

PRODUCT INFORMATION



12(S),13(S)-DiHODE

Item No. 35864

Formal Name: 12S,13S-dihydroxyoctadeca-6Z,9Z-dienoic acid
Synonyms: 12(S),13(S)-dihydroxy-6(Z),9(Z)-ODE,
12(S),13(S)-dihydroxy-*cis*-6,9-Octadecadienoic Acid,
12(S),13(S)-Dihydroxyoctadecadienoic Acid,
FA 18:2;O2

MF: C₁₈H₃₂O₄

FW: 312.4

Purity: ≥95%

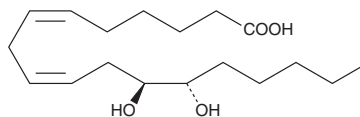
Supplied as: A 100 µg/ml solution in ethanol

Storage: -20°C

Stability: ≥3 years

Special Conditions: Oxygen sensitive

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



Laboratory Procedures

12(S),13(S)-DiHODE 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 purged with an inert gas can be used. 12(S),13(S)-DiHODE is sparingly soluble (1-10 mg/ml) in chloroform.

Description

12(S),13(S)-DiHODE is an oxylipin.

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, 12/01/2025

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