

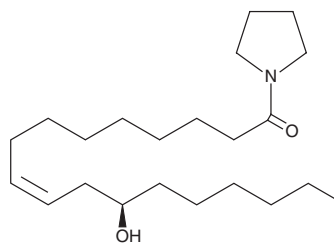
PRODUCT INFORMATION



Pyrrolidine Ricinoleamide

Item No. 17548

CAS Registry No.: 1246776-23-7
Formal Name: 12R-hydroxy-1-(1-pyrrolidinyl)-9Z-octadecen-1-one
MF: C₂₂H₄₁NO₂
FW: 351.6
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 ricinoleamide 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 ricinoleamide in these solvents is approximately 25 and 30 mg/ml, respectively.

Pyrrolidine ricinoleamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of pyrrolidine ricinoleamide should be diluted with the aqueous buffer of choice. Pyrrolidine ricinoleamide has a solubility of 0.25 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

Ricinoleic acid is a naturally occurring 12-hydroxy fatty acid. It constitutes about 90% of the fatty acids in castor oil. Pyrrolidine ricinoleamide is a derivative of the amide of ricinoleic acid that shows potent antiproliferative activity against an array of cancer cell lines, including human glioma U251 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