

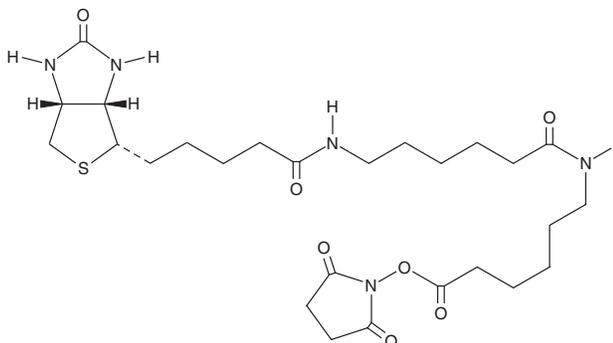
PRODUCT INFORMATION



Biotin-LC-LC-NHS

Item No. 18887

CAS Registry No.: 89889-52-1
Formal Name: (3a*S*,4*S*,6a*R*)-*N*-[6-[[6-[(2,5-dioxo-1-pyrrolidinyl)oxy]-6-oxohexyl]amino]-6-oxohexyl]hexahydro-2-oxo-1*H*-thieno[3,4-*d*]imidazole-4-pentanamide
Synonym: Biotinamidohexanoyl-6-aminohexanoic Acid *N*-hydroxysuccinimide ester
MF: C₂₆H₄₁N₅O₇S
FW: 567.7
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-LC-LC-NHS is supplied as a solid. A stock solution may be made by dissolving the biotin-LC-LC-NHS in the solvent of choice, which should be purged with an inert gas. Biotin-LC-LC-NHS is soluble in the organic solvent DMSO at a concentration of approximately 25 mM.

Description

Biotin-LC-LC-NHS is a biotinylation reagent that contains an *N*-hydroxysuccinimide (NHS) moiety, which activates carboxylic acid groups on biotin to facilitate coupling reactions, and a 30.5 Å spacer to decrease steric hindrance.¹ It has been used to biotinylate small molecules such as paclitaxel (Item No. 10461) and coenzyme Q₁₀ (Item No. 11506) for use in binding site characterization.^{1,2}

References

1. Lis, L.G., Smart, M.A., Luchniak, A., *et al.* Synthesis and biological evaluation of a biotinylated paclitaxel with an extra-long chain spacer arm. *ACS Med. Chem. Lett.* **3(9)**, 745-748 (2012).
2. Murai, M., Yamashita, T., Senoh, M., *et al.* Characterization of the ubiquinone binding site in the alternative NADH-quinone oxidoreductase of *Saccharomyces cerevisiae* by photoaffinity labeling. *Biochemistry* **49(13)**, 2973-2980 (2010).

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, 11/07/2022

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