

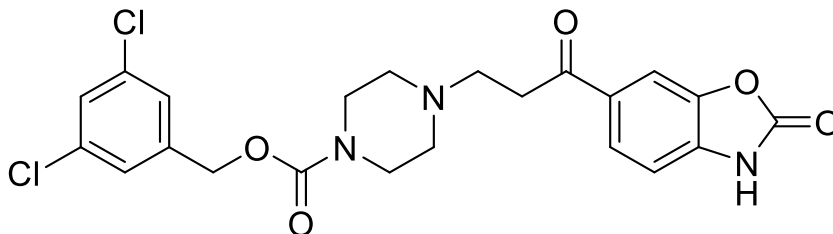
## PF-8380 (ATX Inhibitor)

Catalog number: B-0702

Molecular Formula:  $C_{22}H_{21}Cl_2N_3O_5$

MW: 478.33

CAS: 1144035-53-9



Alternate Name: 3,5-Dichlorobenzyl-[4-[3-oxo-3-(2-oxo-2,3-dihydrobenzoxazol-6-yl)propyl]]piperazine-1-carboxylate

Solubility: DMSO >1 mg/mL

Storage and Handling: Store dry at 4 °C. Stock solutions should be stored frozen (-20 °C or below).

Background: PF-8380 is a potent autotaxin (ATX) inhibitor ( $IC_{50}$  2.8 nM/101 nM purified enzyme/whole blood). Excessive activity of ATX is associated with numerous inflammatory conditions, cancers and other disease. Lysophosphatidic acid (LPA) produced by ATX stimulates cell proliferation through G-protein coupled receptors LPA1-4. PF-8380 is orally bioavailable and effectively inhibits ATX in vivo reducing LPA levels in serum and at the site of inflammation in a rat arthritis model.

References: 1) Gierse, J., *et al.* "A novel autotaxin inhibitor reduces lysophosphatidic acid levels in plasma and the site of inflammation", *J Pharmacol Exp Ther* 334: 310-7, (2010) 2) St. Coeur, *et al.* "PF-8380 and Closely Related Analogs: Synthesis and Structure-Activity Relationship towards Autotaxin Inhibition and Glioma Cell Viability" *Arch. Pharm.* 346: 91-97 (2015)

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.

