

# SMC™ Capture Labeling Kit Instructions

---

Capture Labeling Kit

Catalog #03-0077-02

Kit instructions for the labeling of capture antibody  
and conjugation to microparticles for use with  
**SMC™ Analyte Specific Immunoassays**

**FOR RESEARCH USE ONLY**

**NOT FOR USE IN DIAGNOSTIC PROCEDURES**

**Manufactured & Distributed by:**

---



EMD Millipore Corporation  
3050 Spruce Street  
St. Louis, Missouri 63103  
United States of America  
[emdmillipore.com](http://emdmillipore.com)

The M logo is a trademark of Merck KGaA,  
Darmstadt, Germany.  
© 2013 EMD Millipore Corporation,  
Billerica, MA 01821 USA.

**This page intentionally left blank.**

# TABLE OF CONTENTS

INTRODUCTION.....	2
REAGENTS .....	3
Reagents Provided.....	3
Storage Instructions .....	3
General Supplies Required But Not Provided .....	4
PRECAUTIONS .....	4
CAPTURE ANTIBODY LABELING .....	5
Preparation of Capture Antibody For Labeling .....	5
Label Capture Antibody .....	6
COATING OF MICROPARTICLE BEADS WITH CAPTURE ANTIBODY ...	7
Preparation and coating Microparticle Beads for Conjugation .....	7
WORKSHEET .....	9
ORDERING INFORMATION.....	9

## INTRODUCTION

The SMC™ Immunoassay uses a quantitative fluorescent sandwich immunoassay technique to measure analyte in matrix. This kit is intended for the labeling of **0.3-1.0 mg** of a capture antibody (Ab) and conjugation to magnetic microparticle (MP) beads. The labeled capture antibody is used in conjunction with a sandwich immunoassay to be developed for the SMC Immunoassay system.

This kit should be used in conjunction with 03-0076-02 SMC™ Detection Labeling kit and 03-0078-00 SMC™ Bead Based Immunoassay Development Kit.

**Antibodies used are to be in carrier- and label-free buffer.**

**Note: If labeling proteins other than antibodies of a 150 kD approximate molecular weight, contact MilliporeSigma Technical Support.**

**FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.**

## REAGENTS

The SMC™ Capture Labeling Kit includes all reagents listed in Reagents Provided. Additional reagents and supplies may be required to perform the labeling, as listed in the section titled General Supplies Required But Not Provided. All reagents supplied are for Research Use Only.

### Reagents Provided

Item#	Description	Shipping Conditions	Storage Conditions	Component Part No.	Packaging Details
1	Capture Label	On dry ice	≤ -70°C	02-0557-00	1 x 20 µL
2	Buffer 1	With cold pack	2 - 8°C	02-0552-00	1 x 25 mL
3	Buffer 2	With cold pack	2 - 8°C	02-0553-00	1 x 500 µL
4	Buffer 3	With cold pack	2 - 8°C	02-0554-00	1 x 5 mL
5	Filter Tube Ultrafree	With cold pack	2 - 8°C	02-0555-00	2 pack
6	Ultra 4 Filter Tube with Ultracel-30 membrane	With cold pack	2 - 8°C	02-0556-00	2 pack
7	10X Wash Buffer	With cold pack	2 - 8°C	02-0001-03	1 x 50 mL
8	Coated BeadBuffer	With cold pack	2 - 8°C	02-0558-00	1 x 5 mL
9	Uncoated SMC™ Beads	With cold pack	2 - 8°C	02-0559-00	1 x 4 mL

Note: Additional Ultra-4 Centrifugal Filter Units with Ultracel-30 membrane can be purchased from EMD Millipore (Catalogue # UFC803008).

### Storage Instructions

- The SMC™ Capture Labeling Kit should be stored at 2-8°C.
- The Capture Label should be stored at ≤ -70°C.
- Proper kit performance can only be guaranteed if the materials are stored properly.
- 10X Wash Buffer does not contain preservative. After dilution, the 1X wash buffer may be filter sterilized (Stericup® filter, EMD Millipore PN SCGPU11RE) for storage of up to 1 month at 2 - 8°C. If not filter sterilized, all remaining 1X Wash Buffer should be discarded upon experiment completion.

