

# SANT-2

## Smoothened inhibitor

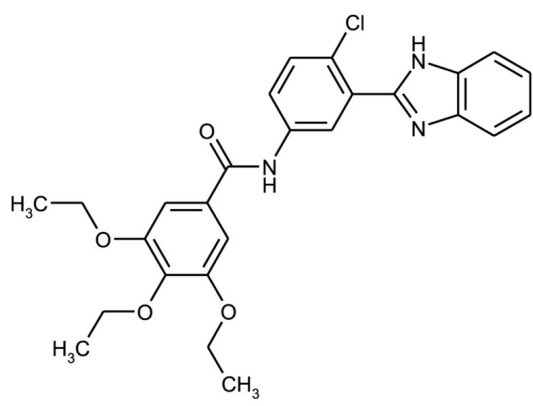
Cell permeable antagonist of the Smoothened (Smo) protein function. Inhibitor of SAG (Prod. No. ALX-270-426) activity. Chemical tool to study the Hedgehog (Hh) signalling pathway and to explore therapeutic uses for inhibitors of Smo.

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Protect from light.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

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

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-be@enzolifesciences.com](mailto:info-be@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)