

PRODUCT INFORMATION



α -Linolenoyl Ethanolamide-d₄

Item No. 9001841

CAS Registry No.: 1451194-68-5

Formal Name: N-(2-hydroxyethyl-1,1',2,2'-d₄)-octadeca-9Z,12Z,15Z-trienamide

MF: C₂₀H₃₁D₄NO₂

FW: 325.5

Chemical Purity: ≥98% (α -linolenoyl ethanolamide)

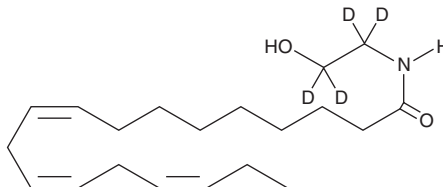
Deuterium

Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀

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

α -Linolenoyl ethanolamide-d₄ is used as an internal standard for the quantification of α -linolenoyl ethanolamide (Item No. 90215) 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).

α -Linolenoyl ethanolamide-d₄ 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 α -linolenoyl ethanolamide-d₄ in these solvents is approximately 30 and 10 mg/ml, respectively.

Description

α -Linolenoyl ethanolamide (Item No. 90215) is an endocannabinoid containing α -linolenic acid in place of the arachidonate moiety of AEA. α -Linolenoyl ethanolamide has been detected in porcine brain, but its specific role and relative importance as a cannabinergic neurotransmitter have not been elucidated.¹

Reference

1. Hanus, L., Gopher, A., Almog, S., *et al.* Two new unsaturated fatty acid ethanolamides in brain that bind to the cannabinoid receptor. *J. Med. Chem.* **36**, 3032-3034 (1993).

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, 02/22/2024

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