

PRODUCT INFORMATION



Biotin-X-NHS

Item No. 13316

CAS Registry No.: 72040-63-2
Formal Name: 6-[[5-[(3aS,4S,6aR)-hexhydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]-2,5-dioxo-hexanoic acid, 1-pyrrolidinyl ester

MF: C₂₀H₃₀N₄O₆S

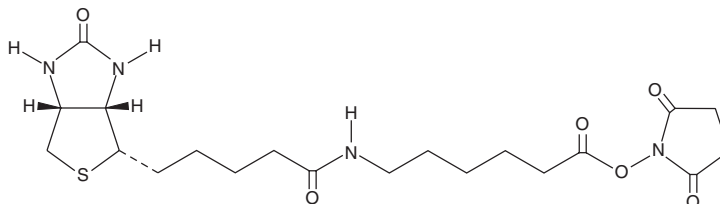
FW: 454.5

Purity: ≥95%

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

Biotin-X-NHS is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, Biotin-X-NHS should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Biotin-X-NHS has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N-Hydroxysuccinimide (NHS) activates carboxylic acid groups (on biotin) to facilitate coupling reactions. Biotin-X-NHS is a compound used to attach biotin to primary amines under alkaline conditions (pH~8-9). For example, lysines on the surface of proteins, including antibodies, are ideal targets for biotinylation with this compound. Biotin-X-NHS is like Biotin-NHS (Item No. 13315), but contains a six-atom spacer arm to minimize steric effects.

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