

FITC and TRITC

46424 46425 46112

2081.1

Number
Description
46424
FITC, 1g
46425
FITC, 100mg

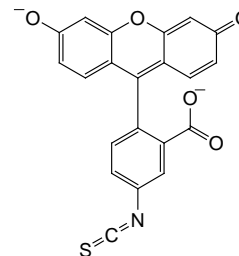
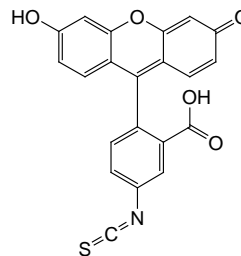
 Chemical name: Fluorescein
 isothiocyanate

Molecular weight: 389.38

 Extinction coefficient: $70,000\text{M}^{-1}\text{cm}^{-1}$
 (in aqueous buffer, pH 9)

Ex/Em wavelength: 494/518nm

CAS # 3326-32-7


46112
TRITC, 10mg

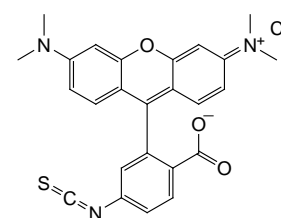
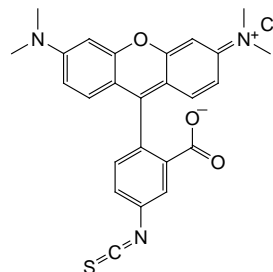
 Chemical name: Tetramethylrhodamine-5-
 (and 6)-isothiocyanate

Molecular weight: 478.97

 Extinction coefficient: $100,000\text{M}^{-1}\text{cm}^{-1}$
 (at 544 nm in methanol)

Ex/Em wavelength: 541/572nm

CAS #6749-36-6



Storage: Upon receipt store at -20°C . Store product in the foil pouch with desiccant to protect from light and moisture. Product is shipped at ambient temperature.

Introduction

Thermo Scientific™ FITC and TRITC are among the most simple and commonly used reagents for protein fluorescent labeling. These isothiocyanates react to amino, sulfhydryl, imidazolyl, tyrosyl or carbonyl groups on proteins; however, only the derivatives of primary and secondary amines generally yield stable products. Reactions are most efficient at pH 8-9 and must be performed in an amine-free buffer such as carbonate/bicarbonate. Antibody and other proteins can be effectively labeled with several fluorophore tags per protein molecule when reacted with a 15- to 20-fold molar excess of isothiocyanate-activated fluorophore. Excess nonreacted and hydrolyzed reagent can be removed by dialysis, gel filtration or dye removal resin.

Procedure for Antibody Labeling with FITC

Materials Required

- Conjugation buffer: 50mM borate buffer, pH 8.5 (Product No. 28384)
- Antibody solution: dissolve ~1mg of antibody in 0.5mL of Conjugation Buffer
- Device to remove excess dye, such as a Dye Removal Columns (Product No. 22858), Thermo Scientific™ Zeba™ Spin Desalting Column (Product No. 89891) or Slide-A-Lyzer™ Dialysis Cassette, 10K MWCO (Product No. 66380)

Procedure

1. Dissolve FITC in DMF at 10mg/mL. Mix well to completely dissolve the FITC.
2. Add 15- to 20-fold molar excess of FITC to 0.5mL of antibody solution and immediately mix the reaction.
3. Incubate for 1 hour at room temperature in the dark.
4. Remove excess and hydrolyzed FITC by gel filtration, dialysis or with a Dye Removal Column.

Procedure for Antibody Labeling with TRITC

This method is adapted from Larsson.¹

Materials Required

- Conjugation Buffer: 100mM carbonate/bicarbonate buffer, pH 9.0
- Antibody solution: dialyze antibody into Conjugation Buffer at 6 mg/ml
- Device to remove excess dye, such as a Dye Removal Column (Product No. 22858), Zeba Spin Desalting Column (Product No. 89891) or Slide-A-Lyzer Dialysis Cassette, 10K MWCO (Product No. 66380)

Procedure

1. Dissolve TRITC in DMSO at 1mg/mL.
2. While stirring, slowly add 35µl of TRITC to 1mL of the 6mg/mL antibody solution and mix thoroughly.
2. Incubate for 2 hours at room temperature in the dark.
3. Remove excess and hydrolyzed TRITC by gel filtration, dialysis or with a Dye Removal Column.

