

SKL2001

Wnt/ β -Catenin pathway activator

SKL2001 is a Wnt/ β -catenin pathway activator that disrupts the Axin/ β -catenin interaction leading to β -catenin stabilization. It suppressed colon cancer spheroid growth via regulation of the E-cadherin/ β -catenin complex. A commonly used reagent to reactivate or rescue the Wnt/ β -Catenin signaling pathway.

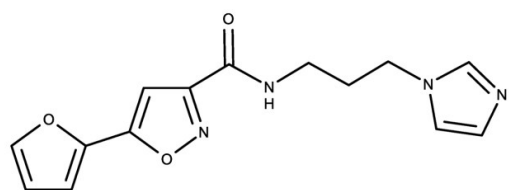
Ordering Information

[Order Online »](#)

ENZ-CHM452-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5-(Furan-2-yl)-N-(3-imidazol-1-ylpropyl)-1,2-oxazole-3-carboxamide

Appearance Beige solid.

CAS 909089-13-0

Couple Type Activator

Formula $C_{14}H_{14}N_4O_3$

Identity Determined by NMR.

MW 286.29

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (>25 mg/ml).

Last modified: January 24, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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