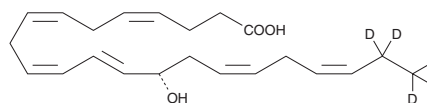


PRODUCT INFORMATION



14(S)-HDHA-d₅ Item No. 9002764

Formal Name:	14S-hydroxy-4Z,7Z,10Z,12E,16Z,19Z-docosa-hexaenoic-21,21,22,22,22-d ₅ acid
Synonyms:	14(S)-HDoHE-d ₅ , 14(S)-hydroxy Docosahexaenoic Acid-d ₅ , FA 22:6-d ₅
MF:	C ₂₂ H ₂₇ D ₅ O ₃
FW:	349.5
Chemical Purity:	≥98% (14(S)-HDHA)
Deuterium Incorporation:	≥99% deuterated forms (d ₁ -d ₅); ≤1% d ₀
Supplied as:	A 100 µg/ml solution in ethanol
Storage:	-20°C
Stability:	≥4 years



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

Laboratory Procedures

14(S)-HDHA-d₅ is intended for use as an internal standard for the quantification of 14(S)-HDHA (Item No. 15253) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Description

14(S)-HDHA is an oxygenation product formed by 12-lipoxygenase (12-LO) or 15-LO processing of docosahexaenoic acid (DHA; Item No. 90310).¹ It is a precursor to the pro-resolving mediator maresin 1 (Item No. 10878) and has been found in peritoneal exudates isolated from a mouse model of zymosan-induced peritonitis.

Reference

1. Serhan, C.N., Yang, R., Martinod, K., *et al.* Maresins: Novel macrophage mediators with potent antiinflammatory and proresolving actions. *J. Exp. Med.* **206**(1), 15-23 (2009).

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, 05/11/2026

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