

PRODUCT INFORMATION



Protonex™ Red 600, SE

Item No. 44340

FW:	953.1
Purity:	≥90%
Ex./Em. Max:	576/597 nm, in acidic environments
Supplied as:	A solid
Storage:	-20°C
Stability:	≥4 years

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Protonex™ Red 600, SE is supplied as a solid. A stock solution may be made by dissolving the Protonex™ Red 600, SE in the solvent of choice, which should be purged with an inert gas. Protonex™ Red 600, SE is soluble in DMSO.

Description

Protonex™ Red 600, SE is a pH-dependent dye that is intended for use in labeling proteins and antibodies. It can be used in conjunction with additional dyes for multi-labeling purposes. Protonex™ Red 600, SE displays excitation/emission maxima of 576/597 nm, respectively, with high fluorescence intensity in acidic pH environments, such as phagosomes, endosomes, and lysosomes, but only weakly fluoresces in the extracellular environment. After conjugation, it can be used for microscopy and flow cytometry applications. It has been used to label anti-CD146 antibodies for immunodetection of cancer cells.¹

Unless otherwise noted, all unused stock solutions should be divided into single-use aliquots and stored at -20°C after preparation. Avoid repeated freeze-thaw cycles.

Protein stock solution (Solution A)

1. Mix 100 µl of a reaction buffer (e.g., 1 M sodium carbonate solution or 1 M phosphate buffer with pH ~9.0) with 900 µl of the target protein solution at a concentration of >2 mg/ml.

Note 1: The pH of the protein stock solution (Solution A) should be 8.5 ±0.5. If the pH of the protein solution is lower than 8.0, adjust the pH to the range of 8.0-9.0 using 1 M sodium bicarbonate solution or 1 M pH 9.0 phosphate buffer.

Note 2: The protein should be dissolved in 1X PBS, pH 7.2-7.4. If the protein is dissolved in Tris or glycine buffer, it should be dialyzed against 1X PBS to remove ammonium salts, such as ammonium sulfate and ammonium acetate, which are widely used for protein precipitation and may precipitate the protein, and free amines.

Note 3: Antibodies used for conjugation should be purified and free of sodium azide and thimerosal, which may interfere with the conjugation reaction. If sodium azide or thimerosal are present, use dialysis or a spin column to remove them. Conjugation of antibodies stabilized with BSA or gelatin will result in poor labeling.

Note 4: For optimal labeling efficiency, the recommended final protein concentration is 2-10 mg/ml.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/03/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Protonex™ Red 600, SE stock solution (Solution B)

1. Add anhydrous DMSO into the vial of Protonex™ Red 600, SE to make a 10 mM stock solution. Mix well by pipetting or vortex.

Note: Prepare the Protonex™ Red 600, SE stock solution (Solution B) before starting the conjugation. Use promptly. Extended storage of the dye stock solution may reduce the dye activity. Solution B can be stored in the freezer for two weeks when protected from light and moisture. Avoid freeze-thaw cycles.

SAMPLE EXPERIMENTAL PROTOCOL

Important Note

This protocol was developed for conjugating Protonex™ Red 600, SE to goat anti-mouse IgG. Additional optimization may be necessary for your specific protein.

Note: Each protein requires a distinct dye/protein ratio, which also depends on the properties of dyes. Over-labeling of a protein could detrimentally affect its binding affinity, while the protein conjugates of a low dye/protein ratio give reduced sensitivity.

Run Conjugation Reaction

1. Use a 10:1 molar ratio of Solution B (dye)/Solution A (protein) as the starting point: Add 5 µl of the dye stock solution (Solution B, assuming the dye stock solution is 10 mM) into the vial of the protein solution (95 µl of Solution A) with effective shaking. The concentration of the protein is ~0.05 mM assuming the protein concentration is 10 mg/ml and the molecular weight of the protein is ~200 kDa.

Note: We recommend using a 10:1 molar ratio of Solution B (dye)/Solution A (protein). If it is too low or too high, determine the optimal dye/protein ratio at 5:1, 15:1, and 20:1, respectively.

2. Continue to rotate or shake the reaction mixture at room temperature for 30-60 minutes.

Purify the Conjugate

The following protocol is an example of dye-protein conjugate purification using a Sephadex G-25 column.

1. Prepare the Sephadex G-25 column according to the manufacturer's instructions.
2. Load the reaction mixture (from "Run conjugation reaction") to the top of the Sephadex G-25 column.
3. Add PBS (pH 7.2-7.4) as soon as the sample runs just below the top resin surface.
4. Add more PBS (pH 7.2-7.4) to the desired sample to complete the column purification. Combine the fractions that contain the desired dye-protein conjugate.

Note 1: For immediate use, the dye-protein conjugate should be diluted with staining buffer and aliquoted for multiple uses.

Note 2: For longer-term storage, the dye-protein conjugate solution should be concentrated or freeze-dried.

Reference

1. Nollet, M., Stalin, J., Moyon, A., et al. A novel anti-CD146 antibody specifically targets cancer cells by internalizing the molecule. *Oncotarget* **8(68)**, 112283-112296 (2017).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM