



Echelon Biosciences Inc.
675 Arapeen Drive, Ste 302
Salt Lake City, UT 84108
Phone: 801-588-0455
Fax: 801-588-0497
Email: echelon@echelon-inc.com
<http://www.echelon-inc.com>

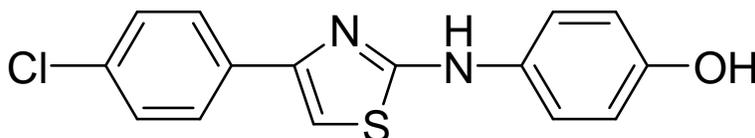
Technical Data Sheet

For research use only

Not intended or approved for
diagnostic or therapeutic use.

Product Name: 2-(*p*-Hydroxyanilino)-4-(*p*-chlorophenyl)thiazole, HCl
Sphingosine Kinase Inhibitor

Product Number: B-0024



Formula: C₁₅H₁₁ClN₂OS·HCl, FW: 339.2

Solubility: water: insoluble, DMSO: 5 mg/mL

Storage: store dry between 2-8°C. Stock solutions should be stored frozen (-20°C).

Field of Interest: 2-(*p*-Hydroxyanilino)-4-(*p*-chlorophenyl)thiazole is a cell permeable, potent, and specific inhibitor of sphingosine kinase (IC₅₀ = 0.5 μM). The compound does not bind to the ATP-binding site and does not affect the kinase activities of hERK2, hPI3K, or hPKCα at concentrations up to 60 μM. The compound showed anti-tumor properties, inducing apoptosis in a number of cell lines including those that express Pgp or MRP1 drug-transport proteins. The product is not sterile.

Reference: French, K.J. *et al. Cancer Res.* **63**, 2003, 5969.

Hazardous Properties and Cautions: The toxicological and pharmacological properties of this compound are not fully known. For further information see the MSDS on request. 2-(*p*-Hydroxyanilino)-4-(*p*-chlorophenyl)thiazole 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.