

# SiteClick™ Antibody Labeling Kits

Catalog Number S10449, S10450, S10451, S10452, S10453, S10454, S10455, S10467, S10469

Pub. No. MAN0007752 Rev. A.0



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from [thermofisher.com/support](http://thermofisher.com/support).

## Product description

The SiteClick™ Antibody Labeling Kits allow you to conjugate your own antibodies to DIBO-modified Qdot™ nanocrystals (525, 565, 585, 605, 625, 655, 705, and 800-nm emission) or DIBO-modified R-Phycoerythrin (R-PE). The SiteClick™ conjugation workflow consists of three steps (antibody carbohydrate domain modification, azide attachment to the antibody, and conjugation with the DIBO-modified label) and relies on copper-free click chemistry to covalently link the label containing the DIBO moiety with the azide-modified antibody without reducing the protein. The antibody concentrators provided in the kits are used to purify and concentrate the antibody at each step of the SiteClick™ antibody labeling workflow (Figure 1 on page 3).

In the first step of SiteClick™ conjugation, terminal galactose residues on the N-linked sugars in the Fc region of the antibody are removed by β-Galactosidase. The azide-containing sugar, GalNAz, is then added to the modified carbohydrate domain of the antibody via the β-1,4-galactosyltransferase (Gal-T)-catalyzed reaction targeting the terminal GlcNAc residues. This specific targeting maintains the integrity of the antigen binding site on the antibody. Finally, the antibody (now containing an azide moiety) is conjugated to the DIBO-modified label (Qdot™ nanocrystals or R-PE) in a copper-free click reaction with simple overnight incubation (Figure 2 on page 4).

Each SiteClick™ Antibody Labeling Kit contains sufficient reagents to perform one conjugation reaction of Qdot™ nanocrystals or R-PE to a primary IgG antibody sample. The protocol in this manual describes a conjugation reaction starting with 100–125 µg of whole IgG from any host species.

**Table 1 Optimal fluorescence excitation and emission maxima of DIBO-modified labels**

| Label           | Excitation (nm) <sup>[1]</sup>  | Emission (nm) <sup>[1]</sup> | Extinction coefficient (ε) (M <sup>-1</sup> cm <sup>-1</sup> ) | Measured at (nm)          | Cat. No.               |
|-----------------|---------------------------------|------------------------------|--|---------------------------|------------------------|
| Qdot™ 525       | <525 nm                         | 525 nm                       | 200,000 M <sup>-1</sup> cm <sup>-1</sup>                       | between 504 nm and 512 nm | <a href="#">S10449</a> |
| Qdot™ 565       | <565 nm                         | 565 nm                       | 300,000 M <sup>-1</sup> cm <sup>-1</sup>                       | between 548 nm and 556 nm | <a href="#">S10450</a> |
| Qdot™ 585       | <585 nm                         | 585 nm                       | 250,000 M <sup>-1</sup> cm <sup>-1</sup>                       | between 572 nm and 580 nm | <a href="#">S10451</a> |
| Qdot™ 605       | <605 nm                         | 605 nm                       | 400,000 M <sup>-1</sup> cm <sup>-1</sup>                       | between 592 nm and 600 nm | <a href="#">S10469</a> |
| Qdot™ 625       | <625 nm                         | 625 nm                       | 500,000 M <sup>-1</sup> cm <sup>-1</sup>                       | between 605 nm and 612 nm | <a href="#">S10452</a> |
| Qdot™ 655       | <655 nm                         | 655 nm                       | 1,700,000 M <sup>-1</sup> cm <sup>-1</sup>                     | 550 nm                    | <a href="#">S10453</a> |
| Qdot™ 705       | <705 nm                         | 705 nm                       | 1,700,000 M <sup>-1</sup> cm <sup>-1</sup>                     | 550 nm                    | <a href="#">S10454</a> |
| Qdot™ 800       | <800 nm                         | 800 nm                       | 1,700,000 M <sup>-1</sup> cm <sup>-1</sup>                     | 550 nm                    | <a href="#">S10455</a> |
| R-Phycoerythrin | 496, 546, 565 nm <sup>[2]</sup> | 578 nm                       | 1,960,000 M <sup>-1</sup> cm <sup>-1</sup>                     | 578 nm                    | <a href="#">S10467</a> |

<sup>[1]</sup> Qdot™ nanocrystals are excitable (Ex, in nm) at any wavelength below their emission maxima (Em, in nm). For most practical applications, they can be excited at wavelengths below 405 nm.

