

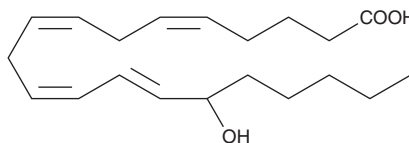
# PRODUCT INFORMATION



## (±)15-HETE

Item No. 34700

**CAS Registry No.:** 71030-36-9  
**Formal Name:** (±)15-hydroxy-5Z,8Z,11Z,13E-eicosatetraenoic acid  
**MF:** C<sub>20</sub>H<sub>32</sub>O<sub>3</sub>  
**FW:** 320.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 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

(±)15-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 (±)15-HETE 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 (±)15-HETE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of (±)15-HETE in PBS (pH 7.2) is approximately 1 mg/ml. Store aqueous solutions of (±)15-HETE on ice and use within 12 hours of preparation.

### Description

(±)15-HETE is one of the six monohydroxy fatty acids produced by the non-enzymatic oxidation of arachidonic acid. The biological activity of (±)15-HETE is similar to that of its constituent enantiomers (Item Nos. 34720 and 34710).

### References

1. Ramis, I., Roselló-Catafau, J., Bulbena, O., *et al.* 15-Hydroxyeicosatetraenoic acid as a major eicosanoid in nasal secretions: Assay by high-performance liquid chromatographic-radioimmunoassay and gas chromatographic-mass spectrometric procedures. *J. Chromatogr.* **496**, 416-422 (1989).
2. Lecomte, M., Laneuville, O., Ji, C., *et al.* Acetylation of human prostaglandin endoperoxide synthase-2 (cyclooxygenase-2) by aspirin. *J. Biol. Chem.* **269**, 13207-13215 (1994).

#### 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, 08/09/2018

#### 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