# Dynabeads<sup>®</sup> Untouched<sup>™</sup> Mouse CD4 Cells

#### Catalog nos. 11415D, 11416D

Store at 2°C to 8°C Rev. Date: March 2012 (Rev. 006)

Wash the cells and add Mouse

incubation.

Depletion Dynabeads® to bind to the

antibody labeled cells during a short

Apply to magnet and transfer the

mouse CD4<sup>+</sup> T cells to a new tube

and discard the bead-bound cells.

supernatant with the untouched

### **Kit Contents**

Kit contents	11415D Volume	11416D Volume*
Mouse Depletion Dynabeads®	2×10 mL	4 mL
Antibody Mix (for Mouse CD4 cells)	2 mL	0.4 mL

\* Sample size

#### Kit capacity

for 11415D:  $\sim 1 \times 10^{9}$  cells for 11416D\*:  $\sim 2 \times 10^{8}$  cells

Mouse Depletion Dynabeads<sup>®</sup> contains  $4 \times 10^8$  beads/mL in phosphate buffered saline (PBS), pH 7.4, containing 0.1% bovine serum albumin (BSA) and 0.02% sodium azide as a preservative. The Antibody Mix for mouse CD4 cells contains a mixture of rat monoclonal antibodies in PBS with 0.02% sodium azide. **Caution:** Sodium azide may react with lead and copper plumbing to form highly explosive metal azides.

# **Product Description**

This product is intended for isolation of untouched mouse CD4<sup>+</sup> T cells from mouse spleen or lymph node cells by depleting CD8<sup>+</sup> T cells, B cells, monocytes/ macrophages, NK cells, dendritic cells, erythrocytes and granulocytes. Isolated untouched mouse CD4<sup>+</sup> T cells are bead- and antibodyfree, and suitable for use in any downstream application (fig. 1).

Add a mixture of monoclonal rat IgG antibodies against the non-CD4<sup>+</sup> T cells to the starting sample to label the unwanted cells. Add Blocking Reagent and Antibody Mix to PBMC. Mix and incubate for 20 minutes at 2°C to 8°C. Wash cells and add Dynabeads\*. Incubate for 15 minutes at 18°C to 25°C. Increase volume and resuspend thoroughly. Apply magnet for 2 minutes.

# Discard beads with unwanted cells.

Figure 1: Simple method for isolating untouched mouse CD4\* T cells.

#### **Downstream Applications**

Isolated mouse CD4<sup>+</sup> T cells can be used in any application, (e.g., studies on CD4<sup>+</sup> T cell proliferation, apoptosis and induction of anergy, studies on antigen-specific T cells, studies on regulation of CD4<sup>+</sup> T cell cytokine expression, flow cytometry/FACS sorting). Isolated cells can be activated/ expanded using Dynabeads<sup>®</sup> Mouse T-Activator CD3/CD28 (polyclonal activation) or Dynabeads<sup>®</sup> Mouse T-Activator CD3/CD28/CD137 (antigenspecific activation).

#### **Required Materials**

- Magnet (DynaMag<sup>™</sup> portfolio). See www.lifetechnologies.com/magnets for recommendations.
- Mixing device with tilting and rotation, e.g. HulaMixer<sup>®</sup> Sample Mixer.
- Heat inactivated Fetal Bovine Serum (FBS)/Fetal Calf Serum (FCS).
- Isolation Buffer: PBS (Ca<sup>2+</sup> and Mg<sup>2+</sup> free) supplemented with 0.1% BSA and 2 mM EDTA.

**Note:** BSA can be replaced by human serum albumin (HSA) or 2% FBS/FCS. EDTA can be replaced by 0.6% sodium citrate.

#### **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 beads do not settle in the tube.
- This product should not be used with the MPC<sup>™</sup>-1 magnet (Cat. no. 12001D).
- Follow the recommended volumes and incubation times.
- Avoid air bubbles (foaming) during pipetting.
- Keep the buffers cold.

