

EZLabel™ Antibody Cy5 Labeling Kit

(Store at 4°C)

6/15

Cat. No. K838-5, contains sufficient reagents to label and purify 5 x 1 mg of antibody

I. Introduction:

BioVision's EZLabel™ Antibody Cy5 Labeling Kit provides an easy way to label antibodies with Cy5 in a user-friendly spin column format. Cy5 is an ideal fluorescent dye for labeling of antibodies with its intense signal in the red region of the spectrum. Each pair of spin columns provided in the kit can be used to purify up to 1 mg of the labeled target protein. The kit provides all of the reagents necessary to perform five labeling reactions using up to 1 mg of antibody per reaction. Cy5-labeled antibody has excitation and emission wavelengths at 650 nm and 670 nm respectively, and can be directly used for multiple downstream applications including ELISA, western blot, Immunohistochemistry, Immunofluorescence, FACS, etc.

II. Applications:

- Cy5 labeled antibodies can be used for ELISA, western blot, Immunohistochemistry, Immunofluorescence, FACS, etc.

III. Kit Contents:

Components	K838-5	Cap Code	Part Number
EZLabel™ Cy5	5 vials	Red	K838-5-1
EZLabel™ Spin Column	10 columns	-	K838-5-2
EZLabel™ Elution Buffer	10 ml	NM	K838-5-3

IV. User Supplied Reagents and Equipment:

- Microcentrifuge, DMSO/DMF, and fresh 0.1 M Sodium Bicarbonate buffer (pH 8.5-9.0).

V. Reagent Preparation and Storage Conditions:

Store the kit at 4°C, protected from light. Read the entire protocol before performing the experiment. Briefly spin small vials prior to opening. Bring the kit components to room temperature before use.

VI. Antibody Cy5 Labeling Protocol:

A. Antibody Solution Preparation: The volume of antibody solution should not exceed 100 µl. For best results, use 100 µl of ~5-10 mg/ml antibody.

Note: Buffers that contain primary amines (e.g. Tris or glycine) interfere with the intended Cy5 conjugation. Dialyze the antibody using **BioVision's DiaEasy™ Dialyzer tubes (Cat #s K1000 to K1022)** against 0.1 M sodium bicarbonate buffer (pH 8.5-9.0) just before labeling experiment is performed to remove the primary amines.

B. Labeling Reaction: Each vial of EZLabel™ Cy5 is sufficient for labeling of 1 mg of antibody. Reconstitute one vial of EZLabel™ Cy5 with 5-10 µl of DMSO or DMF just before use. Dissolve completely by pipetting up and down. Transfer 100 µl of the prepared antibody to a 1.5 ml microcentrifuge tube. Add reconstituted EZLabel™ Cy5 solution and mix well by pipetting up and down. Incubate at room temperature on rotary shaker or mixer for 1 hr. Total volume at this stage should not exceed 110 µl.

Note: If the amount of antibody is less than 1 mg, the amount of EZLabel™ Cy5 also needs to be lowered accordingly to avoid over-labeling of antibody.

C. Purification of Labeled Antibody:

1. During the labeling reaction, snap off the bottom closure of an EZLabel™ Spin Column and place in a fresh microcentrifuge tube. Centrifuge at ~1500 x g for 1 min. to remove the residual storage buffer. Discard the flow through and wash the resin with 110 µl of EZLabel™ Elution Buffer. Close the cap and centrifuge at 1500 x g for 1 min. Discard the flow through. Repeat this washing step for at least a total of three times.
2. Load the labeling reaction mix (max. 110 µl) to the first spin column drop by drop. Centrifuge the column for 2 min. at 1500 x g to collect the eluant.
3. Transfer the eluant onto the second unused spin column drop by drop. Centrifuge the column for 2 min. at 1500 x g to collect the labeled antibody.
4. Optional: Dialyze the labeled antibody in the dark against a desired storage buffer containing 20-30% glycerol and if necessary, add carrier protein (e.g. BSA) after the dialysis. Store the dialyzed antibody in a tube wrapped with aluminum foil at 4°C (for short term) or -20°C (for long term).

