

# SKI-II

## Sphingosine kinase inhibitor

Selective, non-ATP competitive inhibitor of sphingosine kinase ( $IC_{50}$  = 0.5 $\mu$ M GST-hSK;  $IC_{50}$  = 1.1 $\mu$ M intact JC mouse mammary adenovirus carcinoma cells) showing no inhibition of hERK2, hPI3k and hPKC $\alpha$  at concentrations up to 60 $\mu$ M. Cytotoxic effects have been demonstrated against various tumor cell lines including some overexpressing P-glycoprotein or MRP1 (T24,  $IC_{50}$  = 4.6 $\mu$ M; MCF-7,  $IC_{50}$  = 1.2 $\mu$ M; NCI/ADR,  $IC_{50}$  = 1.3 $\mu$ M, MCRF-7/VP,  $IC_{50}$  = 0.9 $\mu$ M). Induced apoptosis in T24 tumor cells.

Citations: 2

[View Online »](#)

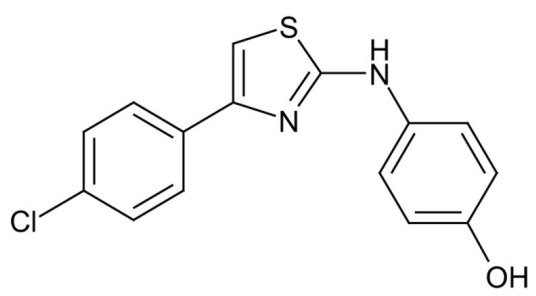
## Ordering Information

[Order Online »](#)

BML-EI406-0010	10mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light and air.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	4-[[4-(4-Chlorophenyl)-1,3-thiazol-2-yl]amino]phenol
Appearance	Off-white solid.
CAS	312636-16-1
Couple Target	Sphingosine kinase
Couple Type	Inhibitor
Formula	$C_{15}H_{11}ClN_2OS$
MW	302.8
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml) or 100% ethanol (>25mg/ml).
Source	Synthetic.



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)