETACHaBEAD®)	100 μL	500 μL
11	Increase cell volume (Isolation Buffer)	2 mL	10 mL
13	Collect residual cells	500 μL × 3	2.5 mL × 3
14*	Remove DETACHaBEAD®	Increase to 10 mL	Increase to 30 mL

\*Transfer the sample to a larger tube that is more appropriate to the volume (e.g. 15 mL tube and 50 mL tube, respectively).

## Description of Materials

Dynabeads<sup>®</sup> CD34 are uniform, superparamagnetic polymer beads (4.5 µm diameter) coated with a primary monoclonal antibody specific for the CD34 membrane antigen predominantly expressed on human hematopoietic progenitor cells and endothelialprogenitor cells. DETACHaBEAD<sup>®</sup> CD34 is a polyclonal anti-Fab antibody specific for the CD34 antibody on the Dynabeads<sup>®</sup> CD34.

## **Related Products**

Product	Cat. no.
DynaMag <sup>™</sup> -5	12303D
DynaMag™-15	12301D
DynaMag™-50	12302D
HulaMixer <sup>®</sup> Sample Mixer	15920D

**REF** on labels is the symbol for catalog number.

## Limited Use Label License

The purchase of this product conveys to the purchaser the limited, nontransferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

Manufactured by Life Technologies AS, Norway. Life Technologies AS complies with the Quality System Standards ISO 9001:2008 and ISO 13485:2003.

SPEC-06167

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners, except where otherwise stated. Life Technologies and/or its affiliate[s] disclaim all warranties with respect to this document, expressed or implied, including but not limited to those of merchantability or fitness for a particular purpose. In no event shall Life Technologies and/or its affiliate[s] disclaim all warranty, or under any statute or on any other basis for special, incidental, indirect, punitive, multiple or consequential damages in connection with or arising from this document, including but not limited to the use thereof.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com



www.lifetechnologies.com