

PRODUCT INFORMATION



Sulfo-NHS-Biotin

Item No. 33362

CAS Registry No.: 194041-65-1

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

Synonym: N-Hydroxysulfosuccinimide-Biotin

MF: $C_{14}H_{18}N_3O_8S_2 \cdot Na$

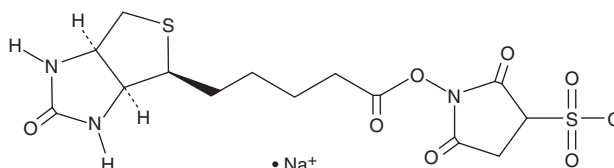
FW: 443.4

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-biotin is supplied as a solid. A stock solution may be made by dissolving the sulfo-NHS-biotin in the solvent of choice, which should be purged with an inert gas. Sulfo-NHS-biotin is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of sulfo-NHS-biotin in these solvents is approximately 20 and 30 mg/ml, respectively.

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 sulfo-NHS-biotin can be prepared by directly dissolving the solid in aqueous buffers. The solubility of sulfo-NHS-biotin in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Sulfo-NHS-biotin is a water-soluble and amine-reactive biotinylation reagent.^{1,2} It is commonly used in the biotin labeling of amine-containing proteins, antibodies, and other molecules.

References

1. Polli, J.W. and Billingsley, M.L. Preparation, characterization and biological properties of biotinylated derivatives of calmodulin. *Biochem J.* **275** (Pt 3), 733-743 (1991).
2. 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).

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, 10/25/2022

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