

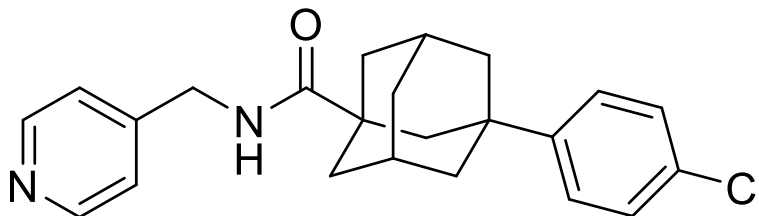
ABC294640 (SPHK2 Inhibitor)

Catalog number: B-0025

Molecular Formula: $C_{23}H_{25}ClN_2O$

MW: 380.91

CAS: 915385-81-8



Solubility: DMSO, > 5mg/mL

Storage and Handling: Store as a solid at -20 °C or below. Stock solutions should be stored at -20 °C.

Background: ABC294640 is an orally bioavailable, selective inhibitor of sphingosine kinase 2 (SPHK2 or SK2). Sphingosine kinases phosphorylate sphingosine producing the bioactive lipid sphingosine 1-phosphate which has roles in cell proliferation and cell motility. Sphingosine kinases are upregulated in numerous forms of cancer and ABC294640, alone or in combination, has been shown to suppress growth in kidney, breast, liver, colon, and lung cancer cell lines.

References: 1) M. B. Thomas, *et al.* "Phase I trial of ABC294640, a first-in-class sphingosine kinase-2 inhibitor" *Mol. Cancer. Ther.* 2013, 12, C62, DOI: 10.1158/1535-7163.TARG-13-C62. 2) K.J French, *et al.* "Pharmacology and antitumor activity of ABC294640, a selective inhibitor of sphingosine kinase-2." *J. Pharmacol. Exp. Ther.* 2010, 333, 129-139. 3) M.M. Gestaut, *et al.* "Inhibition of sphingosine kinase-2 ablates androgen resistant prostate cancer proliferation and survival." *Pharmacol. Rep.* 2014, 66, 174-178.

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.