## General Supplies Required But Not Provided



- De-ionized or distilled water (DI Water)
- Pipettes capable of transferring 5 µL- 5 mL
- Polypropylene Eppendorf tubes
- Centrifuge with swinging buckets (capable of holding 15 mL conical) that can spin at 3900 x g (3900 RCF)
- Spectrophotometer or nanodrop (capable of reading A280)
- Bench top vortex
- Bench top mini-centrifuge
- 15 mL and 50 mL Conical tubes
- Sodium Azide (NaN<sub>3</sub>) (optional)

(Please contact your technical services representative for additional information or assistance selecting required but not provided supplies.)

## PRECAUTIONS

- Use caution when handling biological samples. Wear protective clothing and gloves.
- Components of this reagent kit contain approximately 0.1% sodium azide as a preservative. Sodium azide is a toxic and dangerous compound when combined with acids or metals. Solutions containing sodium azide should be disposed of properly.

## Full Hazardous Label:

Ingredient, Cat #		Full Label	
02-0001-03	10X Wash Buffer		<b>Warning.</b> Causes serious eye irritation. IF IN EYES: Rinse cautiously with water for several minutes. Remove contact lenses, if present and easy to do. Continue rinsing.
02-0557-00	Capture Label		<b>Warning.</b> Combustible liquid.

# CAPTURE ANTIBODY LABELING

## Preparation of Capture Antibody For Labeling

Antibodies need to be unlabeled and free of protein carrier such as BSA as well as supplemental reagents such as gelatin and glycerol to ensure success. Obtain carrier-free antibodies or purify these materials before proceeding.

Some antibody stabilizers (including sodium azide) and amine-containing buffers (including Tris) will significantly reduce capture labeling. Follow the steps below to remove these prior to antibody coupling.

1. Bring all reagents to room temperature.
2. Dilute antibody to approximately 1 mg/mL in Buffer 1.
3. Confirm the antibody concentration by reading the absorbance at 280 nm as follows:

- a. Prepare 100  $\mu$ L of a 1:10 dilution of the 1 mg/mL antibody solution in Buffer 1.
- b. Measure the absorbance (using Buffer 1 as a blank).
- c. Using Beer's Law  $A=\epsilon cl$ , calculate the antibody concentration as shown below (the example assumes a 1 cm path length and a molar extinction coefficient of 1.4):

$$\text{Conc. (mg/mL)} = \frac{A_{280}}{1.4} \times 10$$

4. Use the calculated antibody concentration to calculate the volume of antibody solution needed to label 0.2-1 mg Antibody.
5. Prepare the Ultra 4 30K filter tube:
  - a. Pre-wet the Ultra 4 30K filter tube by adding 4 mL Buffer 1.
  - b. Centrifuge for 5 minutes at 3900 x g (3900 RCF). Discard the flow through.

6. Label two 15 mL conical tubes W1 and W2. Add the antibody volume calculated in step 4 to the Ultra 430K filter tube, W1. Bring the volume up to 4 mL with Buffer 1. Centrifuge for 10 minutes at 3900 x g (3900 RCF). Save the flow through in conical tube W1.

*Note: This is based on a spectrophotometer. [Ab] can be measured neat if using a nanodrop.*

*E.g.: if  $A_{280} = 0.12$  the concentration is  $(0.12/1.4) * 10 = 0.86$  mg/mL*

*E.g.: To label 200  $\mu$ g of Ab:  $0.20 \text{ mg}/0.86 \text{ mg/mL} = 232 \mu\text{L Ab}$ .*

*E.g.:  $0.232 \text{ mL Ab} + 3.76 \text{ mL Buffer 1} = 4 \text{ mL total volume}$*

7. Transfer the filter containing the antibody into conical tube W2, and bring the volume up to 4 mL with Buffer 1. Centrifuge for 10 minutes at 3900 x g (3900 RCF). Save the flow through in conical tube W2.
8. Mix and remove the concentrated antibody from the filter tube and transfer to an Eppendorf tube. Determine the approximate volume remaining.
9. Calculate the volume of Buffer 1 required bringing the antibody volume back to approximate volume at the start of the buffer exchange.
10. Verify the antibody concentration by reading the absorbance at A280 (using Buffer 1 as the blank). *Note: the concentration should be similar to the initial A280 reading.* After the concentration of antibody is confirmed to be close to the initial amount, discard the W1 and W2 flow through tubes. If antibody yield is low, use W1 and W2 to extract the antibody by using new filters

