

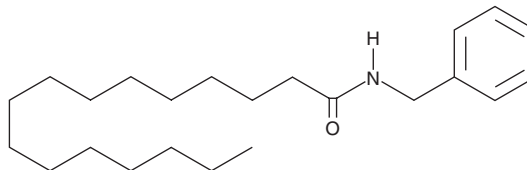
PRODUCT INFORMATION



N-Benzylpalmitamide

Item No. 9002235

CAS Registry No.: 74058-71-2
Formal Name: N-(phenylmethyl)-
hexadecanamide
Synonym: N-Benzylhexadecanamide
MF: C₂₃H₃₉NO
FW: 345.6
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

N-Benzylpalmitamide is supplied as a crystalline solid. A stock solution may be made by dissolving the N-benzylpalmitamide in the solvent of choice, which should be purged with an inert gas. N-Benzylpalmitamide is soluble in the organic solvent ethanol at a concentration of approximately 2 mg/ml.

Description

N-Benzylpalmitamide is a long-chain fatty acid amide (macamide or macaene) isolated from the maca (*L. meyenii*) plant and is structurally related to cannabinoids.¹ N-Benzylpalmitamide has been the most frequently isolated of the 19 macamides currently identified. While many macamides have been identified as potent inhibitors of fatty acid amide hydrolase (FAAH), N-benzylpalmitamide displays only moderate FAAH inhibitory activity (44% inhibition at 500 μM).¹ Additionally, many members of this family demonstrate selective antiproliferative activity against diverse cancer cell lines.²

References

1. Wu, H., Kelley, C.J., Pino-Figueroa, A., *et al.* Macamides and their synthetic analogs: Evaluation of in vitro FAAH inhibition. *Bioorg. Med. Chem.* **21(17)**, 5188-5197 (2013).
2. 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, 05/01/2024

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