D. Calculations (Optional): In some cases, it is advantageous to determine the number of molecules of Cy5 per molecule of antibody (degree of labeling). For that, measure the absorbance of the labeled antibody at 280 nm (A_{280}) and 650 nm (A_{650}). If necessary, dilute the labeled antibody in EZLabel™ Elution Buffer. Calculate the concentration of labeled antibody and degree of labeling using following equations:

$$\text{Concentration of labeled Antibody (M)} = \frac{A_{280} - (A_{650} \times 0.05)}{203000} \times \text{Path Length Correction} \times \text{Dilution Factor}$$

$$\# \text{ of moles Cy5 per mole Antibody} = \frac{A_{650} \times \text{Dilution Factor} \times \text{Path Length Correction}}{250000 \times \text{Antibody Concentration (M)}}$$

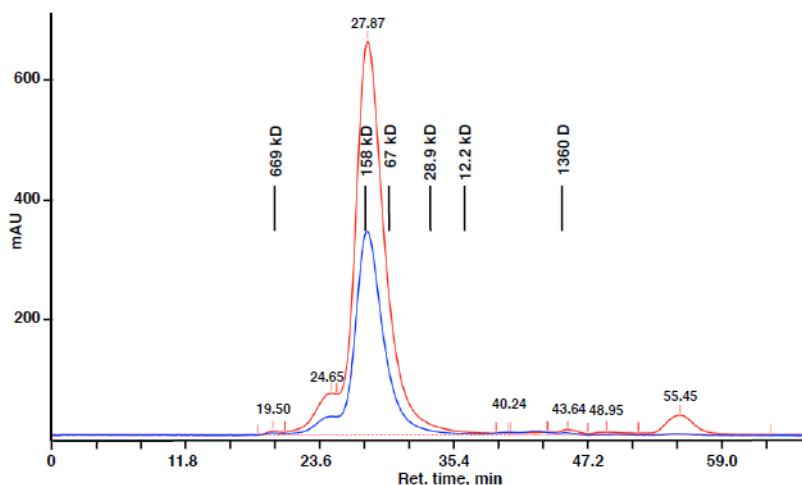


Figure: SEC chromatogram of an Anti-BSA IgG labeled with Cy5 using a Superdex 200 HR 10/30 column at 0.5 ml/min. in 50 mM Tris and 0.25 M NaCl pH 7.5. The absorbance was monitored at 280 nm (Blue line) and 650 nm (Red line). The spin column format ensured that the purification of antibody was fast and there was no unreacted Cy5 left after the antibody was purified according to the kit protocol.

VII. RELATED PRODUCTS:

Annexin V-FITC Apoptosis Kit (K101)	Hi-Bind™ Protein A-Agarose (6520)
Annexin V-FITC Reagent (1001)	Hi-Bind™ Albumin-IgG Depletion Beads (7933)
Human IgG (1296)	Mouse IgG (1265)
Rabbit IgG (1268)	Hi-Bind Ni QR Agarose Beads (6562)
Annexin V-Biotin Apoptosis Kit (K109)	Annexin V-Cy3 Apoptosis Kit (K102)
Annexin V- Cy5 Apoptosis Kit (K103)	Annexin V-EGFP Apoptosis Kit (K104)
Annexin V- PE Apoptosis Detection Kit (K128)	Annexin V-PE-Cy5 Apoptosis Detection Kit (K129)
Red Fluorescent Protein R-PE (R-Phycoerythrin) (6005)	Red Fluorescent Protein Monoclonal Antibody (3984)
EZlabel™ Antibody FITC Labeling Kit (K831)	EZlabel™ Protein FITC Labeling Kit (K832)
EZlabel™ Antibody Biotin Labeling Kit (K834)	EZlabel™ Protein Cy5 Labeling Kit (K839)
EZlabel™ Antibody Cy3 Labeling Kit (K836)	EZlabel™ Protein Cy3 Labeling Kit (K837)
EZ-Desalt™ Spin Desalting Columns (6564)	EZlabel™ Protein Biotin Labeling Kit (K835)

FOR RESEARCH USE ONLY! Not to be used on humans.