### Protocol

#### Wash the Beads

See Table 1 for volume recommendations.

- 1. Resuspend the beads in the vial (i.e. vortex for >30 sec, or tilt and rotate for 5 min).
- 2. Transfer the desired volume of beads 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 beads in the same volume of Isolation Buffer as the initial volume transferred of beads (step 2).

#### Prepare Sample

- Prepare spleen or lymph node cells according to "General Guidelines"
- Resuspend the cells at 1 × 10<sup>8</sup> cells/mL in Isolation Buffer. The protocol might need to be optimized if the cells are isolated from other sources.

#### Isolate Untouched Mouse CD4<sup>+</sup> Cells

This protocol is based on  $5 \times 10^7$  leucocytes, but it is scalable from  $1 \times 10^7 - 1 \times 10^9$  cells, see Table 1.

- 1. Transfer 500  $\mu$ L (5 × 10<sup>7</sup>) leucocytes in Isolation Buffer to a tube.
- 2. Add 100  $\mu$ L heat inactivated FCS/FBS.
- 3. Add 100 µL of Antibody Mix.
- 4. Mix well and incubate for 20 min at 2°C to 8°C.
- Wash the cells by adding 10 mL Isolation Buffer. Mix well by tilting the tube several times and centrifuge at 350 × g for 8 min at 2°C to 8°C. Discard the supernatant.
- 6. Resuspend the cells in 4 mL Isolation Buffer.
- 7. Add 1 mL pre-washed and resupended Mouse Depletion Dynabeads®.
- 8. Incubate for 15 min at 18°C to 25°C with gentle tilting and rotation.
- 9. Add 5 mL Isolation Buffer.
- 10. Resuspend the bead-bound cells by gently pipetting 5 times using a pipette with a narrow tip opening. Avoid foaming.
- 11. Place the tube in the magnet for 2 min and transfer the supernatant containing the untouched CD4<sup>+</sup> T cells to a new tube.

Table 1: Volumes for isolation of mouse CD4\* T cells. This protocol is scalable from 1  $\times$  107 to 3  $\times$  108 leucocytes.

Step	Step description	Volumes per 5 × 10 <sup>7</sup> leucocytes	Volumes per 3 × 10 <sup>8</sup> leucocytes
	Recommended tube	15 mL tubes	50 mL tubes
	Recommended magnet	DynaMag <sup>™</sup> -15	DynaMag <sup>™</sup> -50
1	Cell volume	500 µL	3 mL
2	FCS/FBS	100 µL	600 µL
3	Antibody Mix for mouse CD4 cells	100 µL	600 µL
5*	Wash cells (Isolation Buffer)	~10 mL	~30 mL
6	Resuspend cells (Isolation Buffer)	4 mL	~24 mL
7**	Mouse Depletion Dynabeads®	1 mL	6 mL
9*	Volume added before magnet	~5 mL	~15 mL

\* Adjust the Isolation Buffer volumes to fit to the tube you are using

\*\* When incubating, tilt and rotate 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.

# **Description of Materials**

Mouse Depletion Dynabeads<sup>®</sup> are uniform, superparamagnetic polystyrene beads (4.5 µm diameter) coated with a polyclonal sheep anti-rat IgG antibody. The Antibody Mix for mouse CD4 cells contains a mixture of rat monoclonal antibodies against mouse CD8, CD45R (B220), CD11b (Mac1), Ter-119, and CD16/CD32.

#### **Related Products**

Product	Cat. no.
DynaMag <sup>™</sup> -5	12303D
DynaMag™-15	12301D
DynaMag <sup>™</sup> -50	12302D
HulaMixer <sup>®</sup> Sample Mixer	15920D
Dynabeads <sup>®</sup> Mouse T-Activator CD3/CD28	11452D
Dynabeads® Mouse T-Activator CD3/CD28/CD137	11454D

**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-05651

©2012 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners. 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 EVENTS HALL 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



www.lifetechnologies.com