

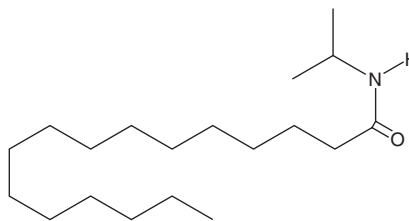
PRODUCT INFORMATION



Palmitoyl N-Isopropylamide

Item No. 91354

CAS Registry No.: 189939-61-5
Formal Name: N-isopropyl-hexadecanamide
Synonyms: PIA
MF: C₁₉H₃₉NO
FW: 297.5
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

Palmitoyl N-isopropylamide (PIA) is supplied as a crystalline solid. A stock solution may be made by dissolving the PIA in the solvent of choice, which should be purged with an inert gas. PIA is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of PIA in these solvents is approximately 20 and 3 mg/ml, respectively.

Description

PIA is a synthetic analog of palmitoyl ethanolamide which incorporates isopropyl amide in place of the native ethanolamide. PIA does not bind appreciably to either CB1 or CB2 receptors.¹ PIA inhibits both the transport (re-uptake) of tritiated AEA into intact cells, as well as the subsequent hydrolysis of AEA to arachidonate and ethanolamine, with 50% inhibition of FAAH-mediated hydrolysis at 10 μM.¹

Reference

1. Jonsson, K.O., Vandevoorde, S., Lambert, D.M., *et al.* Effects of homologues and analogues of palmitoylethanolamide upon the inactivation of the endocannabinoid anandamide. *Br. J. Pharmacol.* **133**(8), 1263-1275 (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, 04/25/2024

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