<sup>[2]</sup> Multiple absorbance peaks

## Contents and storage

**Table 2 SiteClick™ Antibody Labeling Kit contents (Cat. Nos. [S10449](#), [S10450](#), [S10451](#), [S10452](#), [S10453](#), [S10454](#), [S10455](#), [S10467](#), [S10469](#))**

| Component  | Cap color          | Amount                              | Storage <sup>[1]</sup>                          |
|--|--------------------|-------------------------------------|---|
| Antibody preparation buffer (Component A)              | Yellow             | 1.8 mL                              | 2–8°C<br>Do not freeze, and protect from light. |
| Antibody concentrator (small) (Component B)            | N/A <sup>[2]</sup> | each                                |   |
| Collection tube (Component C)                          | N/A                | each                                |   |
| β-Galactosidase (Component D)                          | Green              | 12 μL                               |   |
| UDP-GalNAz (Component E)                               | Blue               | 220 μg                              |   |
| 20X Tris pH 7.0 (Component F)                          | Red                | 1.8 mL                              |   |
| Buffer additive (Component G)                          | Purple             | 30 μL                               |   |
| β-1,4-galactosyltransferase (GalT) (Component H)       | Orange             | 88 μL                               |   |
| Antibody concentrator (large) (Component I)            | N/A                | each                                |   |
| DIBO-modified label (Component J) <sup>[3]</sup>       | Dark Orange        | 55 μL (Qdot™)<br>or<br>80 μL (R-PE) |   |
| Purification concentrator (Component K) <sup>[4]</sup> | N/A                | each                                |   |

<sup>[1]</sup> When stored as directed, this kit is stable for at least 3 months.

<sup>[2]</sup> N/A = not applicable.

<sup>[3]</sup> DIBO-modified Qdot™ nanocrystal or DIBO-modified R-Phycoerythrin (R-PE). See Table 1 for the approximate fluorescence excitation and emission maxima.

<sup>[4]</sup> Only available with SiteClick™ Qdot™ 605 Antibody Labeling Kit (Cat. No. [S10469](#)).

## Required materials not supplied

Unless otherwise indicated, all materials are available through [thermofisher.com](http://thermofisher.com). "MLS" indicates that the material is available from [fisherscientific.com](http://fisherscientific.com) or another major laboratory supplier. Catalog numbers that appear as links open the web pages for those products.

| Item   | Source              |
|--|---------------------|
| <b>Equipment</b>   |                     |
| Centrifuge with fixed angle rotor that can accommodate 1.5-mL centrifuge tubes   | <a href="#">MLS</a> |
| Centrifuge with swinging bucket or fixed angle rotor with 17 mm × 100 mm inserts | <a href="#">MLS</a> |
| <b>Reagents and consumables</b>  |                     |
| 100 to 125 μg of whole IgG antibody <sup>[1]</sup>                               | <a href="#">MLS</a> |
| Distilled water (dH <sub>2</sub> O)  | <a href="#">MLS</a> |
| Centrifuge tubes: 1.5-mL and 15-mL   | <a href="#">MLS</a> |

<sup>[1]</sup> Preferably at a concentration of 2 to 4 mg/mL in a Tris-based buffer, free of carrier proteins and/or azide.

