

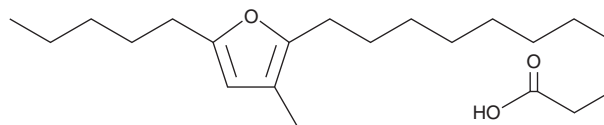
# PRODUCT INFORMATION



## 12,15-epoxy-13-methyl-12,14-Eicosadienoic Acid

Item No. 25862

**CAS Registry No.:** 57818-37-8  
**Formal Name:** 3-methyl-5-pentyl-2-furanundecanoic acid  
**Synonyms:** FA 21:2;O, Furan Fatty Acid F5  
**MF:** C<sub>21</sub>H<sub>36</sub>O<sub>3</sub>  
**FW:** 336.5  
**Purity:** ≥95%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years  
**Special Conditions:** Low melting



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

### Laboratory Procedures

12,15-epoxy-13-methyl-12,14-Eicosadienoic acid is supplied as a low-melting solid. A stock solution may be made by dissolving the 12,15-epoxy-13-methyl-12,14-eicosadienoic acid in the solvent of choice, which should be purged with an inert gas. 12,15-epoxy-13-methyl-12,14-Eicosadienoic acid is slightly soluble in chloroform and methanol.

### Description

12,15-epoxy-13-methyl-12,14-Eicosadienoic acid is a furan fatty acid originally isolated from northern pike (*E. lucius*).<sup>1</sup> Levels of 12,15-epoxy-13-methyl-12,14-eicosadienoic acid are increased in the liver of starving cod.<sup>2</sup>

### References

1. Glass, R.L., Krick, T.P., Sand, D.M., *et al.* Furanoid fatty acids from fish lipids. *Lipids* **10(11)**, 695-702 (1975).
2. Gunstone, F.D., Wijesundera, C., Love, R.M., *et al.* Relative enrichment of furan-containing fatty acids in the liver of starving cod. *J. Chem. Soc. Chem. Commun.* **645(16)**, 630-631 (1976).

#### 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