



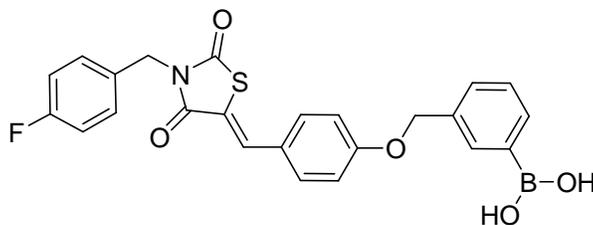
Echelon Biosciences Inc.
675 Arapeen Drive, Suite 302
Salt Lake City, UT 84108
Telephone 866-588-0455
Fax 801-588-0497
echelon@echelon-inc.com
www.echelon-inc.com

Technical Data Sheet

For research use only
Not intended or approved for
diagnostic or therapeutic use.

Product Name: HA130 (Autotaxin Inhibitor)

Product Number: B-0701



Molecular Formula: C₂₄H₁₉BFNO₅S

MW: 463.29

CAS: 1229652-21-4

Solubility: DMSO (>5 mg/mL)

Storage: -20°C or lower

Field of Interest: HA130 is a boronic acid based inhibitor of the lysoPLD activity of autotaxin (ATX). HA130 is a potent (IC₅₀ = 28 nM), reversible inhibitor which interacts with the oxygen in active site residue T210. HA130 is selective against ATX and does not show any activity against NPP1, alkaline phosphatase, phosphodiesterase, at concentrations up to 10 μM. In addition, proteosomal chymotryptic, caspase, and typtic activities were not affected. It also showed minimal toxicity against A2058, HEK293T, and HepG2 cell lines (TD₅₀ = 105, 83, & 2056 μM respectively). Plasma ATX activity was inhibited and circulating LPA levels decreased when administered to mice.

References: 1) H.M.H.G. Albers *et al.* "Boronic acid-based inhibitor of autotaxin reveals rapid turnover of LPA in the circulation" *Proc. Natl. Acad. Sci. USA*, **2010**, *107*, 7257-7262. 2) H.M.H.G. Albers *et al.* "Discovery and Optimization of Boronic Acid Based Inhibitors of Autotaxin" *J. Med. Chem.*, **2010**, *53*, 4958-4967.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. HA130 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.