## Workflow

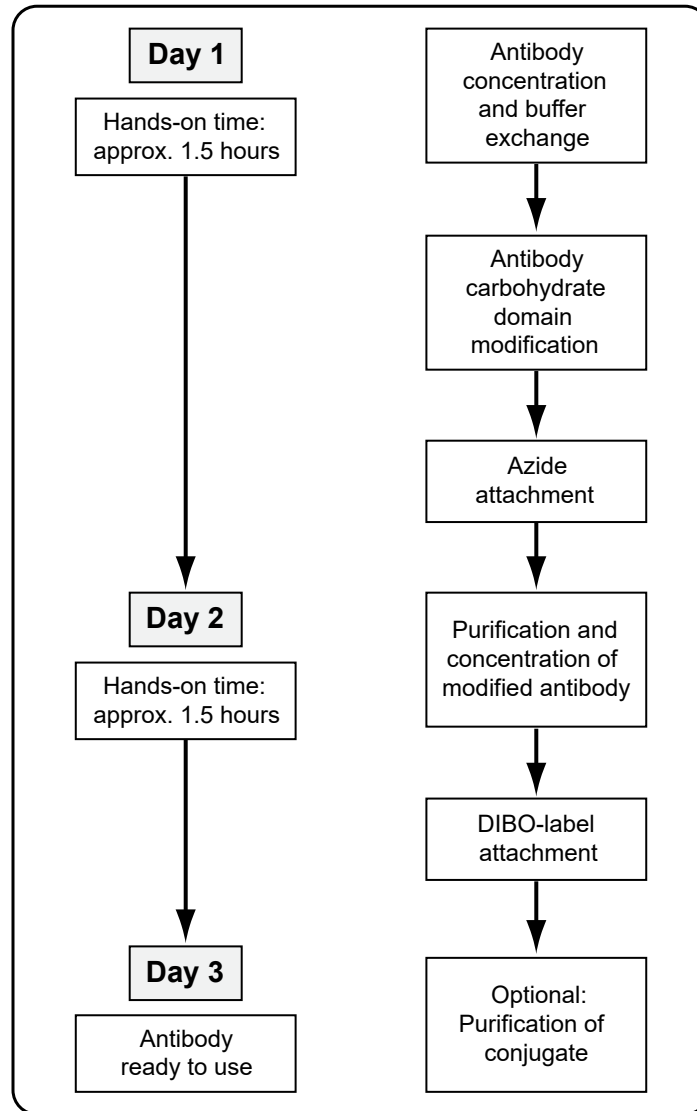


Figure 1 SiteClick™ antibody labeling workflow

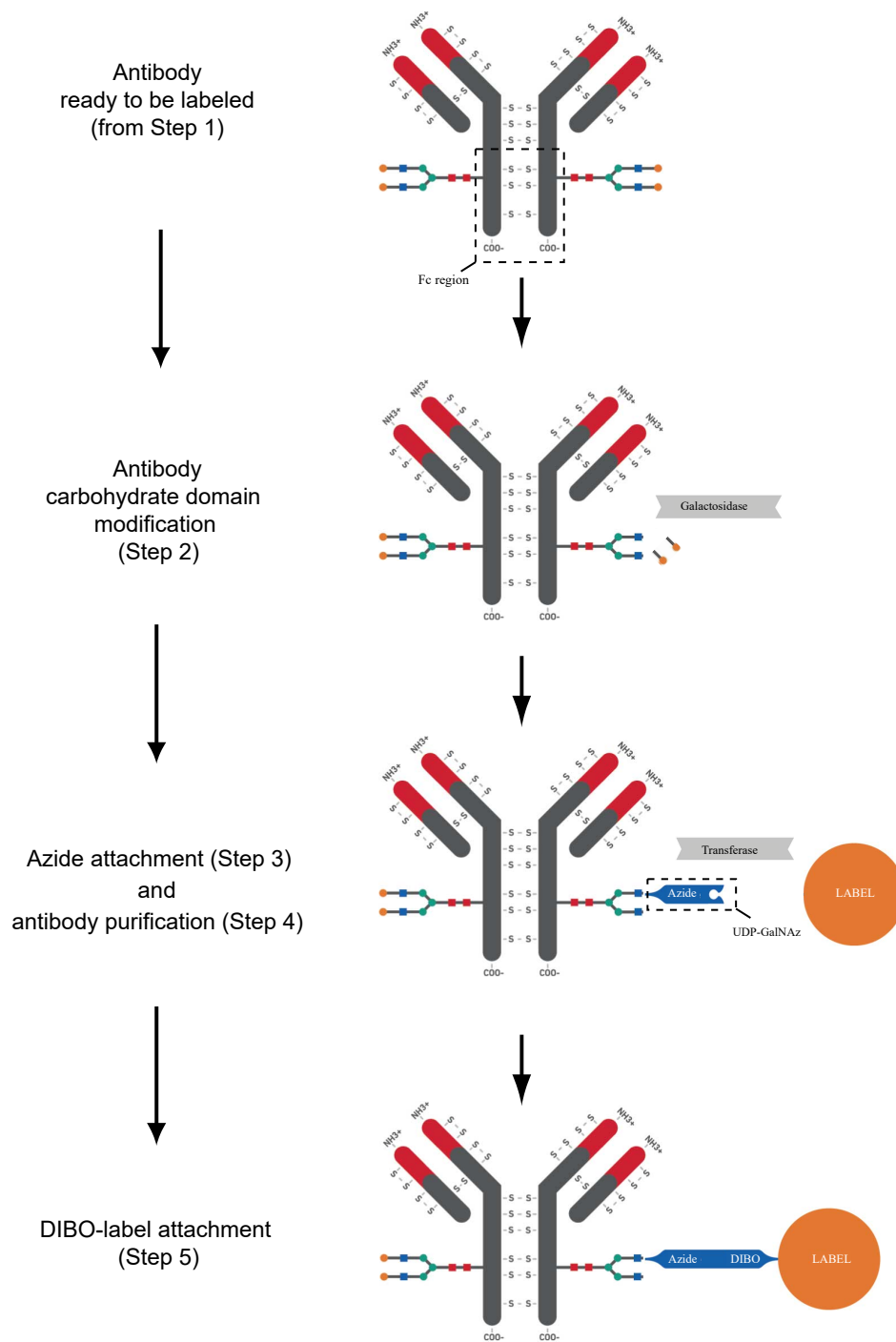


Figure 2 SiteClick™ conjugation reaction

### Caution

- $\beta$ -Galactosidase (Component D) can cause an allergic skin reaction, and it can cause allergy or asthma symptoms or breathing difficulties, if inhaled.
- The Qdot™ conjugate contains cadmium and selenium in an inorganic crystalline form.
- Discard the reagents in compliance with all pertaining local regulations.
- If there is contact with the eyes, rinse the eyes immediately with plenty of water and seek medical advice.
