

PRODUCT INFORMATION



(+)-Biotin-PEG₃-amine

Item No. 30534

CAS Registry No.: 359860-27-8

Formal Name: (3aS,4S,6aR)-N-[2-[2-[2-(2-aminoethoxy)ethoxy]ethoxy]ethyl]hexahydro-2-oxo-1H-thieno[3,4-d]imidazole-4-pentanamide

Synonym: Amine-PEO₃-biotin

MF: C₁₈H₃₄N₄O₅S

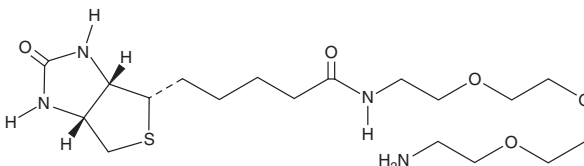
FW: 418.6

Purity: ≥95%

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

(+)-Biotin-PEG₃-amine is supplied as a solid. A stock solution may be made by dissolving the (+)-biotin-PEG₃-amine in the solvent of choice, which should be purged with an inert gas. (+)-Biotin-PEG₃-amine is soluble in organic solvents such as ethanol, DMSO and dimethyl formamide. The solubility of (+)-biotin-PEG₃-amine in these solvents is approximately 20 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of (+)-biotin-PEG₃-amine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of (+)-biotin-PEG₃-amine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(+)-Biotin-PEG₃-amine is a synthetic intermediate.¹ It has been used in the synthesis of biotinylated siderophores for whole-cell bacterial pull-down assays, as well as in the synthesis of protein-polysaccharide conjugate vaccines for *S. pneumoniae*.^{1,2}

References

1. Hu, J., Ghosh, M., Miller, M.J., *et al.* Whole-cell biosensing by siderophore-based molecular recognition and localized surface plasmon resonance. *Anal. Methods* **11**(3), 296-302 (2019).
2. Guo, M., Guo, X., Zhang, C., *et al.* Novel pneumococcal protein-polysaccharide conjugate vaccine based on biotin-streptavidin. *Infect. Immun.* **90**(1), (2022).

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, 08/16/2024

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