Additional Information

- Visit our website for a listing of related products (other fluorophores and convenient labeling kits) and technical resources, such as the following Tech Tips:
 - Tech Tip #43: Protein stability and storage
 - Tech Tip #6: Extinction coefficients guide
 - Tech Tip #31: Calculate dye:protein (F/P) molar ratios
- Fading (photobleaching in tissue sections) sometimes can be reduced by mounting in an alkaline-buffered media (pH 9).¹ There are several reagents that may be used with FITC and/or TRITC derivatives to prevent fading, including n-propyl gallate at 0.1-0.25M dissolved in glycerol.² For FITC derivatives, *o*- or *p*-phenylenediamine added to the mounting buffer from 1µg/mL to 1mg/mL in glycerin also may be used.^{1,2}

Related Thermo Scientific Products

28384	BupH™ Borate Buffer Packs, 40 packs
28372	BupH Phosphate Buffered Saline Packs, 40 packs
22858	Fluorescent Dye Removal Columns
66380	Slide-A-Lyzer Dialysis Cassettes, 10K MWCO
89889	Zeba Spin Desalting Columns, 7K MWCO, 2mL, 5 each
89890	Zeba Spin Desalting Columns, 7K MWCO, 2mL, 25 each
53029	Pierce™ NHS-Fluorescein Antibody Labeling Kit
53031	Pierce NHS-Rhodamine Antibody Labeling Kit
53027	Pierce FITC Antibody Labeling Kit
46409	NHS-Fluorescein, 1g
46410	NHS-Fluorescein, 100mg
46406	NHS-Rhodamine, 25mg
53024	DyLight™ 488 Antibody Labeling Kit
53025	DyLight 488 Microscale Antibody Labeling Kit

Cited References

1. Larsson, L. (1988). *Immunocytochemistry: Theory and Practice*. CRC. Boca Raton, **77-83**:224-5.
2. Goding, J. (1986). *Monoclonal Antibodies: Principles and Practice*, 2nd ed. Academic, London.

General Reference

1. Horisberger, M. (1984). In *Immunolabeling for Electron Microscopy*. Polak, J., Varndel, I. Ed. Elsevier: Amsterdam

Products are warranted to operate or perform substantially in conformance with published Product specifications in effect at the time of sale, as set forth in the Product documentation, specifications and/or accompanying package inserts ("Documentation"). No claim of suitability for use in applications regulated by FDA is made. The warranty provided herein is valid only when used by properly trained individuals. Unless otherwise stated in the Documentation, this warranty is limited to one year from date of shipment when the Product is subjected to normal, proper and intended usage. This warranty does not extend to anyone other than Buyer. Any model or sample furnished to Buyer is merely illustrative of the general type and quality of goods and does not represent that any Product will conform to such model or sample.

NO OTHER WARRANTIES, EXPRESS OR IMPLIED, ARE GRANTED, INCLUDING WITHOUT LIMITATION, IMPLIED WARRANTIES OF MERCHANTABILITY, FITNESS FOR ANY PARTICULAR PURPOSE, OR NON INFRINGEMENT. BUYER'S EXCLUSIVE REMEDY FOR NON-CONFORMING PRODUCTS DURING THE WARRANTY PERIOD IS LIMITED TO REPAIR, REPLACEMENT OF OR REFUND FOR THE NON-CONFORMING PRODUCT(S) AT SELLER'S SOLE OPTION. THERE IS NO OBLIGATION TO REPAIR, REPLACE OR REFUND FOR PRODUCTS AS THE RESULT OF (I) ACCIDENT, DISASTER OR EVENT OF FORCE MAJEURE, (II) MISUSE, FAULT OR NEGLIGENCE OF OR BY BUYER, (III) USE OF THE PRODUCTS IN A MANNER FOR WHICH THEY WERE NOT DESIGNED, OR (IV) IMPROPER STORAGE AND HANDLING OF THE PRODUCTS.

Unless otherwise expressly stated on the Product or in the documentation accompanying the Product, the Product is intended for research only and is not to be used for any other purpose, including without limitation, unauthorized commercial uses, in vitro diagnostic uses, ex vivo or in vivo therapeutic uses, or any type of consumption by or application to humans or animals.

Current product instructions are available at www.thermoscientific.com/pierce. For a faxed copy, call 800-874-3723 or contact your local distributor.

© 2013 Thermo Fisher Scientific Inc. All rights reserved. Triton is a trademark of The Dow Chemical Company. All (other) trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.