

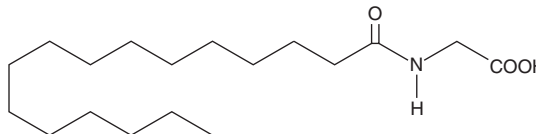
PRODUCT INFORMATION



N-Palmitoyl Glycine

Item No. 10009020

CAS Registry No.: 2441-41-0
Formal Name: N-(1-oxohexadecyl)-glycine
Synonyms: N-Hexadecanoyl-Glycine, PalGly
MF: C₁₈H₃₅NO₃
FW: 313.5
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.

Description

The acyl amides are a family of endogenous lipids that act as potent modulators of pain and inflammation. The best characterized members of this family are the arachidonoyl amides, which includes N-arachidonoyl ethanolamide (AEA; anandamide).

N-palmitoyl glycine (PalGly) contains an 18-carbon saturated fatty acid that is amide-linked to glycine and is structurally similar to the phospholipid-derived N-acyl ethanolamines. Endogenously produced in rat skin and spinal cord, PalGly is present in 100-fold greater amounts in skin and 3-fold greater in brain compared to AEA.¹ Injection of 0.43 μg PalGly in rat hindpaw inhibits heat-induced firing of nociceptive neurons in rat dorsal horn.¹ PalGly treatment induces transient calcium influx in native dorsal root ganglion (DRG) cells and in the PTX-sensitive, DRG-like cell line F-11 (EC₅₀ = 5.5 μM).¹

Reference

1. Rimmerman, N., Bradshaw, H.B., Hughes, H.V., *et al.* N-palmitoyl glycine, a novel endogenous lipid that acts as a modulator of calcium influx and nitric oxide production in sensory neurons. *Mol. Pharmacol.* **74**(1), 213-224 (2008).

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, 12/14/2022

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