

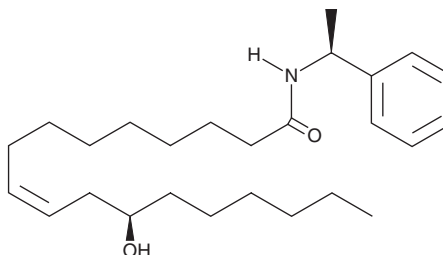
PRODUCT INFORMATION



(S)- α -Methylbenzyl Ricinoleamide

Item No. 17546

CAS Registry No.: 1246776-22-6
Formal Name: 12R-hydroxy-N-[(1S)-1-phenylethyl]-9Z-octadecenamide
MF: C₂₆H₄₃NO₂
FW: 401.6
Purity: \geq 98%
Supplied as: A 10 mg/ml solution in ethanol
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

(S)- α -Methylbenzyl 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 (S)- α -methylbenzyl ricinoleamide in these solvents is approximately 25 and 30 mg/ml, respectively.

(S)- α -Methylbenzyl ricinoleamide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of (S)- α -methylbenzyl ricinoleamide should be diluted with the aqueous buffer of choice. (S)- α -Methylbenzyl 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

(S)- α -Methylbenzyl ricinoleamide is a fatty acid amide derived from ricinoleic acid and methyl benzylamine.¹ It demonstrates potent growth inhibition of glioma (U251), breast (MCF-7), ovarian (NCI-ADR/RES and OVCAR-3), kidney (786-0), non-small cell lung (NCI-H460), and prostate (PC-3) cancer cells with a mean GI₅₀ value of 6.9 μ M.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, 05/26/2026

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