

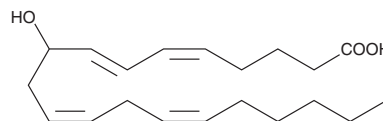
PRODUCT INFORMATION



(±)9-HETE

Item No. 34400

CAS Registry No.: 70968-92-2
Formal Name: (±)-9-hydroxy-5Z,7E,11Z,14Z-eicosatetraenoic acid
Synonym: (±)9-Hydroxyeicosatetraenoic Acid
MF: C₂₀H₃₂O₃
FW: 320.5
Purity: ≥95%
UV/Vis.: λ_{max}: 235 nm
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

(±)9-HETE 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 DMSO and dimethyl formamide purged with an inert gas can be used. (±)9-HETE is miscible in these solvents. The solubility of (±)9-HETE in 0.1 M Na₂CO₃ is approximately 2 mg/ml.

(±)9-HETE is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (±)9-HETE should be diluted with the aqueous buffer of choice. The solubility of (±)9-HETE in PBS (pH 7.2) is approximately 0.8 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

(±)9-HETE is a monohydroxy fatty acid formed by lipid peroxidation of arachidonic acid (Item Nos. 90010 | 90010.1 | 10006607), as well as by rat liver microsomal cytochrome P450 (CYP).^{1,2} Plasma levels of (±)9-HETE are elevated in patients with angiographically defined coronary artery disease.³

References

1. Powell, W.S. and Rokach, J. Biosynthesis, biological effects, and receptors of hydroxyeicosatetraenoic acids (HETEs) and oxoeicosatetraenoic acids (oxo-ETEs) derived from arachidonic acid. *Biochim. Biophys. Acta* **1851**(4), 340-355 (2014).
2. Capdevila, J., Yadagiri, P., Manna, S., *et al.* Absolute configuration of the hydroxyeicosatetraenoic acids (HETEs) formed during catalytic oxygenation of arachidonic acid by microsomal cytochrome P-450. *Biochem. Biophys. Res. Commun.* **141**(3), 1007-1011 (1986).
3. Shishehbor, M.H., Zhang, R., Medina, H., *et al.* Systemic elevations of free radical oxidation products of arachidonic acid are associated with angiographic evidence of coronary artery disease. *Free Rad. Biol. Med.* **41**(11), 1678-1683 (2006).

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, 11/16/2023

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