

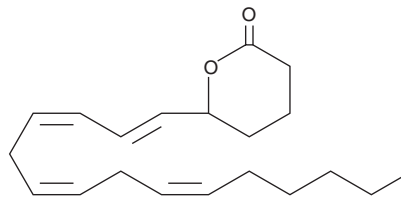
PRODUCT INFORMATION



(±)5-HETE lactone

Item No. 34215

CAS Registry No.: 81517-87-5
Formal Name: tetrahydro-6-(1E,3Z,6Z,9Z)-1,3,6,9-pentadecatetraen-1-yl-2H-pyran-2-one
Synonym: (±)5-Hydroxyeicosatetraenoic Acid lactone
MF: C₂₀H₃₀O₂
FW: 302.5
Purity: ≥98%
UV/Vis.: λ_{max}: 236 nm ε: 27,000
Supplied as: A solution in acetonitrile
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

(±)5-HETE lactone is supplied as a solution in acetonitrile. A stock solution may be made by dissolving the (±)5-HETE lactone in the solvent of choice, which should be purged with an inert gas. (±)5-HETE lactone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of (±)5-HETE lactone in these solvents is approximately 50 mg/ml.

(±)5-HETE lactone is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (±)5-HETE lactone should first be dissolved in acetonitrile and then diluted with the aqueous buffer of choice. The solubility of (±)5-HETE lactone in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(±)5-HETE lactone is a cyclic ester formed by acid-catalyzed nucleophilic addition of the C-5 hydroxyl to the C-1 carboxyl of (±)5-HETE. (±)5-HETE lactone is an inhibitor of 5-lipoxygenase from rat basophilic leukemia cells with an IC₅₀ of 27 μM. The lactone is several times more potent than the linear (±)5-HETE.¹

Reference

1. Kerdesky, F.A.J., Schmidt, S.P., Holms, J.H., *et al.* Synthesis and 5-lipoxygenase inhibitory activity of 5-hydroperoxy-6,8,11,14-eicosatetraenoic acid analogues. *J. Med. Chem.* **30**, 1177-1186 (1987).

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, 11/15/2023

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