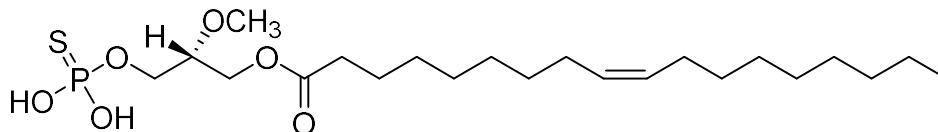


## *D*-sn-1-*O*-oleoyl-2-*O*-methyl-glycerol-3-phosphothionate (2*R*-OMPT)

**Catalog number:** L-9318



**Molecular Formula:** C<sub>22</sub>H<sub>43</sub>O<sub>6</sub>P

**MW:** 466.61

**CAS#:** 645408-60-2

**Solubility:** Methanol > 1 mg/mL

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

**Background:** OMPT is a metabolically-stabilized analog of Lysophosphatidic Acid (LPA) which acts as potent agonist for the LPA3 receptor. Of the two enantiomers, (2*S*)-OMPT is more potent than (2*R*)-OMPT showing 5-20 fold more activity in calcium release from LPA3-transfected Sf9 and rat hepatoma Rh7777 cells. (2*S*)-OMPT is also more potent in inducing calcium release, IL-6 production, and MAPK and Akt phosphorylation in OVCAR3 cells.

**References:** 1) Qian, L., Y. Xu, et al. (2003). "Enantioselective responses to a phosphorothioate analogue of lysophosphatidic acid with LPA3 receptor-selective agonist activity." *J Med Chem* 46(26): 5575-8.  
2) Altman, M. K., V. Gopal, et al. (2010). "Targeting melanoma growth and viability reveals dualistic functionality of the phosphonothionate analogue of carba cyclic phosphatidic acid." *Mol Cancer* 9: 140.

**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.

Technical Data Sheet, Rev 2, 6/27/2024 – **For research use only.** Not intended for diagnostic or therapeutic use.