

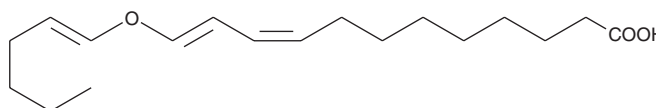
# PRODUCT INFORMATION



## Etheroleic Acid

Item No. 10005119

**CAS Registry No.:** 169217-38-3  
**Formal Name:** 12-[(1E)-1-hexenyloxy]-9Z,11E-dodecadienoic acid  
**Synonym:** FA 18:3;O  
**MF:** C<sub>18</sub>H<sub>30</sub>O<sub>3</sub>  
**FW:** 294.4  
**Purity:** ≥98%  
**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

Etheroleic acid is a divinyl ether oxylipin.<sup>1,2</sup> It is a metabolite of linolenic acid that is formed in plants via 13-lipoxygenase-mediated formation of 13(S)-HpODE (Item No. 48610) or (±)13-HpODE (Item No. 10704) followed by conversion to the divinyl ether by divinyl ether synthase. Etheroleic acid (30 μM) is fungicidal against the powdery mildew *B. graminis* on the second leaves of barley seedlings.<sup>3</sup>

### References

1. Grechkin, A.N., Ilyasov, A.V., and Hamberg, M. On the mechanism of biosynthesis of divinyl ether oxylipins by enzyme from garlic bulbs. *Eur. J. Biochem.* **245(1)**, 137-142 (1997).
2. Hamberg, M. Biosynthesis of new divinyl ether oxylipins in *Ranunculus* plants. *Lipids* **37(4)**, 427-433 (2002).
3. Cowley, T. and Walters, D. Local and systemic effects of oxylipins on powdery mildew infection in barley. *Pest Manag. Sci.* **61(6)**, 572-576 (2005).

#### 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