

# PRODUCT INFORMATION



## N-Oleoyl-DPPE (ammonium salt)

Item No. 39194

**Formal Name:** hexadecanoic acid, [R-(Z)]-1-(3-hydroxy-3-oxido-8-oxo-2,4-dioxo-7-aza-3-phosphapentacos-16-en-1-yl)-1,2-ethanediyl ester, monoammonium salt

**Synonyms:** NAPE(16:0/16:0/18:1), N-Oleoyl-1,2-Dipalmitoyl-3-PE, N-Oleoyl-1,2-Dipalmitoyl-3-Phosphoethanolamine, N-Oleoyl-1,2-Dipalmitoyl-3-Phosphatidylethanolamine

**MF:** C<sub>55</sub>H<sub>105</sub>NO<sub>9</sub>P • NH<sub>4</sub><sup>+</sup>

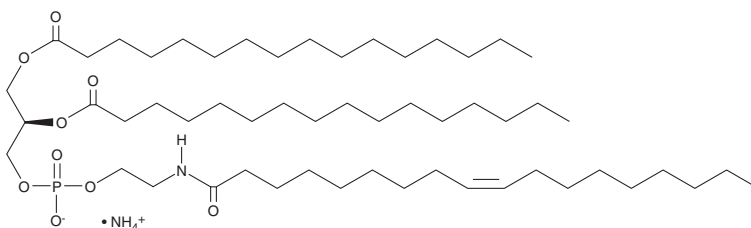
**FW:** 973.5

**Purity:** ≥95%

**Supplied as:** A 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-Oleoyl-DPPE (ammonium salt) is supplied as a solid. A stock solution may be made by dissolving the N-oleoyl-DPPE (ammonium salt) in the solvent of choice, which should be purged with an inert gas. N-Oleoyl-DPPE (ammonium salt) is soluble in the organic solvent chloroform at a concentration of approximately 5 mg/ml.

### Description

N-Oleoyl-DPPE is an N-acyl-phosphatidylethanolamine (NAPE).<sup>1</sup> It is a biosynthetic precursor to the endocannabinoid derivative and peroxisome proliferator-activated receptor  $\alpha$  (PPAR $\alpha$ ) agonist oleoyl ethanolamide (OEA; Item No. 90265).<sup>2</sup>

### References

1. Akoka, S., Tellier, C., Le Roux, C., *et al.* A phosphorus magnetic resonance spectroscopy and a differential scanning calorimetry study of the physical properties of N-acylphosphatidylethanolamines in aqueous dispersions. *Chem. Phys. Lipids* **46(1)**, 43-50 (1988).
2. Orío, L., Alen, F., Pavón, F.J., *et al.* Oleylethanolamide, neuroinflammation, and alcohol abuse. *Front. Mol. Neurosci.* **11**, 490 (2019).

#### 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, 11/07/2023

#### 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