PRODUCT INFORMATION



(S)-(-)-Docosahexaenyl-1'-Hydroxy-2'-Propylamide

Item No. 9001232

Formal Name:	N-((S)-1-hydroxypropan-2-yl)docosa-	
	4Z,7Z,10Z,13Z,16Z,19Z-hexaenamide	НО
MF:	C ₂₅ H ₃₉ NO ₂	
FW:	385.6	
Purity:	≥98%	
Supplied as:	A solution in ethanol	
Storage:	-20°C	
Stability:	≥2 years	
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.		

Laboratory Procedures

(S)-(-)-Docosahexaenyl-1'-hydroxy-2'-propylamide 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 ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of (S)-(-)-docosahexaenyl-1'-hydroxy-2'-propylamide in ethanol and DMSO is approximately 14 mg/ml and 2 mg/ml in DMF.

(S)-(-)-Docosahexaenyl-1'-hydroxy-2'-propylamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (S)-(-)-docosahexaenyl-1'-hydroxy-2'-propylamide should be diluted with the aqueous buffer of choice. (S)-(-)-Docosahexaenyl-1'-hydroxy-2'-propylamide has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N-Acyl ethanolamines (NAEs) have diverse biological actions that are strongly affected by the associated acyl group. Docosahexaenoyl ethanolamide (DHEA) has potential signaling roles in cancer, inflammation, and neurological development and functioning.¹⁻⁴ At least some of DHEA's effects are mediated through cannabinoid (CB) receptors, while some NAEs also act as vanilloid receptor agonists and voltage-gated K⁺ channel blockers.^{1,4,5} (S)-(-)-Docosahexaenyl-1'-hydroxy-2'-propylamide is a homolog of DHEA, characterized by the addition of an (S)- α -methyl group at the methylene carbon adjacent to the amide nitrogen. A similar modification of arachidonoyl ethanolamide (Item No. 90050) to produce S-1 methanandamide (Item No. 90072) results in a diminished affinity for the CB receptor but greatly improved metabolic stability to aminopeptidase hydrolysis.⁶ The physiological actions of this compound have not been evaluated.

References

- 1. Brown, I., Cascio, M.G., Wahle, K.W.J., et al. Carcinogenesis 31(9), 1584-1591 (2010).
- 2. Kim, H.-Y., Moon, H.-S., Cao, D., et al. Biochem. J. 435, 327-336 (2011).
- 3. Balvers, M.G., Verhoeckx, K.C., Plastina, P., et al. Biochim. Biophys. Acta 1801(10), 1107-1114 (2010).
- 4. Poling, J.S., Rogawski, M.A., Salem, N., Jr., et al. Neuropharmacology 35(7), 983-991 (1996).
- 5. Calignano, A., LaRana, G., and Piomelli, D. Eur. J. Pharmacol. 419(2-3), 191-198 (2001).
- 6. Abadji, V., Lin, S., Taha, G., et al. J. Med. Chem. 37, 1889-1893 (1994).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY 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

uyer 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, 02/20/2024

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