- Always wear appropriate laboratory protective clothing and gloves when handling these reagents.

## Step 1. Antibody concentration and/or Buffer exchange (Optional)

**Time required:** 1 hour

Perform the antibody concentration and buffer exchange step if:

- The antibody concentration is less than 2 mg/mL, and/or
- The antibody is in a phosphate-based buffer (e.g. PBS), and/or
- The antibody is in a buffer containing azide.

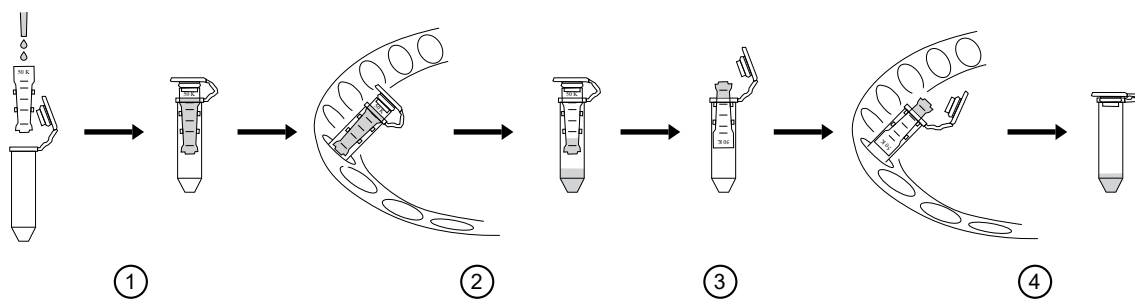
Before you start, briefly centrifuge Components A, C, D, E, F, G, H, and J to ensure that all material is at the bottom of the tubes.

