

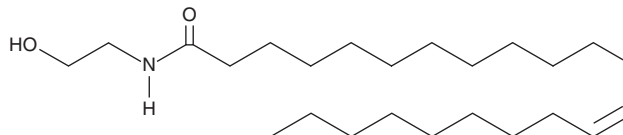
PRODUCT INFORMATION



Docosaenoyl Ethanolamide

Item No. 9001745

CAS Registry No.: 18190-74-4
Formal Name: N-(2-hydroxyethyl)-13Z-docosenamide
MF: $C_{24}H_{47}NO_2$
FW: 381.6
Purity: $\geq 98\%$
Supplied as: A crystalline solid
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

Docosaenoyl ethanolamide is supplied as a crystalline solid. A stock solution may be made by dissolving the docosaenoyl ethanolamide in the solvent of choice, which should be purged with an inert gas. Docosaenoyl ethanolamide is soluble in the organic solvent chloroform at a concentration of approximately 0.3 mg/ml.

Description

Docosaenoyl ethanolamide is a member of the family of fatty N-acyl ethanolamines collectively called endocannabinoids.¹⁻³ This long-chain ethanolamine has been detected at relatively high levels in rat cerebrospinal fluid and has been used to examine the function of voltage-dependent Ca^{2+} channels in skeletal muscle membranes.^{4,5}

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).
5. Oz, M., Alptekin, A., Tchugunova, Y., *et al.* Effects of saturated long-chain N-acyl ethanolamines on voltage-dependent Ca^{2+} fluxes in rabbit T-tubule membranes. *Arch. Biochem. Biophys.* **434**(2), 344-351 (2005).

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, 06/30/2023

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