

# Dynabeads® FlowComp™ Mouse CD8

Catalog no. 11462D

#### Store at 2°C to 8°C

Rev. Date: December 2011 (Rev. 004)

#### Kit Contents

Kit contents	Volume
FlowComp™ Mouse CD8 antibody	1 mL
FlowComp <sup>™</sup> Dynabeads <sup>®</sup>	3 mL
FlowComp™ Release Buffer	2 × 20 mL

#### Kit capacity

 $\sim 2 \times 10^9 \text{ cells}$ 

FlowComp<sup>™</sup> Dynabeads® contains ~1 × 109 (~10 mg) beads/mL in phosphate buffered saline (PBS), pH 7.4, with 0.1% bovine serum albumin (BSA) and 0.02% sodium azide as a preservative. FlowComp<sup>™</sup> Mouse CD8 antibody contains monoclonal anti-mouse CD8 antibody in PBS with 0.5% BSA and 0.02% sodium azide. FlowComp<sup>™</sup> Release Buffer contains modified biotin in 0.1% BSA and 2 mM EDTA. **Caution:** Sodium azide may react with lead and copper plumbing to form highly explosive metal azides.

# **Product Description**

This product is intended for positive magnetic isolation of CD8+T cells from lymphoid organs. The isolated cells are highly pure, viable, and bead-free (fig. 1). In the first step, FlowComp™ Mouse CD8 antibody is added and will bind to the target cells. In the second step, CD8+T cells that have bound the specific antibodies are captured by the Dynabeads®. In the third and last step, the cells are released from the Dynabeads®.

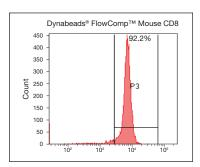


Figure 1: Mouse CD8<sup>+</sup> cells isolated with Dynabeads<sup>®</sup> FlowComp<sup>™</sup> Mouse CD8

## **Downstream Applications**

Isolated cells are bead-free and may be used directly in any downstream application including flow cytometry. The cells readily proliferate in response to Dynabeads® Mouse T Activator CD3/CD28 and can be measured by incorporation of EdU or in a CFSE assay.

# Required Materials

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

  Ca<sup>2+</sup> and Mg<sup>2+</sup> free PBS
  supplemented with
  0.1% BSA and 2 mM EDTA.

  Note: BSA can be replaced by
  human serum albumin (HSA) or
  2% fetal bovine serum (FBS)/fetal
  calf serum (FCS).

- Optional: We recommend using anti-mouse CD8 $\alpha$ -PE or anti-mouse CD3 Alexa Fluor® 488 as primary fluorescent antibody for flow staining of cells after isolation. Avoid using secondary antibodies specific for rat antibodies for flow cytometry staining.
- Optional: Red blood cell lysis buffer.
   See www.lifetechnologies.com/samplepreparation.
- Optional: For viability analysis, SYTOX® Red is recommended.

### 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® do not settle at the bottom of the tube.
- This product should not be used with magnet MPC<sup>™</sup>-1.
- Follow the recommended volumes and incubation times (Table 1).
- Avoid air bubbles (foaming) during pipetting.
- To avoid unspecific labeling of cells during flow staining we recommend using gamma-globulin or Fc blocking reagents prior to staining with primary fluorescent antibody.
- Only DSB-X biotinylated antibodies can be used in the isolation process.
   Standard biotinylated antibodies will not give release of cells after isolation.

### **Protocol**

Approximately 30–35% of mouse spleen cells are T cells, and about 30% of these T cells strongly express the CD8 antigen. This kit isolates highly pure CD8⁺ T cells using Dynabeads® FlowComp™ Mouse CD8.

### Prepare Cells

- Prepare a single cell suspension from lymphoid organs (e.g. lymph nodes or spleen) according to "General Guidelines".
- Resuspend the cells at  $1 \times 10^8$  cells/mL in Isolation Buffer.
- Prepare approximately 10 mL of Isolation Buffer per  $5 \times 10^7$  cells.

## Wash Dynabeads®

See Table 1 for volume recommendations.

- 1. Resuspend the Dynabeads® in the vial (i.e. vortex for >30 sec, or tilt and rotate for 5 min).
- 2. Transfer the desired volume of Dynabeads® to a tube.
- 3. Add the same volume of Isolation Buffer, or at least 1 mL, and resuspend.
- 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® in the same volume of Isolation Buffer as the initial volume of Dynabeads® (step 2).

