

SKI-178 (Sphingosine Kinase 1 Inhibitor)

Catalog number: B-0028

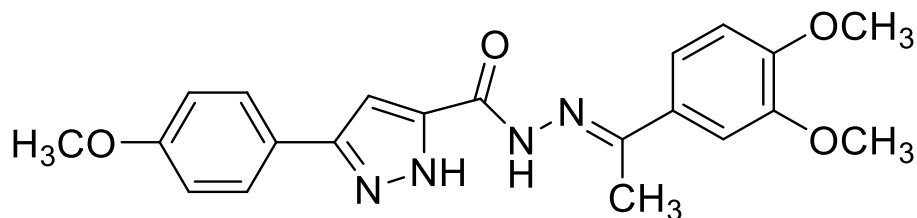
Molecular Formula: C₂₁H₂₂N₄O₄

MW: 394.43

CAS: 1259484-97-3

Solubility: DMSO >5 mg/mL

Storage and Handling: -20 °C or below



Background: SKI-178 is a selective inhibitor of sphingosine kinase 1 which is active *in vitro* and *in vivo*. SKI-178 is a competitive inhibitor of sphingosine binding to SPHK1 ($K_i = 1.3 \mu\text{M}$) and shows no inhibition of SPHK2 up to 25 μM . SKI-178 is cytotoxic to numerous cancer cell lines (MCF-7, A549, Panc-1, etc) at low to sub-micromolar concentrations.

References: 1) J.A. Hengst *et al.* "Development of a sphingosine kinase 1 specific small-molecule inhibitor" *Bioorg. Med. Chem. Lett.* 2010, 20, 7498-7502. doi: 10.1016/j.bmcl.2010.10.005 2) T.E. Dick *et al.* "The apoptotic mechanism of action of the sphingosine kinase 1 selective inhibitor SKI-178 in human acute myeloid leukemia cell lines" *J. Pharmacol. Exp. Ther.* 2015, 352, 494-508. doi: 10.1124/jpet.114.219659

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.

