

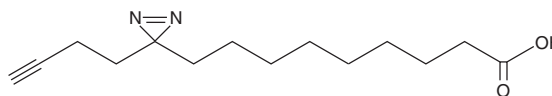
PRODUCT INFORMATION



Photoclick Myristic Acid

Item No. 25363

CAS Registry No.: 2707485-30-9
Formal Name: 3-(3-butyn-1-yl)-3H-diazirine-3-nonanoic acid
Synonyms: Click Tag™ Myristic Acid, YnMyr Diazirine X10, YnMyristic Diazirine X10
MF: C₁₄H₂₂N₂O₂
FW: 250.3
Purity: ≥95%
Supplied as: A crystalline 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 myristic acid is supplied as a crystalline solid. A stock solution may be made by dissolving the photoclick myristic acid in the solvent of choice, which should be purged with an inert gas. Photoclick myristic acid is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of photoclick myristic acid in these solvents is approximately 30 mg/ml.

Photoclick myristic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, photoclick myristic acid should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Photoclick myristic acid has a solubility of approximately 0.04 mg/ml in a 1:20 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Photoclick myristic acid is a derivative of the saturated fatty acid myristic acid (Item No. 13351) that contains a terminal clickable alkyne moiety and a photoactivatable diazirine group at carbon 10. Photoclickable long-chain saturated fatty acids have been used as probes to study protein lipidation.¹

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, 02/16/2024

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