

Product Information

Biotin-PEO3-Amine

Catalog Number: 90067

Unit Size: 10 mg

Storage and Handling

Store desiccated -20°C. The product is highly hygroscopic. Product is stable for at least 12 months from date of receipt when stored as recommended.

Stock solution can be prepared in DMSO and stored at 4°C or -20°C for at least 12 months.

Molecular Information: C₁₆H₃₀N₄O₄S

CAS number: 138529-46-1

Molecular Weight: 375

Color and Form: White solid

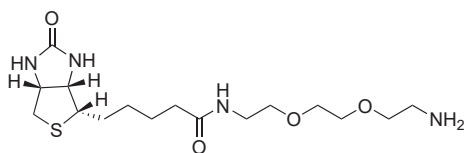


Figure 1. Biotin-PEO3-Amine.

Product Description

Biotin-PEO3-Amine is a water soluble biotin derivative that can be readily coupled to activated carboxylic acids.

Related Products

Cat. No.	Product
92244... 92444	Mix-n-Stain™ Biotin Antibody Labeling Kit
92035... 96010	CF® Dye Amine
90110	Sulforhodamine 101 (Texas Red®)-PEO3-Amine
29031... 29086	Streptavidin Conjugates
90057	Biotin Ethylenediamine, Hydrobromide (Neurobiotin™)
90071	Biotin Ethylenediamine (Free Base)
90075	Biotin Ethylenediamine, Hydrochloride
90079	Biotin Cadaverine, Free Base
90063	Biotin Cadaverine, Trifluoroacetate Salt
90080	Biotin-x Cadaverine, Free Base
90061	Biotin-X Cadaverine, Trifluoroacetate Salt
80018	Lucifer Yellow Cadaverine
80017	Lucifer Yellow Cadaverine Biotin-X, Dipotassium Salt
90108... 91060	ROX Cadaverine, Trifluoroacetate Salt
90109	Sulforhodamine 101 (Texas Red®) Cadaverine
90122	6-TAMRA Cadaverine, Trifluoroacetate Salt
92000	Fluorescein Cadaverine
90082	DMSO, Anhydrous
22014	Bovine Serum Albumin 30% Solution
59002	EDC (EDAC)
22010	Fish Gelatin Blocking Agent (10X)
90083	Biotin-12, SE

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF® Dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Texas Red is a registered trademark of Thermo Fisher Scientific. Neurobiotin is a trademark of Vector Laboratories.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.