

PRODUCT INFORMATION



Sulfo-NHS-LC-Biotin (sodium salt)

Item No. 36579

CAS Registry No.: 191671-46-2

Formal Name: 6-[[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]-hexanoic acid, monosodium salt, 2,5-dioxo-3-sulfo-1-pyrrolidinyl ester

Synonym: N-hydroxysulfosuccinimide-LC-Biotin

MF: $C_{20}H_{29}N_4O_9S_2 \cdot Na$

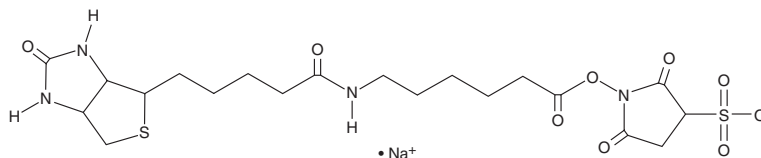
FW: 556.6

Purity: $\geq 90\%$

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

Sulfo-NHS-LC-biotin (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the sulfo-NHS-LC-biotin (sodium salt) in the solvent of choice, which should be purged with an inert gas. Sulfo-NHS-LC-biotin (sodium salt) is soluble in DMSO and methanol.

Description

Sulfo-NHS-LC-biotin is a water soluble and amine-reactive reagent.^{1,2} It contains a spacer arm connecting biotin to the NHS ester, which reduces steric hindrance.² Sulfo-NHS-LC-biotin has commonly been used in the labeling of amine-containing proteins, antibodies, and other molecules.

References

1. Cohen, L. and Walt, D.R. Evaluation of antibody biotinylation approaches for enhanced sensitivity of Single Molecule Array (Simoa) immunoassays. *Bioconjug. Chem.* **29(10)**, 3452-3458 (2018).
2. Elia, G. Biotinylation reagents for the study of cell surface proteins. *Proteomics* **8(19)**, 4012-4024 (2008).

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, 01/29/2025

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