

# Protein Labeling Kits

Pub. No. MAN0009630

Pub. Part. No. 100021421

Revision A.0

Detailed protocol is available at [www.thermofisher.com](http://www.thermofisher.com).

**WARNING!** For every chemical, 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 [www.thermofisher.com/support](http://www.thermofisher.com/support).

Material	Amount	Storage*
Reactive dye (solid) + Stir bar (Component A)	3 vials	<ul style="list-style-type: none"> <li>• 2–8°C</li> <li>• Protect from light</li> </ul>
Sodium bicarbonate (solid) (Component B)	84mg	
Purification resin (slurry) (Component C)	~25 mL	
10X Elution buffer (liquid) (Component D)	~25 mL	
Purification columns, column funnels, foam holders, collection tubes	3 each	
*The kit can be stored under the conditions listed. For optimal storage conditions of individual components, refer to the labels on the vials or bags. <b>Do not freeze Components C and D.</b>		

## Prepare protein

**Protein purity:** For optimal labeling efficiency, the protein must be free of ammonium ions, imidazole, primary amines (e.g. Tris, glycine), and other proteins/peptides (e.g. BSA, gelatin, serum). Purify the protein as needed.

**Protein concentration/volume:** For efficient labeling, protein concentration should be at 2 mg/mL in 0.5 mL of 1X PBS (pH 7.2–8.0) or 0.1 M sodium bicarbonate buffer (pH 8.0–8.3).

## Label protein

- 1. Prepare 1 M solution of sodium bicarbonate buffer:** Add 1 mL of distilled water to the entire vial of Component B (84 mg sodium bicarbonate). Mix to dissolve.
- 2. Dilute protein:** If the protein concentration is >2 mg/mL, dilute the protein to 2 mg/mL (e.g. 1 mg of protein in 0.5 mL of 1X PBS or 0.1 M sodium bicarbonate buffer).
- 3. Adjust pH:** Add 50 µL of 1 M sodium bicarbonate buffer (from step 1) to 0.5 mL of protein in 1X PBS; final pH should be ~7.5–8.3. Skip this step if the protein is already in sodium bicarbonate buffer (pH 8.0–8.3) or PBS (pH 8.0).
- 4. Prepare labeling reaction mixture:** Add the protein solution (from step 3) to 1 vial of reactive dye (Component A). Cap the vial and mix gently by inversion or by pipetting up and down. Stir the reaction mixture. Do not froth the solution.
- 5. Incubate labeling reaction:** Stir the reaction mixture at slow speed on a magnetic stirrer for 1 hour at room temperature. Do not froth the solution.

**Note:** Because the preparation of the purification column takes ~15 minutes, begin pouring the column (see **Purify labeled protein**, page 2) during the labeling reaction.

## Purify labeled protein

- 6. Prepare purification column/resin:** Assemble the purification column as shown in Figure 1. Add purification resin (Component C) up to ~3 cm from top. Check buffer flow-through with 1X PBS. If buffer flow stalls, remove the resin, clear the frit, then repack the column with resin/buffer. Allow the excess buffer to drain before loading the labeled protein.
- 7. Prepare 1X elution buffer:** Dilute 10X elution buffer (Component D) 10-fold in distilled water at room temperature. Typically, less than 10 mL is required for each purification.
- 8. Purify protein:** Remove the funnel. Load the reaction mixture (from step 5) onto the top of the resin in the column; allow the mixture to settle into resin. Elute the protein with 1X elution buffer (from step 7) or 1X PBS. Collect and retain all fractions.
- 9. Calculate DOL (degree of labeling) and store labeled protein:** Refer to the detailed protocol available at [www.thermofisher.com](http://www.thermofisher.com).

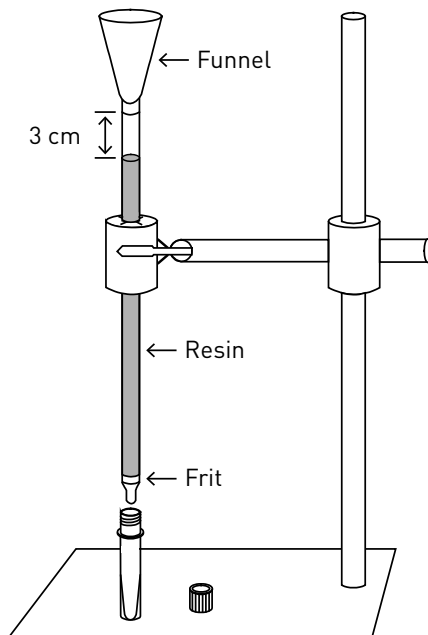


Figure 1

### Documentation and Support

For additional product and technical information, such as Safety Data Sheets (SDS), Certificates of Analysis, etc., visit our website at [www.thermofisher.com](http://www.thermofisher.com). For further assistance, email our Technical Support team at [techsupport@lifetech.com](mailto:techsupport@lifetech.com).

### 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.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 [thermofisher.com/support](http://thermofisher.com/support).

Information in this document 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](http://thermofisher.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)  
[thermofisher.com](http://thermofisher.com)

**ThermoFisher**  
SCIENTIFIC