#### **Isolate Cells**

This protocol is based on  $5 \times 10^7$  cells, but is directly scalable from  $1 \times 10^7$  to  $5 \times 10^8$  cells, according to Table 1. When working with fewer cells than  $1 \times 10^7$ , use the same volumes as for  $1 \times 10^7$ .

- 1. Transfer 500  $\mu$ L (5 × 10<sup>7</sup>) prepared cells to a tube and add 25  $\mu$ L FlowComp<sup>TM</sup> Mouse CD8 antibody.
- 2. Mix well and incubate 10 min at 2°C to 8°C.
- 3. Wash by adding 2 mL Isolation Buffer and centrifuge 8 min at  $350 \times g$ .
- 4. Remove the supernatant, and resuspend in 1 mL Isolation Buffer.
- Add 75 µL washed FlowComp™ Dynabeads® and mix well (e.g. vortex 2–3 seconds).
- 6. Incubate for 15 min at 2°C to 8°C under rolling and tilting.
- 7. Add 1 mL Isolation Buffer, pipet 2–3 times (or vortex 2–3 seconds) and place the tube in a magnet for 2 min.
- 8. While the tube is still in the magnet, carefully remove and discard the supernatant containing the CD8 negative cells.
- 9. Repeat steps 7–8 at least once to wash the bead-bound CD8+ cells. These steps are critical to obtain a high purity of isolated cells.

#### Release Cells

- 10. Resuspend the bead-bound cells in 1 mL Release Buffer.
- 11. Incubate 10 min with rolling and tilting at room temperature.
- 12. Pipet 10 times to efficiently release the cells and place in a magnet for 2 min. Avoid foaming.
- 13. Transfer the supernatant containing the bead-free CD8+ cells to a new tube and again place on the magnet for 1 min to remove any residual beads. Transfer again the supernatant containing the bead-free cells to a new tube.
- 14. Add 2 mL Isolation Buffer followed by centrifugation for 8 min at  $350 \times g$ . Discard the supernatant and resuspend the cell pellet in preferred cell medium.

Keep the cells on 2°C to 8°C until further use in downstream applications.

Table 1: Volumes for mouse CD8+T cells. This protocol is scalable from 1  $\times$   $10^7$  to 5  $\times$   $10^8$  cells.

Step	Step description	Volumes per 5 × 10 <sup>7</sup> cells	Volumes per 5 × 10 <sup>8</sup> cells
	Recommended tube size	5 mL	15 mL
	Recommended magnet	DynaMag <sup>™</sup> -5	DynaMag <sup>™</sup> -15
1	Cell volume	500 μL	5 mL
1	FlowComp™ Human CD8 antibody	25 μL	250 µL
3*	Wash cells (Isolation Buffer)	~2 mL	~10 mL
4	Resuspend cells (Isolation Buffer)	1 mL	10 mL
5**	FlowComp™ Dynabeads®	75 μL	750 μL
7+9	Wash beads (Isolation Buffer)	2 × 1 mL	2 × 10 mL
10	FlowComp™ Release Buffer	1 mL	10 mL
14*	Wash cells (Isolation Buffer)	~2 mL	~20 mL

<sup>\*</sup> Adjust the Isolation Buffer volumes to fit to the tube you are using. For very large volumes use a larger tube than recommended in step 14 to successfully remove the biotin in the sample.

## **Description of Materials**

Dynabeads® FlowComp™ are uniform, superparamagnetic polystyrene beads (2.8  $\mu$ m in diameter) coated with modified streptavidin. FlowComp™ Mouse CD8 antibody contains a DSB-X conjugated monoclonal rat antimouse IgG2a CD8 $\alpha$  that recognizes Ly2.1 and Ly2.2 cells. FlowComp™ Release Buffer contains a modified biotin that out-competes the modified biotin on the antibody to give the cell release from the beads.

### Related Products

Product	Cat. no.
DynaMag™-5	12303D
DynaMag <sup>™</sup> -15	12301D
DynaMag <sup>™</sup> -50	12302D
HulaMixer® Sample Mixer	15920D
Rat anti-mouse CD8 R-PE	MCD0804
Hamster anti-mouse CD3 Alexa Fluor® 488	HM3420
Dynabeads® Mouse T-Activator CD3/CD28	11452D
Phosphate buffered saline	14190
Click-iT®-Edu	A10202
CFSE assay	C34554
SYTOX® Red	S34859

**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.

# **Limited Product Warranty**

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

SPEC-05960

©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 affiliatels] 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) be liable, whether in contract, tort, 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



<sup>\*\*</sup> When incubating, tilt and rotate the vial so the cells and beads are kept in the bottom of the tube. Do not perform end-over-end mixing if the volume is small relative to the tube size.