

SiteClick™ Biotin Antibody Labeling Kit

Cat. No. S20033

Pub. No. MAN0016807

Rev. A

For detailed protocol, visit the product page at www.thermofisher.com.

Storage

Store at 2–8°C. **Do not freeze.** When stored as directed, the product is stable for at least 6 months.

Overview

The SiteClick™ Biotin Antibody Labeling Kit allows you to conjugate your own antibodies to sDIBO-modified biotin in a copper-free click reaction.

Materials needed, but not provided

- 100–250 µg of antibody, preferably at a concentration of 2–20 mg/mL in a Tris-based buffer, free of carrier proteins and/or azide
- PBS or TBS
- Distilled water

For Research Use Only. Not for use in diagnostic procedures.

ThermoFisher
SCIENTIFIC

Day 1 (Hands-on time: ~1.5 hours, then overnight incubation)

1. *Optional*: Concentrate the antibody and exchange buffer.
(Time required: 1 hour)
2. Modify the carbohydrate domain of the antibody.
(Time required: 4 hours, hands-off)
3. Attach azide moiety to the carbohydrate-modified antibody.
(Time required: 5 minutes hands-on, then overnight incubation)

Day 2 (Hands-on time: ~1.5 hours, then overnight incubation)

4. Purify and concentrate the azide-modified antibody.
(Time required: 1 hour)
5. Conjugate the azide-modified antibody with sDIBO-modified biotin.
(Time required: 5 minutes hands-on, then overnight incubation)

Day 3 (Hands-on time: ~1 hour)

6. *Optional*: Purify and concentrate the biotin-antibody conjugate.
(Time required: 1 hour)

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

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES 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.

Corporate entity: Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

© 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit thermofisher.com/support
thermofisher.com

ThermoFisher
SCIENTIFIC