

PRODUCT INFORMATION

(±)9(10)-EpODE

Item No. 40064

CAS Registry No.: 36747-82-7

Formal Name: (2R,3S)-rel-3-(2Z,5Z)-2,5-octadien-1-yl-2-oxiraneoctanoic acid

Synonyms: FA 18:3;O, (±)9(10)-epoxy-12(Z),15(Z)-ODE, (±)9(10)-epoxyoctadecadienoic Acid

MF: C₁₈H₃₀O₃

FW: 294.4

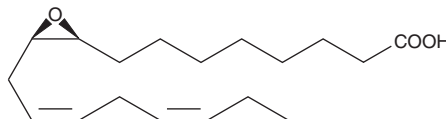
Purity: ≥95%

Supplied as: A 100 µg/ml solution in ethanol

Storage: -20°C

Stability: ≥3 years

Special Conditions: Protect from oxygen



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

Laboratory Procedures

(±)9(10)-EpODE is supplied as a solution in ethanol. To change the solvent, simply evaporate the (±)9(10)-EpODE under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as chloroform and acetonitrile purged with an inert gas can be used. (±)9(10)-EpODE is sparingly soluble (1-10 mg/ml) in chloroform and acetonitrile.

Description

(±)9(10)-EpODE is an oxylipin.¹ It is formed from α-linolenic acid (Item Nos. 90210 | 21910) by the cytochrome P450 (CYP) isoform CYP2B6 in baculosomes containing human CYP2B6.

Reference

- Heintz, M.M., Eccles, J.A., Olack, E.M., *et al.* Human CYP2B6 produces oxylipins from polyunsaturated fatty acids and reduces diet-induced obesity. *PLoS One* **17**(12), e0277053 (2022).

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, 11/18/2025

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