

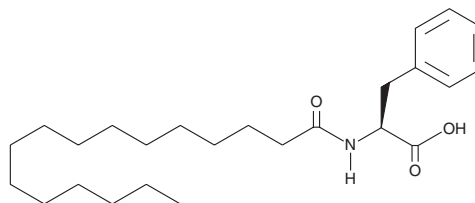
PRODUCT INFORMATION



N-Palmitoyl Phenylalanine

Item No. 20066

CAS Registry No.: 37571-96-3
Formal Name: N-(1-oxohexadecyl)-L-phenylalanine
Synonym: N-Hexadecanoyl Phenylalanine
MF: $C_{25}H_{41}NO_3$
FW: 403.6
Purity: $\geq 95\%$
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

N-Palmitoyl phenylalanine is supplied as a crystalline solid. A stock solution may be made by dissolving the N-palmitoyl phenylalanine in the solvent of choice, which should be purged with an inert gas. N-Palmitoyl phenylalanine is sparingly soluble (1-10 mg/ml) in ethanol.

Description

N-Palmitoyl phenylalanine is an N-acyl amide that contains the long-chain saturated fatty acid palmitic acid (Item No. 10006627) conjugated via an amide bond to the essential amino acid L-phenylalanine (Item No. 31498). It inhibits pre-mRNA splicing in HeLa cell nuclear lysates ($IC_{50} = \sim 150 \mu\text{M}$) but has no splicing activity in intact HeLa cells.¹

Reference

1. Effenberger, K.A., James, R.C., Urabe, V.K., et al. The natural product N-palmitoyl-L-leucine selectively inhibits late assembly of human spliceosomes. *J. Biol. Chem.* **290**(46), 27524-27531 (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, 09/17/2024

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