

SKI-II

Sphingosine kinase inhibitor

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

Citations: 2

[View Online »](#)

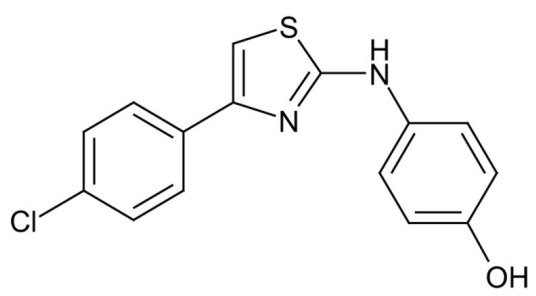
Ordering Information

[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.

Last modified: May 29, 2024



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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)