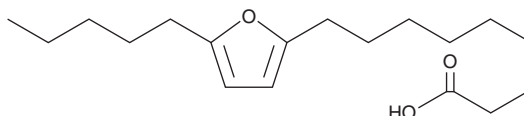


PRODUCT INFORMATION

10,13-epoxy-10,12-Octadecadienoic Acid

Item No. 26767

CAS Registry No.: 4179-43-5
Formal Name: 5-pentyl-2-furannonoic acid
Synonyms: FA 18:2;O, 10,12-furan-FA, 10,12-Furannonoic Acid
MF: C₁₈H₃₀O₃
FW: 294.4
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥4 years



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

Description

10,13-epoxy-10,12-Octadecadienoic acid is a furan fatty acid produced by oxidation of 10(E),12(Z)-conjugated linoleic acid (Item No. 90145).^{1,2} It upregulates the expression of PLIN2, TPD52L2, and SEC13, which are involved in the formation of lipid droplets, and downregulates a variety of genes involved in general cellular processes in Caco-2 cells.² 10,13-epoxy-10,12-Octadecadienoic acid is cytotoxic to Caco-2 and HepG2 cells when used concentrations greater than or equal to 100 μM.

References

1. Yurawecz, M.P., Hood, J.K., Mossoba, M.M., *et al.* Furan fatty acids determined as oxidation products of conjugated octadecadienoic acid. *Lipids* **30**(7), 595-598 (1995).
2. Lengler, I., Buhrke, T., Scharmach, E., *et al.* In-vitro toxicological and proteomic analysis of furan fatty acids which are oxidative metabolites of conjugated linoleic acids. *Lipids* **47**(11), 1085-1097 (2012).

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, 01/16/2026

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