*Tip: Use the Buffer 1 to rinse the membrane to recover more antibody.*

### Label Capture Antibody

1. Calculate the volume of Capture Label required as follows **Note: the volume of capture label is 6 µL per 1 mg antibody**

$$\text{Capture Label } (\mu\text{L}) = 6 \times \text{Ab (mg)}$$

*E.g.: To label 200 µg of antibody requires (6 x 0.20) = 1.2 µL capture label*

2. Add the calculated volume of Capture Label to the antibody tube. Vortex and incubate for **1 hour** at room temperature.
3. During the incubation, prepare 1X Buffer 3 by diluting the 5 mL of the 10X Buffer 3 into 45 mL of DI water.
4. After incubation, calculate the volume of Buffer 2 required to the labeled antibody to quench the reaction. **Note: the volume of Buffer 2 is 5x the volume of Capture Label.** Mix by vortex.

$$\text{Buffer 2 } (\mu\text{L}) = \text{Capture Label } (\mu\text{L}) \times 5$$

*E.g.: Buffer 2: 1.2 µL x 5 µL = 6 µL.*

5. Prepare the Ultra 4 30K filter tube:
  - a. Pre-wet the Ultra 4 30K filter tube by adding 4 mL Buffer 3.
  - b. Centrifuge for 5 minutes at 3900 x g (3900 R.C.F). Discard the flow through,
6. Label four 15 mL conical tubes W1, W2, W3 and W4.

*Tip: Excess can be added to round up (e.g. 10 µL instead of 6 µL)*



7. Add the antibody volume to the Ultra 4 30K filter tube, W1. Bring the volume up to 4 mL with Buffer 3. Centrifuge for 10 minutes at 3900 x g (3900 RCF). Save the flow through in conical tube W1.
8. Repeat the buffer exchange with 1X Buffer 3 three more times and save the flow through in tubes W2, W3, and W4.
9. Remove the labeled antibody from the filter (W4) and dilute to approximately 1 mg/mL in Buffer 3.
10. Verify the antibody concentration by reading the absorbance at A280 (use 1X Buffer 3 as the blank). *Note: the concentration should be similar to the initial A280 reading.* If the concentration is acceptable, discard W1-W4 flow through.
11. Filter the labeled antibody using the Ultrafree filter and a minicentrifuge for two minutes at maximum speed. Label and store the antibody at 4°C.
12. If desired, add NaN<sub>3</sub> to 0.1% as a preservative.

*Tip: To retain most of the Ab, you can use the undiluted Ab to take the measurement.*

## COATING OF MICROPARTICLE BEADS WITH CAPTURE ANTIBODY

This procedure is for the intended purpose of coating magnetic microparticle (MP) beads with the labeled capture antibody.

### Preparation and coating Microparticle Beads for Conjugation

1. Rotate the uncoated MP bead vial at room temperature for 30 minutes to fully resuspend the beads.
2. During the rotation, prepare 1X Wash Buffer by diluting the 5 mL of the 10X Wash Buffer into 45 mL of DI water.
3. Calculate the volume of labeled capture antibody required to coat at 12.5 µg IgG/mg beads and 25 µg IgG/mg bead. **Note: the concentration of uncoated MP beads provided is 10 mg/mL, therefore, 100 µL is 1 mg of MP beads).**

$$\text{Vol}(\mu\text{L}): = \frac{\text{Capture coating } (\mu\text{g})}{\text{Ab. Conc. (mg/mL)}} \times 1 \text{ mg of MP bead}$$

*E.g.: To label 2 mg of MP bead at 12.5 coating: (12.5/0.86) x 2 = 29 µL capture Ab.  
For 25 ug coating: 29 x 2 = 58 µL of capture label*

4. Wash the MP beads as follows:
  - a. Label two Eppendorf tubes, 12.5 and 25

- b. Place the tube(s) containing the beads in a bench top magnetic bead separator.
  - c. Allow beads to sit in magnet for  $\geq 2$  minutes. Visually confirm that all beads have been pulled toward the magnet.
  - d. Aspirate the supernatant with a pipette and discard.
  - e. Add 1 mL of 1X Wash Buffer to the tube(s). Turn the tube in the magnet holder  $180^\circ$  several times to wash the beads.
  - f. Allow beads to sit in magnet for  $\geq 2$  minutes, then remove and discard the Wash Buffer.
  - g. Repeat the wash step two more times using fresh 1X Wash Buffer
  - h. Aspirate and discard the 1X Wash Buffer.
5. Add the calculated volume of capture antibody to the each of the tubes, followed by the calculated volume of wash buffer based on the mg of beads being labeled.

$$\text{Total Vol}(\mu\text{L}) = \text{Capture} (\mu\text{L}) + \text{Wash Buffer} (\mu\text{L})$$

