

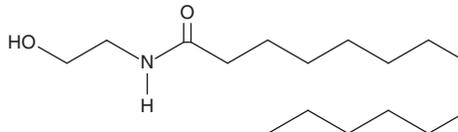
# PRODUCT INFORMATION



## Myristoyl Ethanolamide

Item No. 9001742

**CAS Registry No.:** 142-58-5  
**Formal Name:** N-(2-hydroxyethyl)-  
tetradecanamide  
**Synonyms:** AM3165, Comperlan MM,  
Schercomid MME  
**MF:** C<sub>16</sub>H<sub>33</sub>NO<sub>2</sub>  
**FW:** 271.4  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



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

### Laboratory Procedures

Myristoyl ethanolamide is supplied as a crystalline solid. A stock solution may be made by dissolving the myristoyl ethanolamide in the solvent of choice, which should be purged with an inert gas. Myristoyl ethanolamide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of myristoyl ethanolamide in these solvents is approximately 1, 10, and 20 mg/ml, respectively.

### Description

Myristoyl ethanolamide is a member of the family of fatty N-acyl ethanolamines collectively called endocannabinoids.<sup>1-3</sup> Myristic acid is typically detected at low levels in rat cerebrospinal fluid, however the specific role and relative importance of its ethanolamine metabolite have not been yet determined.<sup>4</sup>

### 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, 04/30/2024

#### 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