

PRODUCT INFORMATION



Palmitoleic Acid Alkyne

Item No. 25362

CAS Registry No.: 2231023-75-7
Formal Name: (9Z)-9-hexadecen-15-ynoic acid
Synonyms: Click Tag™ Palmitoleic Acid Alkyne,
FA 16:3, 9-cis-Hexadecanoic Acid
Alkyne, n-7 Palmitoleate Alkyne,
Palmitoleate Alkyne,
cis-Palmitoleic Acid Alkyne

MF: C₁₆H₂₆O₂

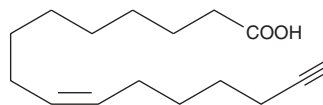
FW: 250.4

Purity: ≥95%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Laboratory Procedures

Palmitoleic 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 and dimethyl formamide purged with an inert gas can be used. The solubility of palmitoleic acid alkyne in these solvents is approximately 100 mg/ml.

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. If an organic solvent-free solution of palmitoleic acid alkyne is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of palmitoleic acid alkyne in PBS, pH 7.2, and 0.15 M Tris-HCl, pH 8.5, is approximately 0.1 and 1 mg/ml, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Palmitoleic acid alkyne is a form of palmitoleic acid (Item No. 10009871) with an ω-terminal alkyne that can be used in click chemistry reactions.¹ It has been used in the study of protein palmitoylation. *cis*-Palmitoleic acid alkyne, but not *trans*-palmitoleic acid alkyne, labels wild-type, but not S209A mutant, Wnt3a in mouse fibroblast L-cells expressing Wnt3a. It also labels secreted wild-type Wnt3a in conditioned media.

Reference

1. Zheng, B., Jarugumilli, G.K., Chen, B., *et al.* Chemical probes to directly profile palmitoylation of proteins. *ChemBioChem* **17**(21), 2022-2027 (2016).

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, 03/06/2024

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