

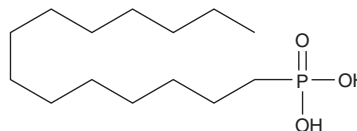
# PRODUCT INFORMATION



## Tetradecyl Phosphonate

Item No. 10007565

**CAS Registry No.:** 4671-75-4  
**Formal Name:** tetradecylphosphonic acid  
**MF:** C<sub>14</sub>H<sub>31</sub>O<sub>3</sub>P  
**FW:** 278.4  
**Purity:** ≥98%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Tetradecyl phosphonate is supplied as a crystalline solid. A stock solution may be made by dissolving the tetradecyl phosphonate in the solvent of choice, which should be purged with an inert gas. Tetradecyl phosphonate is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of tetradecyl phosphonate in these solvents is approximately 30, 2, and 5 mg/ml, respectively.

Tetradecyl phosphonate is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, tetradecyl phosphonate should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. Tetradecyl phosphonate has a solubility of approximately 0.25 mg/ml in a 1:3 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Lysophosphatidic acid (LPA) is a lipid signaling molecule formed by the hydrolysis of lysophosphatidyl choline by lysophospholipase D, also known as autotaxin (ATX).<sup>1</sup> LPA signals through four different G protein-coupled receptors, LPA<sub>1</sub>/EDG-2, LPA<sub>2</sub>/EDG-4, LPA<sub>3</sub>/EDG-7, and LPA<sub>4</sub>/GPR23.<sup>2,3</sup> Activation of peroxisome proliferator-activated receptor  $\gamma$  (PPAR $\gamma$ ) by LPA has also been reported.<sup>4</sup> Tetradecyl phosphonate is a pan-antagonist of LPA<sub>1</sub>, LPA<sub>2</sub>, and LPA<sub>3</sub> receptors with IC<sub>50</sub> values for inhibition of LPA-induced calcium mobilization of 10  $\mu$ M, 5.5  $\mu$ M, and 3.1  $\mu$ M, respectively.<sup>5</sup> At a concentration of 10  $\mu$ M, tetradecyl phosphonate activates a PPAR $\gamma$  reporter construct 4-fold compared to controls and partially inhibits ATX with an IC<sub>50</sub> of approximately 3  $\mu$ M.<sup>5</sup>

### References

1. Tokumura, A., Majima, E., Kariya, Y., *et al.* Identification of human lysophospholipase D, a lysophosphatidic acid-producing enzyme, as autotaxin, a multifunctional phosphodiesterase. *J. Biol. Chem.* **277(42)**, 39436-39442 (2002).
2. Chun, J., Goetzl, E.J., Hla, T., *et al.* International union of pharmacology. XXXIV. Lysophospholipid receptor nomenclature. *Pharmacol. Rev.* **54**, 265-269 (2002).
3. Noguchi, K., Ishii, S., and Shimizu, T. Identification of p2y9/GPR23 as a novel G protein-coupled receptor for lysophosphatidic acid, structurally distant from the Edg family. *J. Biol. Chem.* **278(28)**, 25600-25606 (2003).
4. McIntyre, T.M., Pontsler, A.V., Silva, A.R., *et al.* Identification of an intracellular receptor for lysophosphatidic acid (LPA): LPA is a transcellular PPAR $\gamma$  agonist. *Proc. Natl. Acad. Sci. USA* **100(1)**, 131-136 (2003).
5. Durgam, G.G., Virag, T., Walker, M.D., *et al.* Synthesis, structure-activity relationships, and biological evaluation of fatty alcohol phosphates as lysophosphatidic acid receptors ligands, activators of PPAR $\gamma$ , and inhibitors of autotaxin. *J. Med. Chem.* **48**, 4919-4930 (2005).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/31/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM