

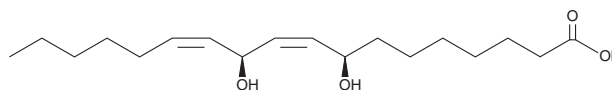
# PRODUCT INFORMATION



## 8(R),11(S)-DiHODE

Item No. 28584

**CAS Registry No.:** 885215-81-6  
**Formal Name:** (8R,9Z,11S,12Z)-8,11-dihydroxy-9,12-octadecadienoic acid  
**Synonym:** 8,11-dihydroxy-9,12-Octadecadienoic Acid  
**MF:** C<sub>18</sub>H<sub>32</sub>O<sub>4</sub>  
**FW:** 312.4  
**Purity:** ≥70%  
**Supplied as:** A solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

8(R),11(S)-DiHODE is supplied as a solid. A stock solution may be made by dissolving the 8(R),11(S)-DiHODE in the solvent of choice, which should be purged with an inert gas. 8(R),11(S)-DiHODE is soluble in ethanol, methanol, and DMSO.

### Description

8(R),11(S)-DiHODE is a fungal oxylipin that has been found in *Aspergillus*.<sup>1,2</sup> It is formed via isomerization of an 8(R)-HpODE intermediate during the oxidation of linoleic acid in several species of *Aspergillus*.

### References

1. Jernerén, F., Garscha, U., Hoffmann, I., *et al.* Reaction mechanism of 5,8-linoleate diol synthase, 10R-dioxygenase, and 8,11-hydroperoxide isomerase of *Aspergillus clavatus*. *Biochim. Biophys. Acta* **1801(4)**, 503-507 (2010).
2. Garscha, U., Jernerén, F., Chung, D., *et al.* Identification of dioxygenases required for *Aspergillus* development. Studies of products, stereochemistry, and the reaction mechanism. *J. Biol. Chem.* **282(48)**, 34707-34718 (2007).

#### 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, 12/02/2022

#### 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