

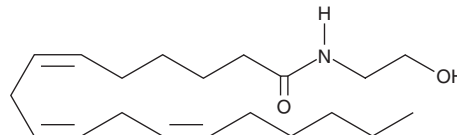
PRODUCT INFORMATION



γ -Linolenoyl Ethanolamide

Item No. 9001747

CAS Registry No.: 150314-37-7
Formal Name: N-(2-hydroxyethyl)-6Z,9Z,12Z-octadecatrienamide
MF: C₂₀H₃₅NO₂
FW: 321.5
Purity: ≥95%
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

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 ethanol, DMSO, and dimethyl formamide (DMF) purged with an inert gas can be used. The solubility of γ -linolenoyl ethanolamide in ethanol is approximately 50 mg/ml and approximately 25 mg/ml in DMSO and DMF.

γ -Linolenoyl ethanolamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of γ -linolenoyl ethanolamide should be diluted with the aqueous buffer of choice. γ -Linolenoyl ethanolamide has a solubility of 0.1 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

γ -Linolenoyl ethanolamide is a member of the family of fatty N-acyl ethanolamines collectively called endocannabinoids.¹⁻⁴ The relative importance of this ethanolamine metabolite has yet to be determined.

References

1. Bachur, N.R. and Udenfriend, S. Microsomal synthesis of fatty acid amides. *J. Biol. Chem.* **241(6)**, 1308-1313 (1966).
2. Doetsch, P.W., Zastawny, T.H., Martin, A.M., *et al.* Monomeric base damage products from adenine, guanine, and thymine induced by exposure of DNA to ultraviolet radiation. *Biochemistry* **34(3)**, 737-742 (1995).
3. Saghatelian, A., Trauger, S.A., Want, E.J., *et al.* Assignment of endogenous substrates to enzymes by global metabolite profiling. *Biochemistry* **43(45)**, 14332-14339 (2004).
4. Buczynski, M.W., Svensson, C.I., Dumlao, D.S., *et al.* Inflammatory hyperalgesia induces essential bioactive lipid production in the spinal cord. *J. Neurochem.* **114(4)**, 981-993 (2010).

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, 02/21/2024

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