

SANT-2

Smoothened inhibitor

Cell permeable antagonist of the Smoothened (Smo) protein function. Inhibitor of SAG (Prod. No. ALX-270-426) activity. Chemical tool to study the Hedgehog (Hh) signalling pathway and to explore therapeutic uses for inhibitors of Smo.

Citations: 1

[View Online »](#)

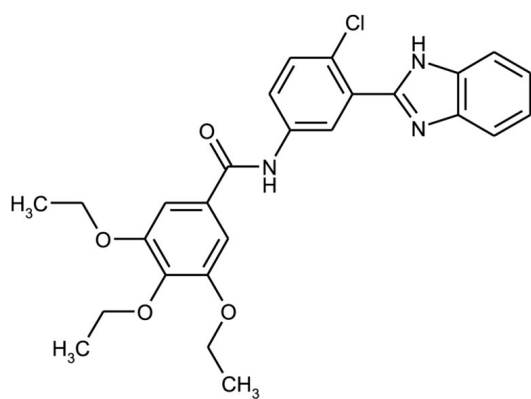
Ordering Information

[Order Online »](#)

ALX-270-436-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 6 months when stored at -20°C in the dark.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-[3-(1H-Benzimidazol-2-yl)-4-chlorophenyl]-3,4,5-triethoxybenzamide
Appearance	White to off-white solid.
CAS	329196-48-7
Couple Target	Smoothened
Couple Type	Inhibitor
Formula	$C_{26}H_{26}ClN_3O_4$
Identity	Identity determined by MS, 1H- and 13C-NMR.
MW	480.0
Purity	≥95%
Solubility	Soluble in DMSO.



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)