

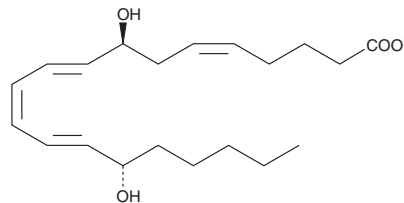
PRODUCT INFORMATION



8(S),15(S)-DiHETE

Item No. 35370

CAS Registry No.: 80234-65-7
Formal Name: 8S,15S-dihydroxy-5Z,9E,11Z,13E-eicosatetraenoic acid
MF: C₂₀H₃₂O₄
FW: 336.5
Purity: ≥95%
UV/Vis.: λ_{max}: 268 nm ε: 40,000
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

8(S),15(S)-DiHETE is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of 8(S),15(S)-DiHETE in these solvents is approximately 50 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of 8(S),15(S)-DiHETE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 8(S),15(S)-DiHETE in PBS (pH 7.2) is approximately 1 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

8(S),15(S)-DiHETE is formed when 15(S)-HETE is subjected to further oxidation by 15-LO.¹ It causes eosinophil chemotaxis with an ED₅₀ value of 1.5 μM but is not chemotactic for neutrophils. 8(S),15(S)-DiHETE antagonizes the hyperalgesic activity of 8(R),15(S)-DiHETE and LTB₄ in the rat hind paw pain model.²

References

1. Morita, E., Schröder, J.M., and Christophers, E. Identification of a novel and highly potent eosinophil chemotactic lipid in human eosinophils treated with arachidonic acid. *J. Immunol.* **144(5)**, 1893-1900 (1990).
2. Levine, J.D., Lam, D., Taiwo, Y.O., et al. Hyperalgesic properties of 15-lipoxygenase products of arachidonic acid. *Proc. Natl. Acad. Sci. USA* **83(14)**, 5331-5334 (1986).

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, 03/29/2024

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