

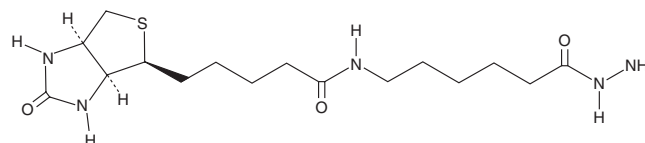
PRODUCT INFORMATION



Biotin LC hydrazide

Item No. 21497

CAS Registry No.: 109276-34-8
Formal Name: 6-[[5-[(3aS,4S,6aR)-hexahydro-2-oxo-1H-thieno[3,4-d]imidazol-4-yl]-1-oxopentyl]amino]-hexanoic acid hydrazide
Synonyms: 6-BACH, Biotinamidocaproyl hydrazide, Biotinamidohexanoic Acid hydrazide
MF: C₁₆H₂₉N₅O₃S
FW: 371.5
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years
Special Conditions: Light and temperature sensitive



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

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

Biotin LC hydrazide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, biotin LC hydrazide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Biotin LC 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 LC hydrazide is a reactive probe that is used to biotinylate carboxyl or vicinal hydroxyl groups on carbohydrates, glycoproteins, or other glycoconjugates.¹ It can also be used to link biotin to antibodies and ribonucleotides.^{1,2} Biotin LC hydrazide contains a long spacer arm to prevent steric hindrances.

References

1. Diamandis, E.P. and Christopoulos, T.K. The biotin-(strept)avidin system: Principles and applications in biotechnology. *Clin. Chem.* **37**(5), 625-636 (1991).
2. Haugland, R.P. and You, W.W. Coupling of antibodies with biotin. *Methods Mol. Bio.* **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, 10/13/2022

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