

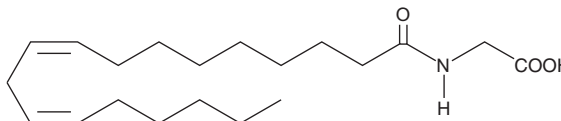
PRODUCT INFORMATION



Linoleoyl Glycine

Item No. 9000326

CAS Registry No.: 2764-03-6
Formal Name: N-9Z,12Z-1-oxo-octadecadien-1-yl-glycine
Synonyms: Glycine Linoleamide, LinGly, N-Linoleoyl Glycine
MF: C₂₀H₃₅NO₃
FW: 337.5
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

Linoleoyl Glycine (LinGly) is supplied as a crystalline solid. A stock solution may be made by dissolving the LinGly in the solvent of choice, which should be purged with an inert gas. LinGly is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of LinGly in these solvents is approximately 25, 15, and 20 mg/ml, respectively.

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. Organic solvent-free aqueous solutions of LinGly can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of LinGly in PBS (pH 7.2) is approximately 2 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

LinGly is an endogenous homolog of linoleoyl ethanolamide, an arachidonyl glycine. It inhibits the hydrolysis of AEA in FAAH-containing N18TG2 cell membranes with an IC₅₀ value of ~25 μM, which is less potent than that of arachidonyl glycine (IC₅₀ = 7 μM).¹ The biological significance of LinGly has not yet been determined.

Reference

1. Huang, S.M., Bisogno, T., Petros, T.J., *et al.* Identification of a new class of molecules, the arachidonyl amino acids, and characterization of one member that inhibits pain. *J. Biol. Chem.* **276(46)**, 42639-42644 (2001).

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, 04/10/2024

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