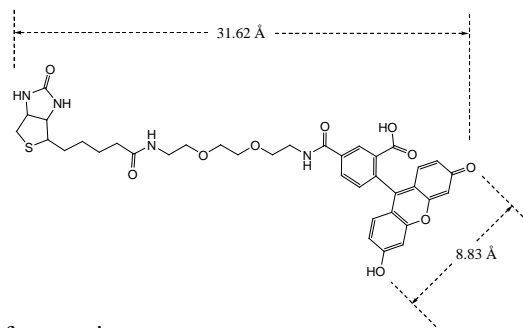


# Biotin-Fluorescein

22030

0745.3

Number	Description
22030	<b>Biotin-Fluorescein</b> [(+)-biotinyl-fluorescein], 5mg
	Formula: C <sub>37</sub> H <sub>40</sub> N <sub>4</sub> O <sub>10</sub> S
	Molecular Weight: 732.80
	Maximum Solubility: > 20mg/mL in DMF or DMSO
	Ex/Em Wavelength: 491/518nm
	Extinction Coefficient: 76,600M <sup>-1</sup> cm <sup>-1</sup>



**Storage:** Upon receipt store product at 4°C, protected from moisture.  
 Product is shipped at room temperature

## General Information

Thermo Scientific Biotin-Fluorescein is a heterobifunctional reagent containing a biotin moiety and a fluorescein group separated by a long-chain hydrophilic polyethylene glycol (PEG) spacer arm. Biotin-Fluorescein is used to label biotin-binding protein molecules including avidin, Thermo Scientific NeutrAvidin Protein and streptavidin, which each have four biotin-binding sites. The resulting fluorescent biotin-binding protein can then amplify a signal generated in assays that use avidin-biotin detection systems.

**Note:** Biotin-Fluorescein is light-sensitive. Avoid exposure to direct sunlight and prolonged exposure to ambient light. Fluorescein does not fluoresce at pH ≤ 6; therefore, use buffers at pH > 6.

## Related Thermo Scientific Products

21120	<b>Hydrazide-Activated Streptavidin, 2mg</b>
21125	<b>Streptavidin, 5mg</b>
31007	<b>Maleimide Activated NeutrAvidin<sup>®</sup> Protein, 5mg</b>
31000	<b>NeutrAvidin Protein, 10mg</b>
21121	<b>Avidin, 10mg</b>
31242	<b>Mouse Anti-Fluorescein Antibody, 1mg</b>
31510	<b>Mouse Anti-Fluorescein, Texas Red-conjugated Antibody, 0.5mg</b>
31687	<b>Mouse Anti-Fluorescein, Rhodamine-conjugated Antibody, 0.5mg</b>

This product ("Product") is 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") and to be free from defects in material and workmanship. Unless otherwise expressly authorized in writing, Products are supplied for research use only. 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 the original purchaser of the Product ("Buyer").

**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 replacement of or refund for the non-conforming Product(s).**

There is no obligation to replace 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.

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

© 2012 Thermo Fisher Scientific Inc. All rights reserved. Unless otherwise indicated, all trademarks are property of Thermo Fisher Scientific Inc. and its subsidiaries. Printed in the USA.