

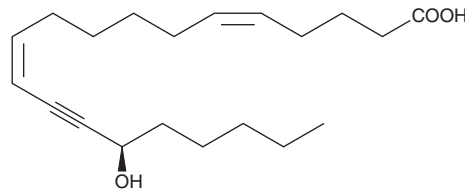
PRODUCT INFORMATION



(5Z,11Z,15R)-15-Hydroxyeicosa-5,11-dien-13-ynoic Acid

Item No. 19557

CAS Registry No.: 339534-01-9
Formal Name: 15R-hydroxy-5Z,11Z-eicosadien-13-ynoic acid
Synonyms: HETE Analog 1,
Hydroxyeicosatetraenoic Acid Analog 1
MF: C₂₀H₃₂O₃
FW: 320.5
Purity: ≥98%
UV/Vis.: λ_{max}: 227 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Laboratory Procedures

(5Z,11Z,15R)-15-Hydroxyeicosa-5,11-dien-13-ynoic acid 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 (DMF) purged with an inert gas can be used. The solubility of (5Z,11Z,15R)-15-hydroxyeicosa-5,11-dien-13-ynoic acid in these solvents is approximately 20 mg/ml.

(5Z,11Z,15R)-15-Hydroxyeicosa-5,11-dien-13-ynoic acid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (5Z,11Z,15R)-15-hydroxyeicosa-5,11-dien-13-ynoic acid should be diluted with the aqueous buffer of choice. (5Z,11Z,15R)-15-Hydroxyeicosa-5,11-dien-13-ynoic acid has a solubility of 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

(5Z,11Z,15R)-15-Hydroxyeicosa-5,11-dien-13-ynoic acid is a stable isomer of 15(S)-HETE (Item No. 34720), a major arachidonic acid metabolite from the 15-lipoxygenase pathway.¹ (5Z,11Z,15R)-15-Hydroxyeicosa-5,11-dien-13-ynoic acid elicits concentration-dependent contraction of isolated pulmonary arteries from rabbits and inhibits the proliferation and migration of hormone-independent prostate carcinoma PC-3 cells.¹

Reference

1. Pfister, S. L., Klimko, P. G., and Conrow, R. E. (5Z,11Z,15R)-15-Hydroxyeicosa-5,11-dien-13-ynoic acid: A stable isomer of 15(S)-HETE that retains key vasoconstrictive and antiproliferative activity. *Prostaglandins Other Lipid Mediat.* **Apr.** (2016).

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, 05/06/2022

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