

PRODUCT INFORMATION



Biotin-XX hydrazide

Item No. 13304

CAS Registry No.: 211237-33-1
Formal Name: 6-[[6-[[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]-1-oxohexyl]amino]-hexanoic acid hydrazide

MF: C₂₂H₄₀N₆O₄S

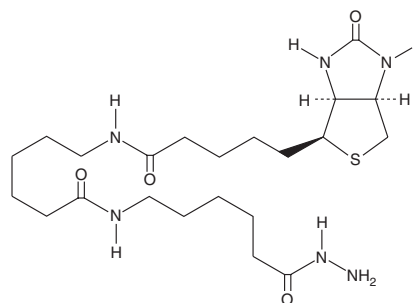
FW: 484.7

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-XX hydrazide is supplied as a crystalline solid. A stock solution may be made by dissolving the biotin-XX hydrazide in the solvent of choice. Biotin-XX hydrazide is soluble in the organic solvent DMSO, which should be purged with an inert gas. The solubility of biotin-XX hydrazide in DMSO is approximately 5 mg/ml.

Biotin-XX hydrazide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, biotin-XX hydrazide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Biotin-XX hydrazide 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

Biotin-XX hydrazide is a biotylating reagent for carbohydrates and carbonyl compounds that contains two aminohexanoic spacers to increase the efficiency of avidin-binding by alleviating steric hindrance.¹

Reference

1. Haugland, R.P., and You, W.W. Coupling of antibodies with biotin. *Methods Mol. Biol.* **418**, 13-23 (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, 12/12/2022

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