

# 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 $\alpha$ , PKC $\delta$ , PKA, AKT1, ERK1, EGFR, CDK2, IKK $\beta$  or CamK2 $\beta$ . 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 »](#)

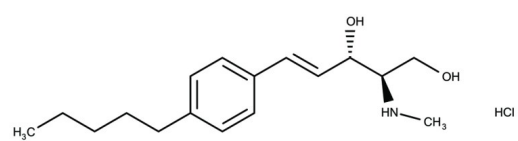
## Ordering Information

[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 <sub>17</sub> H <sub>27</sub> NO <sub>2</sub> . HCl
Identity	Determined by NMR.
MW	313.9
Purity	≥98% (HPLC)
Solubility	Soluble in water or DMSO (100 mM).



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)