

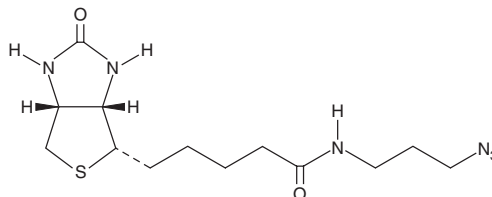
PRODUCT INFORMATION



Biotin-azide

Item No. 13040

CAS Registry No.: 908007-17-0
Formal Name: (3a*S*,4*S*,6a*R*)-*N*-(3-azidopropyl) hexahydro-2-oxo-1*H*-thieno[3,4-*d*]imidazole-4-pentanamide
Synonyms: Azido-biotin, Click Tag™ Biotin-azide
MF: C₁₃H₂₂N₆O₂S
FW: 326.4
Purity: ≥98%
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-azide is supplied as a crystalline solid. A stock solution may be made by dissolving the biotin-azide in the solvent of choice, which should be purged with an inert gas. Biotin-azide is soluble in the organic solvent DMSO, at a concentration of approximately 5 mg/ml.

Biotin-azide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, biotin-azide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Biotin-azide 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-azide is a form of biotin with a terminal azide group. It is used to add biotin to other molecules that bear either an alkyne group, through click chemistry, or a phosphine group, using Staudinger ligation.^{1,2} Biotin-azide has been commonly used to biotin-tag alkynylated lipids, particularly those associated with proteins through post-translational modification.^{1,3,4} It can also be used to biotin-tag proteins, lipids, carbohydrates, and nucleic acids that have been modified with alkyne or phosphine groups.⁵⁻⁸

References

1. Yap, M.C., Kostiuk, M.A., Martin, D.D.O., *et al. J. Lipid Res.* **51**, 1566-1580 (2010).
2. Zhang, J., Li, S., Zhang, D., *et al. Org. Lett.* **12(18)**, 4208-4211 (2010).
3. Kim, H.-Y.H., Tallman, K.A., Liebler, D.C., *et al. Mol. Cell Proteomics* **8(9)**, 2080-2089 (2009).
4. Yount, J.S., Zhang, M.M., and Hang, H.C. *Curr. Protoc. Chem. Biol.* **3(2)**, 65-79 (2011).
5. Ma, Y., Zhang, H., Gruzdys, V., *et al. Langmuir* **27(21)**, 13097-13103 (2011).
6. Crump, C.J., am Ende, C.W., Ballard, T.E., *et al. Bioorg. Med. Chem. Lett.* **22(8)**, 2997-3000 (2012).
7. Sohn, C.H., Agnew, H.D., Lee, J.E., *et al. Anal. Chem.* **84(6)**, 2662-2669 (2012).
8. Al-Shareffi, E., Chaubard, J.-L., Leonhard-Melief, C., *et al. Glycobiology* **23(2)**, 188-198 (2013).

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, 10/14/2022

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