



AnaTag™ 5-ROX Protein Labeling Kit

Catalog #	71005
Unit Size	1 Kit
Kit Size	5 Conjugation Reactions

This kit is optimized to conjugate 5-ROX (5-Carboxy-X-rhodamine) to proteins (e.g., IgG). It provides ample materials to perform five protein conjugations and purifications. One conjugation reaction can label up to 5 mg protein. The entire process only takes about half an hour.

- **Convenient Format:** Complete kit includes all the assay components.
- **Optimized Performance:** Optimal conditions for conjugation and purification.
- **Enhanced Value:** Less expensive than the sum of individual components.
- **High Speed:** Minimal hands-on time.
- **Assured Reliability:** Detailed protocol and references are provided.

USA and Canada Ordering Information

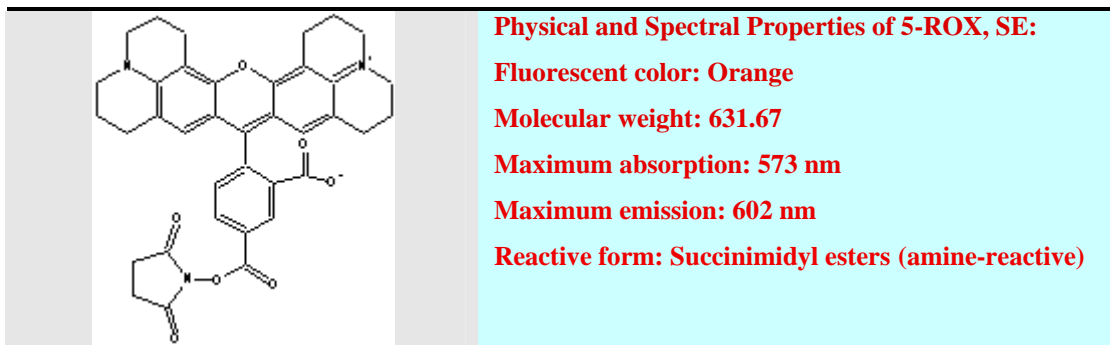
AnaSpec Corporate Headquarter

2149 O'Toole Ave.
San Jose, CA 95131
Toll-Free: 800-452-5530
Tel: 408-452-5055
Fax: 408-452-5059
E-mail: service@anaspec.com
Internet: www.anaspec.com

International Ordering Information

A list of international distributors is available at www.anaspec.com.

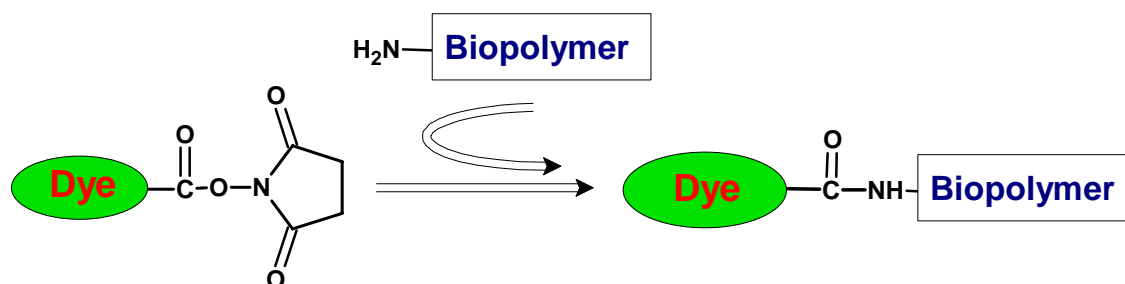
Introduction



The AnaTag™ 5-ROX Protein Labeling Kit provides a convenient way to label proteins with 5-ROX (5-Carboxy-X-rhodamine). 5-ROX is the purified single isomer of 5-(and 6)-ROX that provides reproducibility for critical biological applications and avoids the inconsistency brought by minor positional difference between 5-ROX and 6-ROX.

The succinimidyl ester (SE) reactive form of 5-ROX reacts with the amine group on the protein and forms stable carboxamide bonds, which is identical to, and is as stable as the natural peptide bond (Scheme 1). 5-ROX-protein conjugates are very stable and suitable for immunofluorescent staining, fluorescence *in situ* hybridization, flow cytometry and other biological applications.

5-ROX SE is generally stable if it is properly stored, and shows good reactivity and selectivity towards aliphatic amines. It has a very low reactivity towards aromatic amines, alcohols, phenols (including tyrosine) and histidine. The kit has all the essential components for performing the conjugation reaction and for purifying the conjugate.



KIT COMPONENTS, STORAGE AND HANDLING

Note: Store component A at -20 °C with desiccant. Store all the rest kit components at room temperature.

Component	Function	Quantity
A. 5-ROX, SE	Amino-reactive dye	5 vials
B. Reaction buffer	For pH adjustment of the conjugate reaction	0.5 mL
C. Desalting column	Purify dye-protein conjugate	5 Pre-packed columns
D. DMSO	Solvent for preparing dye solution	1 mL
E. 10X Elution buffer	Solution for eluting dye-protein conjugates	30 mL
F. Buffer reservoir	Hold buffer for desalting column	1

Standard Operating Protocol (SOP)

1. Preparing the protein solution

Add reaction buffer (component B) at 1/10 (v/v) ratio to your target protein (e.g. antibody) solution (2-10 mg/mL is the recommended concentration range).

Note 1: The protein can be dissolved in phosphate, carbonate, borate, triethanolamine or MOPS buffer, pH 7.2-7.5, without reducing reagents (e.g. DTT) or protein stabilizers (e.g. BSA). If the protein is dissolved in Tris or glycine buffer, it should be dialyzed against 0.01 M phosphate buffer saline, pH 7.2-7.4 to get rid of free amines. Ammonium salts (such as ammonium sulfate and ammonium acetate) that are widely used for protein precipitation must also be removed before performing the dye conjugations.

