

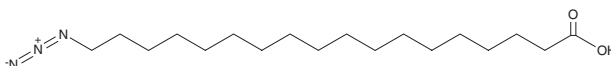
PRODUCT INFORMATION



18-azido Stearic Acid

Item No. 26693

CAS Registry No.: 1529763-58-3
Formal Name: 18-azido-octadecanoic acid
Synonyms: C18:0 Azide, FA 18:0 Azide, Octadecanoic Acid Azide, ω -azido Stearate
MF: $C_{18}H_{35}N_3O_2$
FW: 325.5
Purity: $\geq 95\%$
Supplied as: A 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

18-azido Stearic acid is supplied as a solid. A stock solution may be made by dissolving the 18-azido stearic acid in the solvent of choice, which should be purged with an inert gas. 18-azido Stearic acid is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

18-azido Stearic acid is a clickable form of stearic acid (Item No. 10011298) functionalized with azide, which reacts with alkynes, such as dibenzocyclooctyne (DBCO), to form a stable conjugate.¹ It has been used as a probe to study S-acylation by zinc finger DHHC domain-containing (zDHHC) palmitoyltransferases *in vitro*.² 18-azido Stearic acid has also been used as a building block in the synthesis of a Cy7-labeled lysophosphatidylcholine (LPC) derivative to evaluate brain tumor penetration in an orthotopic GL261 murine glioma model.¹

References

1. Yu, J., Tu, Z., Fan, J., *et al.* Phosphatidylcholine-derived carriers facilitate brain tumor delivery and enhance glioblastoma therapy. *J. Control. Release* **385:113999**, (2025).
2. Greaves, J., Munro, K.R., Davidson, S.C., *et al.* Molecular basis of fatty acid selectivity in the zDHHC family of S-acyltransferases revealed by click chemistry. *Proc. Natl. Acad. Sci. USA* **114(8)**, E1365-E1374 (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, 08/04/2026

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