

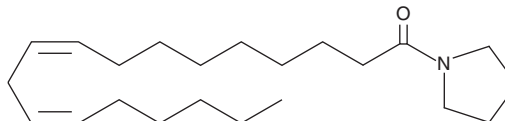
PRODUCT INFORMATION



Pyrrolidine Linoleamide

Item No. 17539

CAS Registry No.: 3140-51-0
Formal Name: 1-(1-pyrrolidinyl)-9Z,12Z-octadecadien-1-one
MF: C₂₂H₃₉NO
FW: 333.6
Purity: ≥98%
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

Pyrrolidine linoleamide is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of pyrrolidine linoleamide in these solvents is approximately 25 and 30 mg/ml, respectively.

Pyrrolidine linoleamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of pyrrolidine linoleamide should be diluted with the aqueous buffer of choice. Pyrrolidine linoleamide has a solubility of 0.25 mg/ml mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Linoleic acid (Item No. 90150) is an essential fatty acid and one of the most abundant polyunsaturated fatty acids in the western diet. Pyrrolidine linoleamide is a derivative of the amide of linoleic acid that shows potent antiproliferative activity against an array of cancer cell lines, including human glioma U251 cells.¹ It is greater than 4-fold more effective against cancer cells than non-cancer cells.¹

Reference

1. dos Santos, D.S., Piovesan, L.A., D'Oca, C.R.M., *et al.* Antiproliferative activity of synthetic fatty acid amides from renewable resources. *Bioorg. Med. Chem.* **23(2)**, 340-347 (2015).

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, 02/02/2024

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