### Wash the antibody concentrator

1. Add 450  $\mu\text{L}$  of  $\text{dH}_2\text{O}$  to the small antibody concentrator (Component B), then cap the device (see Figure 3).
2. Centrifuge for 6 minutes at  $5000 \times g$ .  
**Note:** Ensure that the cap strap and one membrane panel of the concentrator faces the center of the rotor.
3. Discard the flow-through.

### Concentrate the antibody and the exchange buffer

1. Add a sufficient volume of antibody solution to contain 100–125  $\mu\text{g}$  of antibody to the small antibody concentrator. For example, if the antibody concentration is 1 mg/mL, add 125  $\mu\text{L}$ .
2. Dilute the added antibody to 500  $\mu\text{L}$  using the antibody preparation buffer (Component A).
3. Centrifuge for 6 minutes at  $5000 \times g$ . Ensure that the cap strap and one membrane panel of the concentrator faces the center of the rotor.
4. Discard the flow-through.
5. Add 450  $\mu\text{L}$  of antibody preparation buffer (Component A) to the small antibody concentrator (Component B), then centrifuge for 6 minutes at  $5000 \times g$ . Ensure that the cap strap and one membrane panel of the concentrator faces the center of the rotor.  
**Note:** If antibody volume in concentrator is greater than 50  $\mu\text{L}$  after Step 5, centrifuge for an extra 3 minutes at  $5000 \times g$  or until the appropriate volume is achieved.
6. Invert the small antibody concentrator (Component B) into the collection tube (Component C) (see Figure 3).
7. Centrifuge for 3 minutes at  $1000 \times g$  to collect the concentrated antibody. After collection, you should have approximately 50  $\mu\text{L}$  of concentrated antibody in the collection tube.



**Figure 3** Antibody concentration and/or buffer exchange

- ① Add antibody and antibody preparation buffer to antibody concentrator
- ② Centrifuge

- ③ Invert the antibody concentrator into collection tube
- ④ Centrifuge to collect the concentrated antibody

## Step 2. Modification of the antibody carbohydrate domain

**Time required:** 4 hours, hands-off

### Add $\beta$ -galactosidase

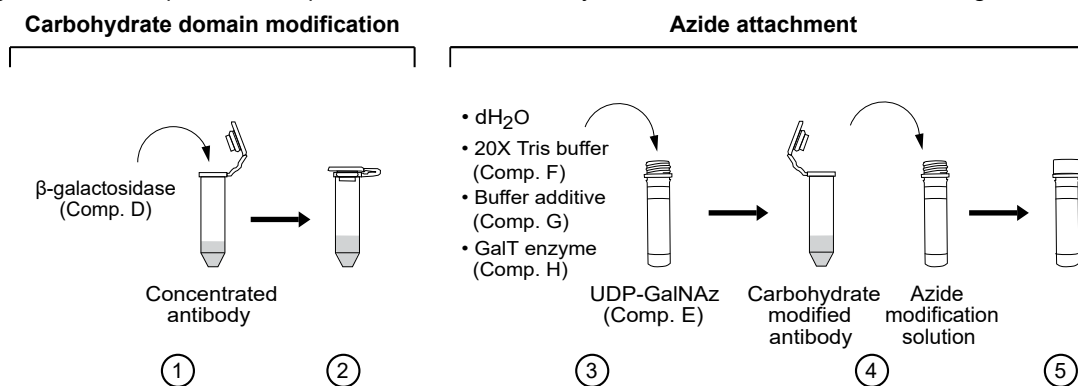
1. Add 10  $\mu$ L of  $\beta$ -galactosidase (Component D) to the antibody collected in step 7 on page 5 (see Figure 4).
2. Wrap the tube cap with Parafilm™ laboratory film or similar, then incubate for 4 hours at 37°C.

