

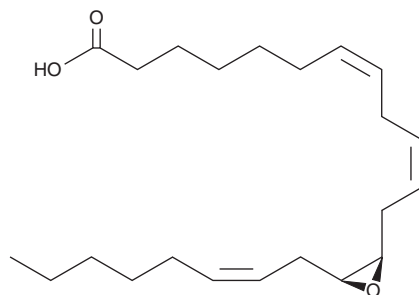
PRODUCT INFORMATION



(±)13,14-EDT

Item No. 38937

CAS Registry No.: 355016-19-2
Formal Name: *rel*-12-[(2R,3S)-3-(2Z)-2-octen-1-yl-2-oxiranyl]-7Z,10Z-dodecadienoic acid
Synonyms: (±)-Dihomo-13,14-EET, (±)13,14-EpDoTrE, (±)13,14-Epoxydocosatrienoic Acid
MF: C₂₂H₃₆O₃
FW: 348.5
Purity: ≥98%
Supplied as: A 100 µg/ml 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

(±)13,14-EDT is an oxylipin and a metabolite of adrenic acid (Item No. 90300) formed *via* the cytochrome P450 (CYP) pathway.^{1,2} It activates large-conductance calcium-activated potassium channels (K_{Ca}1.1/BK) in isolated rat coronary small arterial smooth muscle cells when used at a concentration of 50 nM and induces dilation of isolated porcine arterioles (EC₅₀ = 12 pM).³ (±)13,14-EDT also induces relaxation of isolated bovine coronary arteries precontracted with the TP receptor agonist U-46619 (Item No. 16450).²

References

1. Singh, N., Barnych, B., Wagner, K.M., *et al.* Adrenic acid-derived epoxy fatty acids are naturally occurring lipids and their methyl ester prodrug reduces endoplasmic reticulum stress and inflammatory pain. *ACS Omega* **6**(10), 7165-7174 (2021).
2. Yi, X.-Y., Gauthier, K.M., Cui, L., *et al.* Metabolism of adrenic acid to vasodilatory 1α,1β-dihomo-epoxyeicosatrienoic acids by bovine coronary arteries. *Am. J. Physiol. Heart Circ. Physiol.* **292**(5), H2265-H2274 (2007).
3. Zhang, Y., Oltman, C.L., Lu, T., *et al.* EET homologs potently dilate coronary microvessels and activate BK_{Ca} channels. *Am. J. Physiol. Heart Circ. Physiol.* **280**(6), H2430-H2440 (2001).

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, 07/21/2026

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