

User Guide

# SMC® Capture Labeling Kit Instructions

## Microparticle Assay

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

**03-0077-02**

Introduction.....	3	Worksheet .....	13
Reagents.....	4	Pre-Label .....	13
Reagents Provided.....	4	Labeling .....	14
Storage Instructions .....	5	Post-label .....	14
General Supplies Required .....	5	MP Conjugation .....	15
Precautions.....	6	Product Ordering .....	15
Capture Antibody Labeling.....	7	Notice .....	16
Preparation of Capture Antibody		Technical Assistance.....	16
for Labeling .....	7	Terms and Conditions of Sale .....	16
Label Capture Antibody .....	9	Safety Data Sheets (SDS).....	16
Prepare the Ultra-4 30K		Contact Information .....	16
filter tube .....	10		
Coating of Microparticle Beads with			
Capture Antibody .....	11		
Preparation and Coating			
Microparticle Beads for Conjugation			
.....	11		

---

## Introduction

The SMC<sup>®</sup> Immunoassay System 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<sup>®</sup> Immunoassay System.

This kit should be used in conjunction with (03-0076-02) SMC<sup>®</sup> Detection Labeling Kit and (03-0078-00) SMC<sup>®</sup> 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 [SigmaAldrich.com/techservice](https://www.sigmaaldrich.com/techservice).

**This kit is for Research Use Only. Not for Use in Diagnostic Procedures.**

## Reagents

The SMC<sup>®</sup> 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 (Not Provided). All reagents supplied are for Research Use Only.

### Reagents Provided

All products are available at [SigmaAldrich.com](https://www.sigmaaldrich.com) unless otherwise noted.

Description	Storage Conditions	Packaging Details	Component Number
Capture Label	≤ -70 °C	1 x 20 µL	02-0557-00
Buffer 1	2-8 °C	1 x 25 mL	02-0552-00
Buffer 2	2-8 °C	1 x 500 µL	02-0553-00
Buffer 3	2-8 °C	1 x 5 mL	02-0554-00
Filter Tube Ultrafree <sup>®</sup>	2-8 °C	2 pack	02-0555-00
Amicon <sup>®</sup> Ultra-4 Filter Tube with Ultracel <sup>®</sup> -30 membrane	2-8 °C	2 pack	02-0556-00
10X Wash Buffer	2-8 °C	1 x 50 mL	02-0001-03
Coated Bead Buffer	2-8 °C	1 x 5 mL	02-0558-00
Uncoated SMC <sup>®</sup> Beads	2-8 °C	1 x 4 mL	02-0559-00

**Note:** Additional Amicon<sup>®</sup> Ultra-4 Centrifugal Filter Units with Ultracel<sup>®</sup>-30 membrane can be purchased from [SigmaAldrich.com](https://www.sigmaaldrich.com) (Cat. No. UFC803008).

---

## Storage Instructions

- The SMC<sup>®</sup> 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 with Stericup<sup>®</sup> Filter 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.

## Additional Supplies Required (Not Provided)

- De-ionized or distilled water (DI Water)
- Pipettes capable of transferring 5 µL-5 mL
- Polypropylene Eppendorf<sup>®</sup> 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 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. No.	Label	
10X Wash Buffer	02-0001-03		<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.
Capture Label	02-0557-00		<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:
  - Prepare 100  $\mu$ L of a 1:10 dilution of the 1 mg/mL antibody solution in Buffer 1.
  - Note:** This is based on a spectrophotometer. Ab can be measured neat if using a nanodrop.
  - Measure the absorbance (using Buffer 1 as a blank).
  - Using Beer's Law  $A = \epsilon Lc$ , calculate the antibody concentration as shown below. The example assumes a 1 cm path length and a molar extinction coefficient of 1.4:

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

(For example, if  $A_{280} = 0.12$  the concentration is  $(0.12/1.4) * 10 = 0.86$  mg/mL)

4. Use the calculated antibody concentration to calculate the volume of antibody solution needed to label 0.2-1 mg Antibody. (For example, to label 200  $\mu$ g of Ab:  $0.20 \text{ mg} / 0.86 \text{ mg/mL} = 232 \mu\text{L Ab}$ .)
5. Prepare the Ultra-4 30K filter tube:
  - Pre-wet the Ultra-4 30K filter tube by adding 4 mL Buffer 1.
  - Centrifuge for 5 minutes at  $3900 \times g$  (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-4 30K filter tube, W1. Bring the volume up to 4 mL with Buffer 1. Centrifuge for 10 minutes at  $3900 \times g$  (3900 RCF). Save the flow through in conical tube W1.

(For example,  $0.232 \text{ mL Ab} + 3.76 \text{ mL Buffer 1} = 4 \text{ mL total volume}$ .)

- 
8. 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.
  9. Mix and remove the concentrated antibody from the filter tube and transfer to an Eppendorf® Tube. Determine the approximate volume remaining.  
**Tip:** Use the Buffer 1 to rinse the membrane to recover more antibody.
  10. Calculate the volume of Buffer 1 required bringing the antibody volume back to approximate volume at the start of the buffer exchange.
  11. 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.



