

PRODUCT INFORMATION

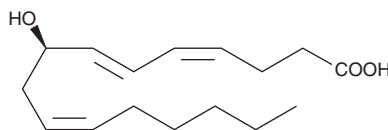


tetranor-12(R)-HETE

Item No. 34565

CAS Registry No.: 135271-51-1
Formal Name: 8R-hydroxy-4Z,6E,10Z-hexadecatrienoic acid
Synonyms: 8(R)-HHxTrE,

2,3,4,5-tetranor-12(R)-HETE,
tetranor-12(R)-Hydroxyeicosatetraenoic Acid
MF: $C_{16}H_{26}O_3$
FW: 266.4
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 234 nm
Supplied as: A 100 $\mu\text{g/ml}$ solution in ethanol
Storage: -80°C
Stability: ≥ 1 year



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

Laboratory Procedures

tetranor-12(R)-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. The solubility of tetranor-12(R)-HETE in these solvents is approximately 25 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 tetranor-12(R)-HETE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of tetranor-12(R)-HETE in PBS (pH 7.2) is approximately 0.50 mg/ml. For greater aqueous solubility, tetranor-12(R)-HETE can be directly dissolved in 0.1 M Na_2CO_3 (2 mg/ml) and then diluted with PBS (pH 7.2) to achieve the desired concentration or pH. We do not recommend storing the aqueous solution for more than one day.

Description

tetranor-12(R)-HETE is a metabolite of 12(R)-HETE (Item No. 34560) formed via β -oxidation.¹

Reference

1. Nishimura, M., Schwartzman, M.L., Falck, J.R., *et al.* Metabolism of 12(R)-hydroxy-5,8,10,14-eicosatetraenoic acid (12(R)-HETE) in corneal tissues: Formation of novel metabolites. *Arch. Biochem. Biophys.* **290**(2), 326-335 (1991).

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, 01/14/2026

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