

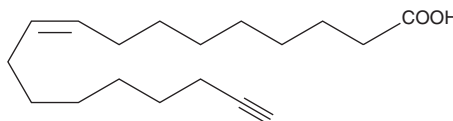
# PRODUCT INFORMATION



## Oleic Acid Alkyne

Item No. 9002078

**CAS Registry No.:** 151333-45-8  
**Formal Name:** (9Z)-9-octadecen-17-ynoic acid  
**Synonyms:** Click Tag™ Oleic Acid Alkyne,  
FA 18:3  
**MF:** C<sub>18</sub>H<sub>30</sub>O<sub>2</sub>  
**FW:** 278.4  
**Purity:** ≥95%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

Oleic acid alkyne is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO, dimethyl formamide (DMF), and 0.1 M Na<sub>2</sub>CO<sub>3</sub> purged with an inert gas can be used. The solubility of oleic acid alkyne in 0.1 M Na<sub>2</sub>CO<sub>3</sub> is approximately 1 mg/ml and approximately 100 mg/ml in DMSO and DMF.

### Description

Oleic acid alkyne is a form of oleic acid (Item No. 90260) with an ω-terminal alkyne. The terminal alkyne group can be used in click chemistry linking reactions, to tag oleic acid with fluorescent or biotinylated labels for analysis of its metabolism and biological activity.<sup>1-4</sup>

### References

1. Kolb, H.C. and Sharpless, K.B. The growing impact of click chemistry on drug discovery. *Drug Discov. Today* **8(24)**, 1128-1137 (2003).
2. Lutz, J.-F. and Zarafshani, Z. Efficient construction of therapeutics, bioconjugates, biomaterials and bioactive surfaces using azide-alkyne “click” chemistry. *Adv. Drug Deliv. Rev.* **60(9)**, 958-970 (2008).
3. Vila, A., Tallman, K.A., Jacobs, A.T., et al. Identification of protein targets of 4-hydroxynonenal using click chemistry for ex vivo biotinylation of azido and alkynyl derivatives. *Chem. Res. Toxicol.* **21(2)**, 432-444 (2008).
4. Jiang, H., Khan, S., Wang, Y., et al. SIRT6 regulates TNF-α secretion through hydrolysis of long-chain fatty acyl lysine. *Nature* **496(7443)**, 110-113 (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, 02/15/2024

#### 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