

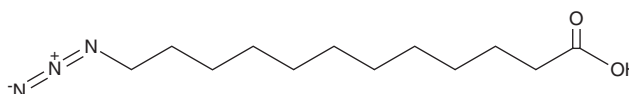
PRODUCT INFORMATION



12-azido Lauric Acid

Item No. 46216

CAS Registry No.: 80667-36-3
Formal Name: 12-azido-dodecanoic acid
Synonyms: C12:0 Azide, FA 12:0 Azide, NSC 667271
MF: $C_{12}H_{23}N_3O_2$
FW: 241.3
Purity: $\geq 95\%$
Supplied as: A low melting 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

12-azido Lauric acid is supplied as a low melting solid. A stock solution may be made by dissolving the 12-azido lauric acid in the solvent of choice, which should be purged with an inert gas. 12-azido Lauric acid is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 12-azido lauric acid can be prepared by directly dissolving the low melting solid in aqueous buffers. 12-azido Lauric acid is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

12-azido Lauric acid is a clickable form of lauric acid (Item Nos. 10006626 | 45444). It has been used as a probe to study protein fatty acylation and in the synthesis of proteolysis-targeting chimeras (PROTACs).^{1,2} 12-azido Lauric acid is active against *C. albicans*, *C. neoformans*, and *S. cerevisiae* (MICs = 1.3, 0.16, and 0.3 mM, respectively) and inhibits the replication of HIV-1 *in vitro*.^{3,4}

References

1. Hang, H.C., Geutjes, E.-J., Grotenberg, G., *et al.* Chemical probes for the rapid detection of fatty-acylated proteins in mammalian cells. *J. Am. Chem. Soc.* **129**(10), 2744-2745 (2007).
2. Cross, J.M., Coulson, M.E., Smalley, J.P., *et al.* A 'click' chemistry approach to novel entinostat (MS-275) based class I histone deacetylase proteolysis targeting chimeras. *RSC Med. Chem.* **13**(12), 1634-1639 (2022).
3. Parang, K., Knaus, E.E., Wiebe, L.I., *et al.* Synthesis and antifungal activities of myristic acid analogs. *Arch. Pharm. (Weinheim)* **329**(11), 475-482 (1996).
4. Devadas, B., Lu, T., Katoh, A., *et al.* Substrate specificity of *Saccharomyces cerevisiae* myristoyl-CoA: Protein N-myristoyltransferase. Analysis of fatty acid analogs containing carbonyl groups, nitrogen heteroatoms, and nitrogen heterocycles in an *in vitro* enzyme assay and subsequent identification of inhibitors of human immunodeficiency virus I replication. *J. Biol.Chem.* **267**(11), 7224-7239 (1992).

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, 09/09/2026

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