

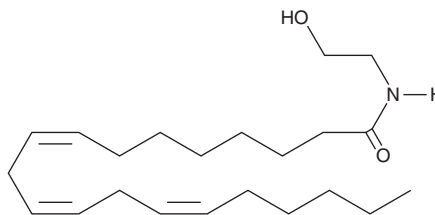
PRODUCT INFORMATION



Dihomo- γ -Linolenoyl Ethanolamide

Item No. 90235

CAS Registry No.: 50314-34-4
Formal Name: N-(2-hydroxyethyl)-8Z,11Z,14Z-
eicosatrienamide
MF: $C_{22}H_{39}NO_2$
FW: 349.6
Purity: $\geq 98\%$
Supplied as: A solution in ethanol
Storage: $-20^{\circ}C$
Stability: ≥ 2 years



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

Laboratory Procedures

Dihomo- γ -linolenoyl ethanolamide 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 or dimethyl formamide purged with an inert gas can be used. The solubility of dihomogamma-linolenoyl ethanolamide is approximately 30 and 10 mg/ml, respectively.

Dihomo- γ -linolenoyl ethanolamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of dihomogamma-linolenoyl ethanolamide should be diluted with the aqueous buffer of choice. Dihomogamma-linolenoyl ethanolamide has a solubility of approximately 8 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

Dihomo- γ -linolenoyl ethanolamide is an endocannabinoid containing dihomogamma-linoleate in place of the arachidonate moiety of AEA. Dihomogamma-linolenoyl ethanolamide binds to recombinant human CB_1 and CB_2 receptor with K_i values of 857 and 598 nM, respectively.¹ Its specific role and relative importance as a cannabinergic neurotransmitter have not been elucidated.²

References

1. Felder, C.C., Joyce, K.E., Briley, E.M., *et al.* Comparison of the pharmacology and signal transduction of the human cannabinoid CB_1 and CB_2 receptors. *Mol. Pharmacol.* **48**, 443-450 (1995).
2. 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, 02/29/2024

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