



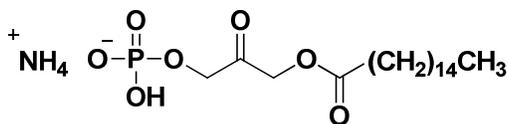
Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Product Name: Palmitoyl dihydroxyacetone phosphate, ammonium salt
(palmitoyl-DHAP)



Product Number: L-0310

Molecular Formula: C₁₉H₄₀NO₇P **MW:** 425.50

Solubility: CHCl₃:MeOH:H₂O (1:2:0.8) ; tert-Butyl Alcohol:H₂O (1:1)

Storage: -20 °C

Field of Interest: Ether lipids have been implicated in cell membrane function and can act as signaling molecules; however their roles in cell physiology are still not completely understood. The key step for ether lipid biosynthesis is the conversion of acyldihydroxyacetone phosphate (acyl-DHAP) to alkyl dihydroxyacetone phosphate (alkyl-DHAP) by the enzyme alkyl dihydroxyacetone phosphate synthase (ADPS, alternately referred to as AGPS). Recently, elevated levels of ADPS have been found in human cancer cells and primary tumors which lead to increased levels of oncogenic signaling lipids.

References: 1) Benjamin, D.I.; Cozzo, A. *et al.* Ether lipid generating enzyme AGPS alters the balance of structural and signaling lipids to fuel cancer pathogenicity. *Proc. Natl. Acad. Sci. USA* **110**, 2013. 14912-7

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. **Palmitoyl-DHAP** 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.