

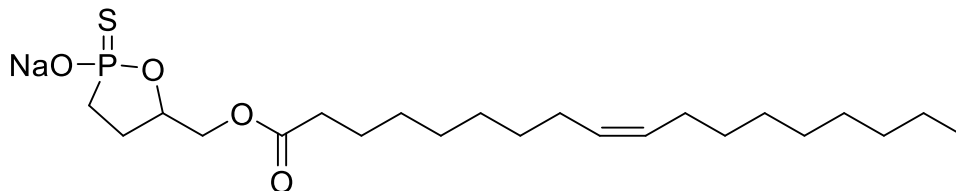
Thio-ccPA

Catalog number: L-7118

Molecular Formula: $C_{22}H_{40}NaO_4PS$

MW: 454.6

CAS: 779333-56-1



Alternate Name: 3-Hydroxy-4-oleoyloxybutane 1,3-Cyclic Phosphonothioate

Solubility: Methanol 1 mg/mL

Storage and Handling: Store dry at $-20^{\circ}C$. Stock solutions should be stored frozen ($-20^{\circ}C$ or below). Storage in basic (pH > 9.0) or acidic (pH < 4.0) buffers may cause decomposition.

Background: Thio-ccPA is a stabilized analogue of carbacyclic phosphatidic acid (ccPA). It inhibits the lysoPLD activity of Autotaxin (ATX) and acts as an LPA1/3 receptor antagonist. Thio-ccPA decreases the viability of B16F10, A375, and MeWo melanoma cell lines. Addition of endogenous LPA did not reverse the effects of Thio-ccPA in B16F10 cells. Thio-ccPA decreased the number of pulmonary and non-pulmonary metastases in a B16F10 metastatic melanoma mouse model.

References: 1) M.K. Altman, *et al.* "Targeting melanoma growth and viability reveals dualistic functionality of the phosphonothionate analogue of carbacyclic phosphatidic acid" *Molecular Cancer*, 2010, 9, 140.

2) Y. Xu, *et al.* "Phosphonothioate and fluoromethylene phosphonate analogues of cyclic phosphatidic acid: novel antagonists of lysophosphatidic acid receptors" *J. Med. Chem.*, 2006, 49, 5309.

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.

