

PRODUCT INFORMATION



13-OxoODE Alkyne

Item No. 9004372

Formal Name: 13-oxooctadeca-9Z,11E-dien-17-ynoic acid
Synonyms: ClickTag™ 13-OxoODE Alkyne, FA 18:3;O Alkyne,
13-KODE Alkyne,
13-keto-9(Z),11(E)-Octadecadienoic Acid Alkyne,
13-keto-9(Z),11(E)-ODE Alkyne,
13-oxo-cis-9-trans-11-Octadecadienoic Acid Alkyne,
13-oxo-cis-9-trans-11-ODE Alkyne

MF: C₁₈H₂₆O₃

FW: 290.4

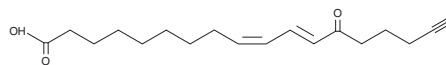
Purity: ≥95%

Supplied as: A solution in ethanol

Storage: -80°C

Stability: ≥2 years

Special Conditions: Oxygen sensitive



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

Laboratory Procedures

13-OxoODE 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 chloroform and dichloromethane purged with an inert gas can be used. 13-OxoODE alkyne is sparingly soluble (1-10 mg/ml) in these solvents.

Description

13-OxoODE alkyne is a clickable form of the oxylipin 13-OxoODE (Item No. 38620) containing an ω-terminal alkyne.

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, 03/27/2024

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