## Step 3. Azide attachment

**Time required:** 5 minutes hands-on, then overnight incubation

### Add GalT enzyme

1. Prepare the azide modification solution by adding the following components to the tube containing UDP-GalNAz (Component E) (see Figure 4):
  - 75  $\mu$ L of dH<sub>2</sub>O
  - 10  $\mu$ L of 20X Tris buffer, pH 7.0 (Component F)
  - 25  $\mu$ L of buffer additive (Component G)
  - 80  $\mu$ L of GalT enzyme (Component H)
2. Vortex the reaction components, then add the modified antibody (from “Add  $\beta$ -galactosidase”, step 2 on page 6) to the tube.
3. Briefly centrifuge the tube, wrap the tube cap with Parafilm™ laboratory film or similar, then incubate overnight at 30°C.



**Figure 4 Modification of antibody carbohydrate domain and azide attachment**

- ① Add  $\beta$ -galactosidase to the concentrated antibody
- ② Incubate for 4 hours
- ③ Prepare the azide modification solution
- ④ Add the azide modification solution to the modified antibody
- ⑤ Incubate overnight

## Step 4. Purification and concentration of the azide-modified antibody

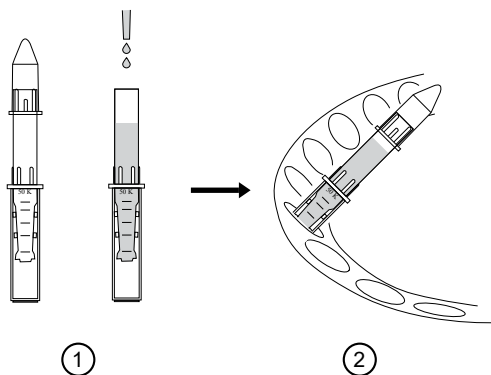
**Time required:** 1 hour

### Wash the antibody concentrator

1. Prepare 10 mL of 1X Tris, pH 7.0 by adding 500  $\mu$ L of 20X Tris, pH 7.0 (Component F) to 9.5 mL of dH<sub>2</sub>O in a 15-mL conical tube. Vortex briefly to mix.  
**Note:** You can also use TBS for the purification and collection of the modified antibody. 20X Tris, pH 7.0 is provided for convenience.
2. Remove the conical collection tube from the large antibody concentrator (Component I) (see Figure 5).
3. Add 1 mL of 1X Tris, pH 7.0 to the large antibody concentrator (Component I), then centrifuge for 10 minutes at 1200  $\times$  g. Ensure that one membrane panel of the concentrator faces the center of the rotor.
4. Discard the flow-through

## Purify the azide-modified antibody

1. Add 1.75 mL of 1X Tris, pH 7.0 and 250  $\mu$ L of the azide-modified antibody (from “Add GalT enzyme”, step 3 on page 6) to the large antibody concentrator (Component I) (see Figure 5).
2. Centrifuge for 6 minutes at 1200  $\times$  *g*. Ensure that one membrane panel of the concentrator faces the center of the rotor.
3. Discard the flow-through.
4. Add 1.8 mL of 1X Tris, pH 7.0 to the large antibody concentrator (Component I), then centrifuge for 10 minutes at 1200  $\times$  *g*. Ensure that one membrane panel of the concentrator faces the center of the rotor.
5. Discard the flow-through, then repeat step 4.



