

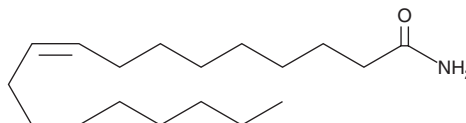
Product Information



9-Octadecenamide

Catalog No. 90375

CAS Registry No.: 301-02-0
Formal Name: 9Z-octadecenamide
Synonym: Oleamide
MF: C₁₈H₃₅NO
FW: 281.5
Purity: ≥98%
Stability: ≥1 year at -20°C
Supplied as: A crystalline solid



Laboratory Procedures

For long term storage, we suggest that 9-octadecenamide be stored as supplied at -20°C. It should be stable for at least one year.

9-Octadecenamide is supplied as a crystalline solid. Stock solutions of 9-octadecenamide can be made by dissolving the crystalline compound in the solvent of choice. Solvents such as ethanol, DMSO, or dimethyl formamide purged with an inert gas can be used. The solubility of 9-octadecenamide in these solvents is approximately 14 mg/ml.

9-Octadecenamide is only sparingly soluble in aqueous buffers. Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Store aqueous solutions of 9-octadecenamide on ice and use within 12 hours of preparation. Although the aqueous solutions of 9-octadecenamide may be stable for more than 12 hours, we strongly recommend using a fresh preparation each day.

9-Octadecenamide is the amide of oleic acid. It was first identified in the cerebrospinal fluid of sleep-deprived cats. It has also been detected in the cerebrospinal fluid of rats and humans. It induces physiological sleep when injected into rats intraperitoneally at 5 to 50 mg doses.¹

Reference

1. Cravatt, B.F., Prospero-Garcia, O., Siuzdak, G., *et al.* Chemical characterization of a family of brain lipids that induce sleep. *Science* **268**, 1506-1509 (1995).

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 01/28/2010