

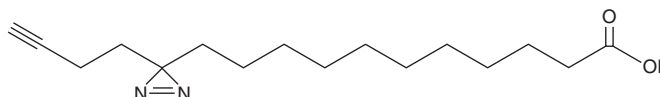
# PRODUCT INFORMATION



## Photoclick Palmitic Acid

Item No. 37969

**CAS Registry No.:** 2988174-40-7  
**Formal Name:** 3-(3-butyn-1-yl)-3H-diazirine-3-undecanoic acid  
**Synonyms:** Click Tag™ Palmitic Acid, 12-Diazirinhexadec-15-yneic Acid, YnPalm Diazirine X12, YnPalmic Diazirine X12



**MF:**  $C_{16}H_{26}N_2O_2$   
**FW:** 278.4  
**Purity:**  $\geq 95\%$   
**Supplied as:** A crystalline solid  
**Storage:**  $-20^{\circ}C$   
**Stability:**  $\geq 4$  years

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

### Laboratory Procedures

Photoclick palmitic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the photoclick palmitic acid in the solvent of choice, which should be purged with an inert gas. Photoclick palmitic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of photoclick palmitic acid in these solvents is approximately 30 mg/ml.

### Description

Photoclick palmitic acid is a derivative of the long-chain saturated fatty acid palmitic acid (Item No. 10006627) that contains a terminal clickable alkyne moiety and a photoactivatable diazirine group at carbon 12. 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, 09/30/2025

#### 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](http://WWW.CAYMANCHEM.COM)