

PRODUCT INFORMATION



Fluorescein Biotin (triethylamine salt)

Item No. 25574

Formal Name: (3aS,4S,6aR)-N-[6-[[[5-[[[(3',6'-dihydroxy-3-oxospiro[isobenzofuran-1(3H),9'-[9H]xanthen]-5-yl)amino]thioxomethyl]amino]pentyl]amino]-6-oxohexyl]hexahydro-2-oxo-1H-thieno[3,4-d]imidazole-4-pentanamide, ditriethylamine salt

Synonym: FITC Biotin

MF: $C_{42}H_{50}N_6O_8S_2 \cdot 2C_6H_{15}N$

FW: 1,033.4

Purity: $\geq 90\%$

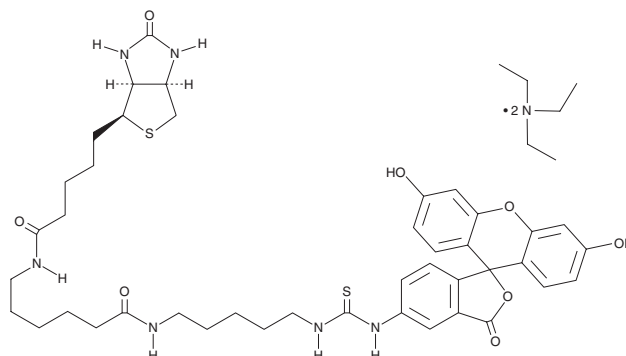
UV/Vis.: λ_{max} : 222 nm

Ex./Em. Max: 485/535 nm

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

Fluorescein biotin (triethylamine salt) is supplied as a solid. A stock solution may be made by dissolving the fluorescein biotin in the solvent of choice, which should be purged with an inert gas. Fluorescein biotin (triethylamine salt) is soluble in methanol, DMSO, and dimethyl formamide.

Fluorescein biotin (triethylamine salt) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Fluorescein biotin (FITC biotin) is a fluorescent biotin derivative and avidin and streptavidin ligand that contains a FITC fluorophore which displays excitation/emission maxima of 485/535 nm, respectively.¹ FITC fluorescence is quenched upon binding to avidin or streptavidin and decreases in fluorescence intensity have been used to quantify biotin binding stoichiometry.

Reference

1. Chu, Y.H., Lees, W.J., Stassinopoulos, A., *et al.* Using affinity capillary electrophoresis to determine binding stoichiometries of protein-ligand interactions. *Biochemistry* **33**(35), 10616-10621 (1994).

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, 05/05/2025

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