

PRODUCT INFORMATION



Tricosanoyl Ethanolamide

Item No. 9001743

CAS Registry No.: 171022-15-4

Formal Name: N-(2-hydroxyethyl)-tricosanamide

MF: $C_{25}H_{51}NO_2$

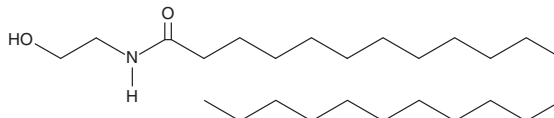
FW: 397.7

Purity: $\geq 98\%$

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: ≥ 4 years



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

Laboratory Procedures

Tricosanoyl ethanolamide is supplied as a crystalline solid. A stock solution may be made by dissolving the tricosanoyl ethanolamide in the solvent of choice, which should be purged with an inert gas. Tricosanoyl ethanolamide is soluble in the organic solvent chloroform at a concentration of approximately 1 mg/ml.

Description

Tricosanoyl 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, 11/14/2025

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