

SK1-I

Sphingosine kinase 1 inhibitor exclusive to Enzo Life Sciences

SK1-I is a sphingosine analog and a sphingosine competitive inhibitor specific for sphingosine kinase 1 (SK1), with $K_i \sim 10 \mu\text{M}$ and excellent water solubility. It is not to be confused with SKI-I, 5-naphthalen-2-yl-2*H*-pyrazole-3-carboxylic acid (2-hydroxy-naphthalen-1-ylmethylene)-hydrazide, CAS 306301-68-8, a noncompetitive inhibitor of both SK1 and SK2 with poor water solubility (K.J. French, et al., 2006; N.J. Pyne and S. Pyne, 2010). SK1-I does not inhibit SK2, PKC α , PKC δ , PKA, AKT1, ERK1, EGFR, CDK2, IKK β or CamK2 β . Not only does it decrease sphingosine-1-phosphate levels, it also causes an accumulation of its proapoptotic precursor ceramide. Inhibits tumor cell growth *in vitro* and *in vivo*.

Citations: 14

[View Online »](#)

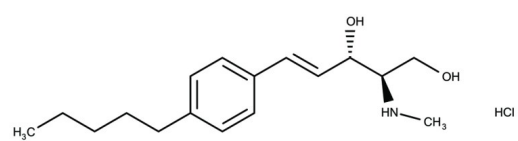
Ordering Information

[Order Online »](#)

BML-EI411-0005	5mg
BML-EI411-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C.
Handling	Protect from light and air.
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(2R,3S,4E)-N-methyl-5-(4'-pentylphenyl)-2-aminopent-4-ene-1,3-diol . HCl
Appearance	White solid.
CAS	1072443-89-0
Couple Target	Sphingosine kinase
Couple Type	Inhibitor
Formula	$C_{17}H_{27}NO_2 \cdot HCl$
Identity	Determined by NMR.
MW	313.9
Purity	≥98% (HPLC)
Solubility	Soluble in water or DMSO (100 mM).

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-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)