Note 2: The conjugation efficiency is poor when the concentration of protein is less than 2 mg/mL.

2. Preparing the dye solution

Add 20 µL of DMSO (component D) to one vial of 5-ROX SE (component A) to get 10 mg/mL dye solution. Completely dissolve all the dye contents by vortexing.

Note: Dye solution must be prepared fresh for each conjugation reaction. Extended storage of the dye solution may reduce dye activity. Any solutions containing the dye should be kept away from light.

3. Performing the conjugation reaction

3.1 Add the dye solution to the IgG solution at the molar ratio 10:1 (dye: IgG). Table 1 gives a quick reference for labeling IgG. Mix the reaction mixture completely.

Table 1. The volume of dye solution needed for different amount of IgG.

Ig G	Dye solution	Ig G	Dye solution
0.5 mg	2.1 µL	3 mg	12.6 µL
1 mg	4.2 µL	3.5 mg	14.7 µL
1.5 mg	6.3 µL	4 mg	16.8 µL
2 mg	8.4 µL	4.5 mg	19.9 µL
2.5 mg	10.5 µL	5 mg	21.1 µL

Note 1: The molecular weight of IgG is 150 kDa.

Note 2: The dye: protein molar ratio 10:1 is optimized for labeling IgG with 5-ROX SE. For proteins other than IgG, the optimal dye/protein molar ratio needs to be experimentally determined. The desired dye/protein molar ratio usually should be between 2:1 and 20:1.

3.2 Keep the reaction mixture away from light and shake for 15 min at room temperature on a rotator or a shaker.

4. Purify dye-protein conjugates

Note: Desalting column (component C) is best suited for purifying proteins of MW>5,000. For smaller proteins, we recommend using Sephadex LH-20 or dialysis. HPLC may also be used to purify the smaller protein conjugates.

4.1 Dilute 10X elution buffer (component E) to 1X in deionized water.

4.2 Hold the desalting column (component C) upright. Remove the top cap of the column, and then cut its bottom tip. Place the buffer reservoir (component F) on the top of the column.

4.3 Add 25 mL 1X elution buffer into the buffer reservoir to pre-equilibrate the column.

4.4 As soon as the liquid runs just below the top solid surface, load the column with the reaction mixture (directly from step 3.2.).

4.5 As soon as the reaction mixture runs just below the top gel surface, add 10 mL 1X elution buffer into the column.

4.6 When the reaction mixture runs down the column, you should see it separated into two bands. The faster-running band (lower band) contains the desired dye-labeled protein, while the slower-running band (upper band) contains the free dye.

4.7 Collect the faster-running band only. Avoid the slower-running band, which will contaminate your conjugate.

4.8 The degree of substitution (DOS) of the conjugate can be determined according to [Appendix I](#).

Appendix I. Characterizing the Dye-Protein Conjugate

The Degree of substitution (DOS) is the most important factor for characterizing dye-labeled protein. Proteins of lower DOS usually have weaker fluorescence intensity, but proteins of higher DOS (e.g. DOS>8) tend to have reduced fluorescence. The optimal DOS recommended for most antibodies is between 2 and 8.

1. Read the absorbance of dye-protein conjugate at 280 nm (A_{280}) and A_{max}

For most spectrophotometers, dilute a small portion of conjugate solution in phosphate buffered saline so that the absorbance readings are in the 0.1 to 0.9 range. The maximal absorption of protein is at 280 nm (A_{280}). The maximal absorption of 5-ROX amide (A_{max}) is approximately at 573 nm.

2. Calculating the DOS using the following equations for IgG labeling

Molar concentration of dye:

$$[\text{Dye}] = (A_{max} / \epsilon_{5\text{-ROX}}) \times \text{dilution factor}, \quad \epsilon_{5\text{-ROX}} = 82,000 \text{ cm}^{-1}\text{M}^{-1}$$

Molar concentration of protein:

$$[\text{Protein}] = [(A_{280\text{nm}} - 0.16 \times A_{max}) / \epsilon_{\text{protein}}] \times \text{dilution factor} \quad \epsilon_{\text{IgG}} = 203,000 \text{ cm}^{-1}\text{M}^{-1}$$

$$\text{DOS} = [\text{Dye}] / [\text{Protein}] \quad \text{where } \epsilon \text{ is the extinction coefficient}$$

Protein concentration (mg/mL):

$$\text{IgG (mg/mL)} = [\text{IgG}] \times 150,000 \quad \text{MW}_{\text{IgG}} = 150,000$$

For effective labeling, the degree of substitution should fall within 2-8 moles of 5-ROX to one mole of protein.

The Storage of Dye-protein Conjugates

The dye labeled protein should be stored at > 0.5 mg/mL or in the presence of a carrier protein (e.g., 0.1% Bovine Serum Albumin). We recommend adding preservative (e.g. 0.1% sodium azide) to the solution. The dye labeled protein can be stored at 4°C for two months without significant changes if kept from light. For extended storage, it should be divided into aliquots and stored at -20°C in the dark. Avoid multiple thaw-freeze cycles.

Reference

1. Hermanson GT (1996). *Bioconjugate Techniques*, Academic Press, New York.
2. Haugland RP (1995). Coupling of monoclonal antibodies with fluorophores. *Methods Mol Biol* **45**, 205-21.
3. Brinkley M (1992). A brief survey of methods for preparing protein conjugates with dyes, haptens, and cross-linking reagents. *Bioconjug Chem* **3**, 2-13.