

# PRODUCT DATA SHEET



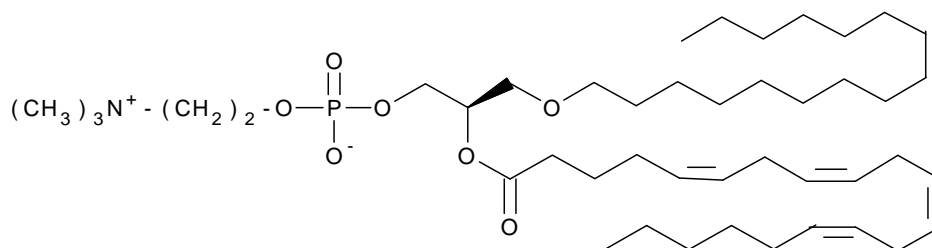
PRODUCT: 2-AA-PAF  
1-O-Hexadecyl-2-O-arachidonoyl-sn-glycero-3-phosphorylcholine

CAS NO: 86288-11-1

CATALOG NO.: L-102

LOT#: temp

STRUCTURE:



PHYSICAL APPEARANCE: pale yellow waxy solid in ethanol (25 mg/mL)

MOLECULAR FORMULA:  $\text{C}_{44}\text{H}_{82}\text{NO}_7\text{P}$

MOLECULAR WEIGHT: 768.2

PURITY: 98% (TLC)

STORAGE: Store, as supplied, unopened at  $-80^\circ\text{C}$  for up to 1 year. Store opened product under an inert atmosphere at  $-80^\circ\text{C}$  for up to 3 months.

APPLICATION NOTES: A biosynthetic precursor of PAF via lipid remodeling pathway.<sup>1</sup> Open ampule and aliquot into suitable tubes or vials. Product is subject to oxidation. Protect from air - store under an inert atmosphere. Store at  $-80^\circ\text{C}$  or colder. For experiments with cells or tissues, aliquots of ethanol solutions may be added directly to aqueous media

REFERENCES: 1. F.H.Chilton *et al.* J.Biol.Chem. 1984 **259** 12014

**The pharmacological and toxicological properties of this product have not been fully investigated. Exercise caution in use and handling. This product must not be used in humans.**