

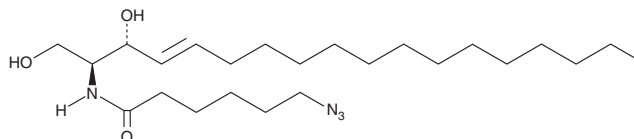
PRODUCT INFORMATION



ω -azido-C6 Ceramide

Item No. 37650

CAS Registry No.: 2108102-18-5
Formal Name: 6-azido-N-[(1S,2R,3E)-2-hydroxy-1-(hydroxymethyl)-3-heptadecen-1-yl]-hexanamide
Synonyms: ω -azido-Cer(18:1/6:0), ω -azido-Ceramide (d18:1/6:0), Click Tag™ ω -azido-C6 Ceramide, ω -azido-N-hexanoyl-D-erythro-Sphingosine, ω -N₃-C6-Cer, ω -N₃-C6 Ceramide
MF: C₂₄H₄₆N₄O₃
FW: 438.7
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

ω -azido-C6 Ceramide is supplied as a crystalline solid. A stock solution may be made by dissolving the ω -azido-C6 ceramide in the solvent of choice, which should be purged with an inert gas. ω -azido-C6 Ceramide is soluble in the organic solvent ethanol at a concentration of approximately 5 mg/ml.

Description

ω -azido-C6 Ceramide is a clickable form of C6 ceramide (Item No. 62525).¹ It is active against *N. meningitidis* and *N. gonorrhoeae* (MIC = 2 μ g/ml for both).² ω -azido-C6 Ceramide has been used to conjugate alkynylated fluorescent probes to liposomes containing ω -azido-C6 ceramide via a click reaction.¹

References

1. Eder, S., Hollmann, C., Mandasari, P., *et al.* Synthesis and characterization of ceramide-containing liposomes as membrane models for different T cell subpopulations. *J. Funct. Biomater.* **13**(3), 111 (2022).
2. Becam, J., Walter, T., Burgert, A., *et al.* Antibacterial activity of ceramide and ceramide analogs against pathogenic *Neisseria*. *Sci. Rep.* **7**(1), 17627 (2017).

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, 11/14/2025

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