

SAG

Smoothened agonist

Smoothened (Smo) is a transmembrane receptor that relays the Hedgehog (Hh) signal and plays an essential role during embryonic development and tumorigenesis. SAG is a potent cell-permeable benzothiophene compound that modulates the coupling of Smo with its downstream effector by interacting with the Smo heptahelical domain ($K_D = 59 \text{ nM}$). SAG was shown to induce Hedgehog pathway activation independently of the Hh binding receptor Patched (Ptch) ($EC_{50} = \sim 3 \text{ nM}$ in NIH 3T3-derived Shh-LIGHT2 cells) and counteract inhibition of Smo by Cyclopamine (Prod. No. BML-GR334). Smo is reported to act as an activator at low concentrations and as an inhibitor at very high concentrations.

Citations: 60

[View Online »](#)

Ordering Information

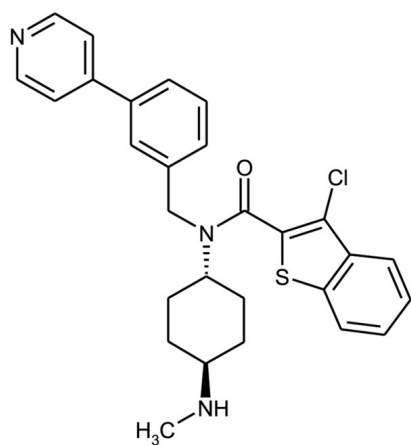
[Order Online »](#)

ALX-270-426-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)

- Potent Smoothened (Smo) receptor agonist
- Binds directly to Smo and antagonizes Cyclopamine
- Enhancer of neuronal differentiation of iPSCs into dopaminergic neurons



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions in DMSO are stable for at least 1 month when stored at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-chloro-N-[trans-4-(methyamino)cyclohexyl]-N-[[3-(4-pyridinyl)phenyl]methyl]-benzo[b]thiophene-2-carboxamide
Appearance	White to off-white solid.
CAS	912545-86-9
Couple Target	Smoothened
Couple Type	Activator, Ligand
Formula	$C_{28}H_{28}ClN_3OS$
Identity	Identity determined by MS and NMR.
MW	490.1
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (10mg/ml), sparingly soluble in water (acidification with HCl increases water solubility).



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)