

# PRODUCT INFORMATION



## Photoclick 3-hydroxy Stearic Acid

Item No. 9004901

**Formal Name:** 13-(3-(but-3-yn-1-yl)-3H-diazirin-3-yl)-3-hydroxytridecanoic acid

**Synonyms:** ClickTag™ 3-hydroxy Stearic Acid, Photoclick 3-HODA, Photoclick 3-Hydroxyoctadecanoic Acid

**MF:** C<sub>18</sub>H<sub>30</sub>N<sub>2</sub>O<sub>3</sub>

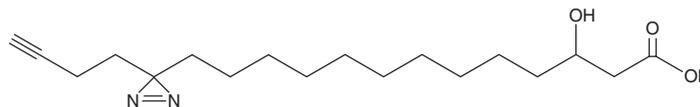
**FW:** 322.4

**Purity:** ≥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

Photoclick 3-hydroxy stearic acid is supplied as a solid. A stock solution may be made by dissolving the photoclick 3-hydroxy stearic acid in the solvent of choice, which should be purged with an inert gas. Photoclick 3-hydroxy stearic acid is soluble in chloroform.

### Description

Photoclick 3-hydroxy stearic acid is a derivative of 3-hydroxy stearic acid (Item No. 10011298) that contains a terminal clickable moiety and a photoactivatable diazirine group at carbon 14. Photoclickable long-chain saturated fatty acids have been used as probes to study protein lipidation.<sup>1</sup>

### Reference

1. Peng, T. and Hang, H.C. Bifunctional fatty acid chemical reporter for analyzing S-palmitoylated membrane protein-protein interactions in mammalian cells. *J. Am. Chem. Soc.* **137**(2), 556-559 (2015).

#### 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, 01/22/2026

#### 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