*E.g.: To label 2 mg of MP bead required at 12.5 (200 – 29) = 171  $\mu\text{L}$  Wash Buffer plus 29  $\mu\text{L}$  capture Ab totaling 200  $\mu\text{L}$*

6. Incubate for **60 minutes** at room temperature with gentle inversion on a rotator.
7. After incubation, wash the beads by placing in a bead magnetic separator and repeat wash step 4, five times.
8. After final wash resuspend antibody coated MP beads in Coated Bead Buffer as follows:
  - a. After final wash in above step allow beads to sit in magnet for  $\geq 2$  minutes.
  - b. Visually confirm that all MPs have been pulled to the magnet.
  - c. Aspirate the supernatant with a pipette and discard.
  - d. Add equal amount of Coated Bead Buffer to resuspend and block the coated beads at a 10 mg/mL.
  - e. Either block the coated beads for 1hr at RT or overnight at  $4^\circ\text{C}$  rotating.
9. Store the blocked beads at  $4^\circ\text{C}$  until use.

*Eg: 2 mg (200  $\mu\text{L}$ ) of beads requires 200  $\mu\text{L}$  of Coated Bead Buffer*

# WORKSHEET

See preceding protocol for more detailed information on each step.

## Pre-Label

Step	Result
Dilute Antibody to ~1 mg/mL in Buffer 1	Vol. added: _____ mL
Read A280 of antibody	A280: _____ Conc.: (C) _____ mg/mL
Determine amount and volume of antibody to label	Ab Amount (A): _____ mg (up to 1 mg) Volume.: (A/C) _____ mL
Rinse Ultra-4 30K filter tube with 4 mL Buffer 1, 5 minutes at 3900 RCF	
Add Antibody to Ultra-4 30K filter tube	
Wash #1 with Buffer 1 (10 min.)	Ab. Vol.: _____ mL + _____ mL Buffer = 4 mL
Wash #2 with Buffer 1 (10 min.)	Up to 4 mL Buffer 1
Dilute to ~1.0 mg/mL in Buffer 1	Recovered Vol: _____ mL Buffer 1 added: _____ mL
Read A280 of antibody	A280: _____ Conc.: (C) _____ mg/mL

## Labeling

Step	Result
Add capture Label. Mix by vortexing.	Capture label added = _____ $\mu$ L
Incubate 1 hour at RT.	Start time: _____ End Time: _____

## Post-label

Step	Result
Rinse new Ultra 4 30K filter tube with 4 mL 1x Buffer 3, 5 minutes at 3900 RCF	
Add labeled Ab to Ultra-4 30K filter tube	Ab vol: _____ mL
Wash #1 with 1X Buffer 3 (10 min.)	Ab vol: _____ mL + _____ mL Buffer 3 = 4mL
Wash #2, #3, #4 with 1X Buffer 3, (10 min.)	Up to 4 mL Buffer 3
Dilute the antibody to approx. 1 mg/mL with 1X Buffer 3	Buffer 3 added: _____ mL
Read A280 of antibody	A280: _____ Conc: (C) _____ mg/mL

## MP Conjugation

Step	Result
Wash uncoated beads with 1x Wash Buffer 3 times.	
Add calculated volume of Antibody to coat at 12.5 and 25 µg/IgG/mg bead.	Ab vol. added to coat 12.5 µg: _____ µL Ab vol. added to coat 25 µg: _____ µL
Incubate 1 hour at RT, rotating.	Start time: _____ End Time: _____
Wash beads with 1x Wash Buffer 5 times	
Add calculated amount of coated bead buffer to block beads at RT for 1hr, or overnight at 4°C rotating.	Start time: _____ End Time: _____

## ORDERING INFORMATION

To place an order or to obtain additional information about SMC™ products, please contact your Customer Service or Technical Support Specialist.

Contact information for each region can be found on our website:

[emdmillipore.com/contact](http://emdmillipore.com/contact)

### **Conditions of Sale**

For Research Use Only. Not for Use in Diagnostic Procedures.

### **Safety Data Sheets (SDS)**

Safety Data Sheets for EMD Millipore products may be ordered by fax or phone or through our website at [emdmillipore.com/msds](http://emdmillipore.com/msds)



EMD Millipore Corporation  
3050 Spruce Street  
St. Louis, Missouri 63103  
United States of America

Toll-Free US: (800) 645-5476  
Fax: (800) 645-5439  
[emdmillipore.com](http://emdmillipore.com)