**Figure 5 Purification and concentration of azide-modified antibody**

- ① Add azide-modified antibody to the large antibody concentrator      ② Centrifuge

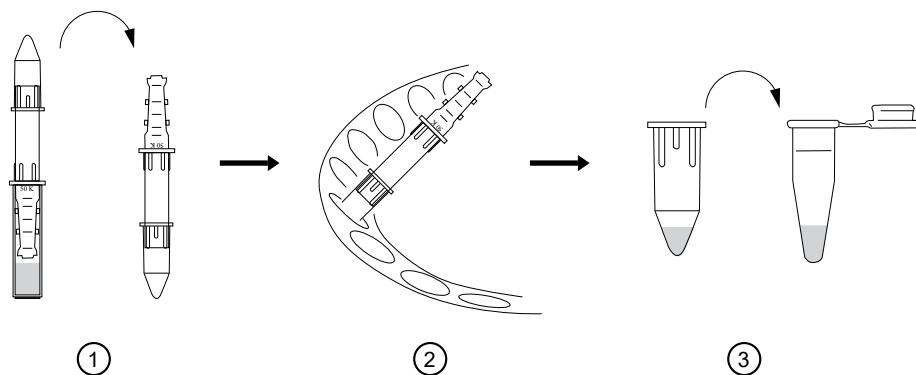
## Concentrate and collect the purified azide-modified antibody

1. Add 1.8 mL of 1X Tris, pH 7.0 to the large antibody concentrator (Component I), then centrifuge for 10 minutes at 1400  $\times$  *g*. Discard the flow-through. The final volume in the concentrator should be approximately 80–120  $\mu$ L.

**Note:** If the antibody volume in the concentrator is greater than 100  $\mu$ L, centrifuge for an extra 5 minutes at 1400  $\times$  *g* or until the appropriate volume is achieved.

2. Invert the antibody concentrator into the conical collection tube (see Figure 6).
3. Centrifuge for 3 minutes at 1000  $\times$  *g* to collect the concentrated antibody.
4. Transfer the antibody from the conical collection tube to a 1.5-mL centrifuge tube. If the final collected volume is less than 100  $\mu$ L, dilute the antibody to 100  $\mu$ L with 1X Tris, pH 7.0.

**Note:** At this stage, you can store the antibody at 2–8°C and attach the DIBO-modified label at a later time.



**Figure 6 Collection of the purified and concentrated azide-modified antibody**

- ① Invert the antibody concentrator      ③ Collect the concentrated antibody  
② Centrifuge

## Step 5. Conjugation with DIBO-modified label

**Time required:** 5 minutes hands-on, then overnight incubation

### Add DIBO-modified label

1. Add the DIBO-modified label (Component J) to the azide-modified antibody in the 1.5-mL centrifuge tube:
  - If using DIBO-modified Qdot™ nanocrystal, add 50  $\mu$ L of Qdot™ DIBO (Component J).
  - If using DIBO-modified R-PE, add 75  $\mu$ L of R-PE DIBO (Component J).
2. Vortex the reaction mixture, briefly centrifuge, then incubate overnight at 25°C.
3. Store the antibody conjugate at 2–8°C, protected from light (see “Antibody conjugate storage” on page 9).

**Note:** (*Optional*) If desired, you can further purify the antibody conjugate of excess unconjugated antibody (“Step 6. Purification and concentration of the antibody conjugate (optional)” on page 8).

## Step 6. Purification and concentration of the antibody conjugate (*optional*)

**Time required:** 1 hour

The purification step removes any excess antibody that has not been conjugated with the Qdot™ DIBO-modified label as well as unconjugated Qdot™ nanocrystals. Typical yield from this step is approximately 80%.

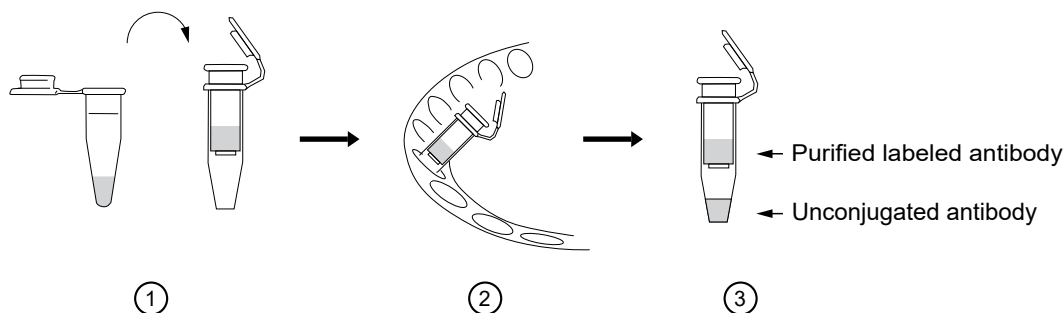
**Note:** The flow-through from the final concentration and purification step can be colored, and this color can vary. This is normal and does not affect the quality of the final purified conjugate.