## Label Capture Antibody

Calculate the volume of Capture Label required as follows:

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

**Note:** The volume of capture label is 6  $\mu\text{L}$  per 1 mg antibody.

For example, to label 200  $\mu\text{g}$  of antibody requires  $(6 \times 0.20) = 1.2 \mu\text{L}$  capture label.

1. Add the calculated volume of Capture Label to the antibody tube. Vortex and incubate for 1 hour at room temperature.
2. During the incubation, prepare 1X Buffer 3 by diluting the 5 mL of the 10X Buffer 3 into 45 mL of DI water.
3. 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. (For example, Buffer 2:  $1.2 \mu\text{L} \times 5 \mu\text{L} = 6 \mu\text{L}$ )

**Tip:** Excess can be added to round up (For example, 10  $\mu\text{L}$  instead of 6  $\mu\text{L}$ ).

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

---

## Prepare the Ultra-4 30K filter tube

1. Pre-wet the Ultra-4 30K filter tube by adding 4 mL Buffer 3.
2. Centrifuge for 5 minutes at  $3900 \times g$  (3900 RCF). Discard the flow through.
3. Label four 15 mL conical tubes W1, W2, W3 and W4.
4. 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  $3900 \times g$  (3900 RCF). Save the flow through in conical tube W1.
5. Repeat the buffer exchange with 1X Buffer 3 three more times and save the flow through in tubes W2, W3, and W4.
6. Remove the labeled antibody from the filter (W4) and dilute to approximately 1 mg/mL in Buffer 3.
7. 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.

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

8. 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.
9. If desired, add  $\text{NaN}_3$  to 0.1% as a preservative.

# 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  $\mu\text{g}$  IgG/mg beads and 25  $\mu\text{g}$  IgG/mg bead.

**Note:** the concentration of uncoated MP beads provided is 10 mg/mL, therefore, 100  $\mu\text{L}$  is 1 mg of MP beads.

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

(For example: To label 2 mg of MP bead at 12.5  $\mu\text{g}$  coating,  $(12.5/0.86) \times 2 = 29 \mu\text{L}$  capture Ab. For 25  $\mu\text{g}$  coating,  $29 \times 2 = 58 \mu\text{L}$  of capture label.)

Wash the MP beads as follows:

4. Label two Eppendorf® Tubes, 12.5  $\mu\text{g}$  and 25  $\mu\text{g}$ .
5. Place the tube(s) containing the beads in a bench top magnetic bead separator.
6. Allow beads to sit in magnet for  $\geq 2$  minutes. Visually confirm that all beads have been pulled toward the magnet.
7. Aspirate the supernatant with a pipette and discard.
8. Add 1 mL of 1X Wash Buffer to the tube(s). Turn the tube in the magnet holder 180° several times to wash the beads.
9. Allow beads to sit in magnet for  $\geq 2$  minutes, then remove and discard the Wash Buffer.
10. Repeat the wash step two more times using fresh 1X Wash Buffer.
11. Aspirate and discard the 1X Wash Buffer.

12. Add the calculated volume of capture antibody to each of the tubes, followed by the calculated volume of wash buffer based on the mg of beads being labeled.

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

(For example: To label 2 mg of MP bead at 12.5  $\mu\text{g}$ , 200  $\mu\text{L}$  MP bead – 29  $\mu\text{L}$  capture antibody = 171  $\mu\text{L}$  Wash Buffer.)

13. Incubate for 60 minutes at room temperature with gentle inversion on a rotator.
14. After incubation, wash the beads by placing in a bead magnetic separator and repeat wash step 4, five times.

After final wash resuspend antibody coated MP beads in Coated Bead Buffer as follows:

15. After final wash in above step allow beads to sit in magnet for  $\geq 2$  minutes.
16. Visually confirm that all MPs have been pulled to the magnet.
17. Aspirate the supernatant with a pipette and discard.
18. Add equal amount of Coated Bead Buffer to resuspend and block the coated beads at a 10 mg/mL. For example, 2 mg (200  $\mu\text{L}$ ) of beads requires 200  $\mu\text{L}$  of Coated Bead Buffer.
19. Either block the coated beads for 1hr at RT or overnight at 4  $^{\circ}\text{C}$  rotating.
20. Store the blocked beads at 4  $^{\circ}\text{C}$  until use.

## 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

For research use only. Not for use in diagnostic procedures.

## 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 = 4 mL
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 hr 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 1 hr, or overnight at 4 °C rotating.	Start time: _____ End Time: _____

## Product Ordering

Products are available for online ordering at [SigmaAldrich.com](https://www.sigmaaldrich.com).

---

## Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

### Technical Assistance

Visit the tech service page at [SigmaAldrich.com/techservice](https://SigmaAldrich.com/techservice).

### Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/terms](https://SigmaAldrich.com/terms).

### Safety Data Sheets (SDS)

Safety Data Sheets are available on the product page at [SigmaAldrich.com](https://SigmaAldrich.com).

### Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/offices](https://SigmaAldrich.com/offices).

The life science business of Merck KGaA, Darmstadt, Germany operates as MilliporeSigma in the U.S. and Canada.

MilliporeSigma, Amicon, Ultrafree, Ultracel, Stericup, Millipore and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.

© 2021-2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

For research use only. Not for use in diagnostic procedures.