

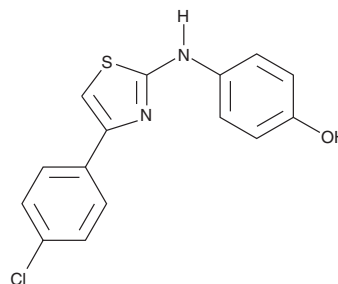
# PRODUCT INFORMATION



## Sphingosine Kinase Inhibitor 2

Item No. 10009222

**CAS Registry No.:** 312636-16-1  
**Synonyms:** SKI II, SPHK I2  
**Formal Name:** 4-[[4-(4-chlorophenyl)-2-thiazolyl]amino]-phenol  
**MF:** C<sub>15</sub>H<sub>11</sub>ClN<sub>2</sub>OS  
**FW:** 302.8  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 259, 282 nm  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

SPHK I2 is supplied as a crystalline solid. A stock solution may be made by dissolving the SPHK I2 in an organic solvent purged with an inert gas. SPHK I2 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of SPHK I2 in these solvents is approximately 20 mg/ml.

SPHK I2 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SPHK I2 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. SPHK I2 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Sphingosine kinase inhibitor 2 is an inhibitor of sphingosine kinase 1 (SPHK1) with an IC<sub>50</sub> value of 0.5 μM for non-ATP-competitive inhibition of human recombinant SPHK1.<sup>1</sup> It is selective for SPHK1 over ERK2, PI3K, and PKCα at concentrations up to 60 μM. Sphingosine kinase inhibitor 2 inhibits proliferation of several cancer cell lines, including T-24, MCF-7, NCI/ADR, and MCF-7/VP (IC<sub>50</sub>s = 0.9-4.6 μM).

### Reference

1. French, K.J., Schrecengost, R.S., Lee, B.D., et al. Discovery and evaluation of inhibitors of human sphingosine kinase. *Cancer Res.* **63(18)**, 5962-5969 (2003).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/25/2024

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

**PHONE:** [800] 364-9897  
[734] 971-3335

**FAX:** [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
[WWW.CAYMANCHEM.COM](http://WWW.CAYMANCHEM.COM)