**Note:** For many applications, purification of the antibody conjugate is not necessary, because the efficiency of the SiteClick™ conjugation reaction yields a minimal quantity of free antibody.

**Note:** The purification concentrator (Component K) is only available in SiteClick™ antibody purification kits containing DIBO-modified Qdot™ nanocrystals with emission maxima of 605, 625, 655, 705, and 800 nm. Antibody conjugates labeled with R-PE or with Qdot™ nanocrystals that emit at 525, 565, or 585 nm are too small to be retained by the membrane in the purification concentrator.

### Purify and concentrate the antibody conjugate

1. Add 500  $\mu$ L of dH<sub>2</sub>O to the purification concentrator (Component K), then centrifuge for 5 minutes at 1500  $\times$  g. Discard the flow-through.
2. Add 150  $\mu$ L of the labeled antibody conjugate (from step 2 on page 8) to the purification concentrator (Component K) (see Figure 7).
3. Add buffer (TBS or PBS) to the purification concentrator to bring the volume to 500  $\mu$ L, then centrifuge for 10 minutes at 1500  $\times$  g. Discard the flow-through.
4. Add 500  $\mu$ L of buffer (TBS or PBS) to the purification concentrator, then centrifuge for 10 minutes at 1500  $\times$  g.
5. Discard the flow-through, then repeat step 4.
6. Collect the purified labeled antibody conjugate from the top of the concentrator, then transfer to a new 1.5-mL centrifuge tube. You can store the conjugated antibody at 2–8°C, protected from light (see “Antibody conjugate storage” on page 9).



**Figure 7** Optional purification of the conjugated antibody

- ① Add the labeled antibody conjugate to purification concentrator
- ② Centrifuge
- ③ Collect the purified labeled antibody conjugate



## Storage and use of the antibody conjugates

### Antibody conjugate storage

Store the antibody conjugate labeled with the Qdot™ nanocrystal or R-PE at 2–8°C, in the dark, until needed. **Do not freeze.** You can add sodium azide or thimerosal at this stage to a final concentration of 0.02% (w/v) for long term storage, if preferred.

### Antibody conjugate use

You can determine the concentration of the label in the final preparation by measuring the optical density of the label at the specified wavelength, then using the formula  $A = \epsilon cL$ , where A is the absorbance,  $\epsilon$  is the molar extinction coefficient (Table 1 on page 1), c is the molar concentration, and L is the pathlength.

For example, for the Qdot™ 605 nanocrystal, if the material eluting from the final column has  $A = 0.80$  measured at peak between 592 nm and 600 nm in a cuvette with 1 cm pathlength, then  $c = A/\epsilon = 0.80/400,000 = 2 \mu\text{M}$  of label, based on nanocrystal absorbance.

Determine the optimal concentration of any labeled antibody empirically for a particular application or experiment. Determine the optimal working concentrations by performing a titration series for the application of interest.

## 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 at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



Life Technologies Corporation | 29851 Willow Creek Road | Eugene, Oregon 97402 USA

For descriptions of symbols on product labels or product documents, go to [thermofisher.com/symbols-definition](http://thermofisher.com/symbols-definition).

Revision history: Pub. No. MAN007752 A.0

| Revision | Date         | Description   |
|----------|--------------|---|
| A.0      | 20 June 2023 | Correct the concentration of Tris buffer needed to dilute the antibody in the concentration step, and reformat the user guide. This version of the user guide (Rev. A.0) replaces the Rev. 2.0. |

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Parafilm™ is the trademark of Bemis Company, Inc.