

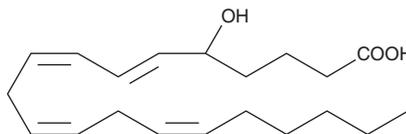
PRODUCT INFORMATION



(±)5-HETE

Item No. 34210

CAS Registry No.: 71030-39-2
Formal Name: (±)5-hydroxy-6E,8Z,11Z,14Z-eicosatetraenoic acid
Synonym: (±)5-Hydroxyeicosatetraenoic Acid
MF: C₂₀H₃₂O₃
FW: 320.5
Purity: ≥95%
UV/Vis.: λ_{max}: 236 nm ε: 27,000
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



Special Conditions: Oxygen and light sensitive

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

Laboratory Procedures

(±)5-HETE is supplied as a solution in ethanol. To change the solvent, evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as dimethyl formamide or DMSO purged with an inert gas can be used. The solubility of (±)5-HETE in these solvents is approximately 50 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made just 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. Organic solvent-free solutions of (±)5-HETE can be prepared using concentrated basic buffers (pH > 8.0 and ionic strength ≥ 0.1 M). Add 400 μl of cold buffer (0°C) per mg of (±)5-HETE and vortex vigorously until completely dissolved. Store aqueous solutions of (±)5-HETE on ice and use within twelve hours.

Description

(±)5-HETE is one of the six monohydroxylated products resulting from the non-enzymatic oxidation of arachidonic acid. It contains an equal mixture of (S)- and (R)-isomers. 5(S)-HETE is the 5-lipoxygenase metabolite of arachidonic acid. (±)5-HETE is a potent neutrophil-aggregating agent, inducing a half-maximal response at 200 nM.¹

Reference

1. O'Flaherty, J.T., Thomas, M.J., Lees, C.J., *et al.* Neutrophil-aggregating activity of monohydroxyeicosatetraenoic acids. *Am. J. Pathol.* **104**, 55-62 (1981).

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