

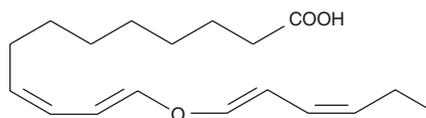
PRODUCT INFORMATION



Etherolenic Acid

Item No. 10005123

CAS Registry No.: 2247927-22-4
Formal Name: 12-(1,3-hexadien-1-yloxy)-9,11-dodecadienoic acid
Synonyms: (9Z,11E,1'E,3'Z)-12-(1',3'-hexadienyloxy)-9,11-Dodecadienoic Acid, FA 18:4;O
MF: C₁₈H₂₈O₃
FW: 292.4
Purity: ≥95% (mixture of isomers)
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.

Description

Etherolenic acid is a divinyl ether oxylipin.^{1,2} It is a metabolite of linolenic acid that is formed in plants via 13-lipoxygenase-mediated formation of 13(S)-HpOTrE (Item No. 45220) followed by conversion to the divinyl ether by divinyl ether synthase.

References

1. Grechkin, A.N., Fazliev, F.N., and Mukhtarova, L.S. The lipoxygenase pathway in garlic (*Allium sativum* L.) bulbs: Detection of the novel divinyl ether oxylipins. *FEBS Letters* **371**, 159-162 (1995).
2. Hamberg, M. Biosynthesis of new divinyl ether oxylipins in Ranunculus plants. *Lipids* **37(4)**, 427-433 (2002).

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, 01/30/2024

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