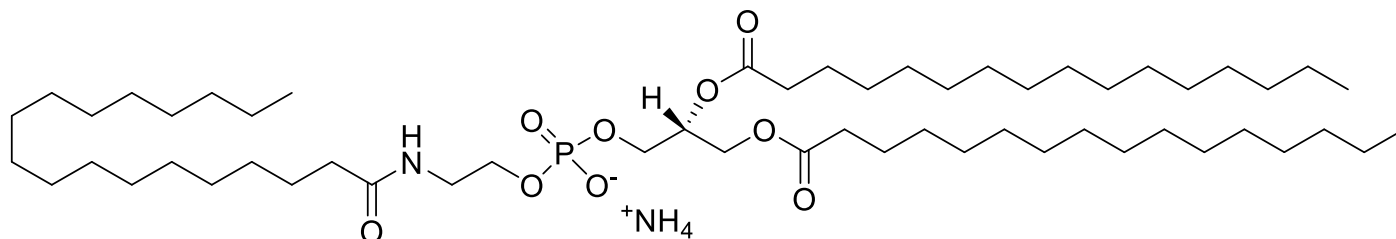


## N-Stearoyl-DPPE (1,2-Dipalmitoyl-sn-glycero-3-phosphoethanolamine-N-stearoyl) (ammonium salt)

Catalog number: L-2418



Molecular Formula: C<sub>55</sub>H<sub>111</sub>N<sub>2</sub>O<sub>9</sub>P

MW: 975.47

CAS: 113701-56-7

Alternate Name: Stearoyl-NAPE, 16:0-PE-N-18:0

Solubility: CHCl<sub>3</sub> (5 mg/mL)

Storage and Handling: Store dry at -20 °C. Stock solutions should be stored frozen (-20 °C or below).

Background: N-acylphosphatidylethanolamines (NAPEs) are important intermediates in the eicocannabinoid biosynthesis pathway, producing anandamide through hydrolysis by NAPE-specific phospholipase D. NAPEs are also produced in the small intestine, travel through the circulatory system to the brain, accumulate in the hypothalamus, and suppress appetite. Injected NAPEs cause a reduction in food intake and weight loss in rats.

References: 1) J. Wang, N. Ueda "Biology of the eicocannabinoid synthesis system" *Prostaglandins and other Lipid Mediators*, 2009, *89*, 112-119.

2) M. P Gillum et al. "N-acylphosphatidylethanolamine, a gut-derived circulating factor induced by fat ingestion, inhibits food intake" *Cell*, 2008, *135*, 813-824.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. This product is manufactured and shipped only in small quantities, intended for research and development in a laboratory utilizing prudent procedures for handling chemicals of unknown toxicity, under the supervision of persons technically qualified to evaluate potential risks and authorized to enforce appropriate health and safety measures. As with all research chemicals, precautions should be taken to avoid unnecessary exposures or risks.

Warranty and Disclaimer: Echelon warrants the product conforms to the specifications stated herein. In the event of nonconformity, Echelon will replace products or refund purchase price, at its sole option, and Echelon shall not be responsible for any other loss or damage, whether known or foreseeable to Echelon. No other warranties apply, express or implied, including but not limited to warranty of fitness for any purpose or implied warranty of merchantability. Purchaser is solely responsible for all consequences of its use of the product and Echelon assumes no responsibility therefore, including success of purchaser's research and development. or health or safety of any uses of the product.

