

Echelon Biosciences Inc.

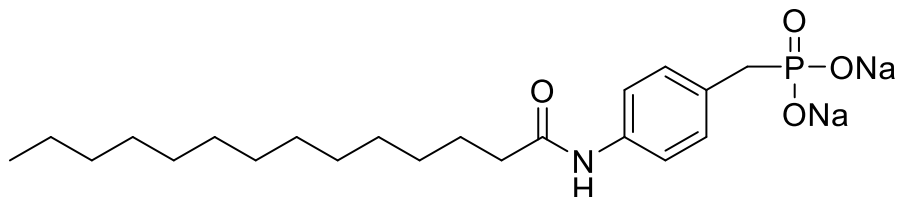
S32826

Catalog number: L-3282

Molecular Formula: $C_{21}H_{34}NNa_2O_4P$

MW: 441.20

CAS: 1103672-43-0



Alternate Name: [4-(Tetradecanoylamino)benzyl]phosphonic acid disodium salt

Solubility: water, 1 mg/mL, may require sonication for complete dissolution

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

Background: S32826 is a potent inhibitor of the enzyme autotaxin, with an IC_{50} in the nanomolar range¹. S32826 is suitable for *in vitro* studies using purified components or cultured cells. Due to its short circulation half-life¹, S32826 is not recommended for use in animal models.

References: 1) Ferry, G., Moulharat, N., Pradere, J.P., Desos, P., Try, A., Genton, A., Giganti, A., Beucher-Gaudin, M., Lonchamp, M., Bertrand, M., Saulnier-Blache, J.S., Tucker, G.C., Cordi, A., & Boutin, J.A., S32826, a nanomolar inhibitor of autotaxin: discovery, synthesis and applications as a pharmacological tool. *J Pharmacol Exp Ther* 327 (3), 809-819 (2008).

2) Samadi, N., Gaetano, C., Goping, I.S., & Brindley, D.N., Autotaxin protects MCF-7 breast cancer and MDA-MB-435 melanoma cells against Taxol-induced apoptosis. *Oncogene* 28 (7), 1028-1039 (2009).

3) Conrotto, P, *et al.* Knock-down of SOX11 induces autotaxin-dependent increase in proliferation *in vitro* and more aggressive tumors *in vivo* *Mol. Oncology* 5 (7), 527 (2011).

4) Li, S. Xiong, C., Zhang, J. ATX and LPA receptor 3 are coordinately up-regulated in lipopolysaccharide-stimulated THP-1 cells through PKR and SPK1-mediated pathways. *FEBS Lett.* 586 (6), 792 (2012).

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 4, 1-31-19 – For research use only. Not intended for diagnostic or therapeutic use.



Echelon Biosciences Inc.

Ph: 866-588-0455

Fax: 801-588-0497

Echelon-inc.com