

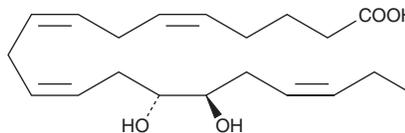
# PRODUCT INFORMATION



## (±)14(15)-DiHETE

Item No. 10006998

<b>Formal Name:</b>	(±)14,15-dihydroxy-5Z,8Z,11Z,17Z-eicosatetraenoic acid
<b>Synonym:</b>	(±)14,15-dihydroxy-eicosa-5,8,11,17-Tetraenoic Acid
<b>MF:</b>	C <sub>20</sub> H <sub>32</sub> O <sub>4</sub>
<b>FW:</b>	336.5
<b>Purity:</b>	≥98%
<b>Supplied as:</b>	A solution in ethanol
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥2 years



NOTE: Relative stereochemistry shown in chemical structure

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

### Laboratory Procedures

(±)14(15)-DiHETE 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 (±)14(15)-DiHETE in these solvents is at least 30 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 (±)14(15)-DiHETE is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of (±)14(15)-DiHETE in PBS (pH 7.2) is at least 0.5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

(±)14(15)-DiHETE is an oxylipin and a metabolite of the ω-3 fatty acid eicosapentaenoic acid (EPA; Item Nos. 90110 | 90110.1 | 21908),<sup>1,2</sup> It is formed via cytochrome P450-mediated epoxidation of EPA to (±)14(15)-EpETE (Item No. 10173), followed by conversion of the epoxide to a diol by soluble epoxide hydrolase.

### References

1. Jónasdóttir, H.S., Ioan-Facsinay, A., Kwekkeboom, J., *et al.* An advanced LC-MS/MS platform for the analysis of specialized pro-resolving lipid mediators. *Chromatographia* **78**(5-6), 391-401 (2014).
2. Shearer, G.C., Harris, W.S., Pedersen, T.L., *et al.* Detection of omega-3 oxylipins in human plasma and response to treatment with omega-3 acid ethyl esters. *J. Lipid. Res.* **51**(8), 2074-2081 (2010).

#### 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, 05/17/2022

#### 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