

GANT61

GLI inhibitor

Inhibitor of Hedgehog (Hh) signalling pathway. Blocks the GLI function which constitutes the final step in the Hh pathway. Interferes with GLI1 DNA binding in living cells. Inhibits cell proliferation of tumor cells *in vitro* in a GLI-dependent manner. Blocks cell growth in an *in vivo* xenograft model using human prostate cancer cells.

Citations: 18

[View Online »](#)

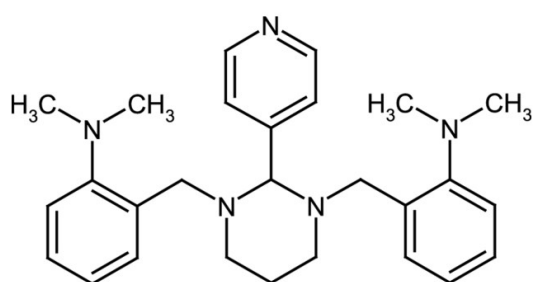
Ordering Information

[Order Online »](#)

ALX-270-482-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	NSC 136476, 2,2'-[[Dihydro-2-(4-pyridinyl)-1,3(2H,4H)-pyrimidinediyl]bis(methylene)]bis[N,N-dimethyl]-benzenamine
Appearance	White to off-white solid.
CAS	500579-04-4
Couple Target	GLI
Couple Type	Inhibitor
Formula	$C_{27}H_{35}N_5$
MW	429.6
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (5mg/ml).

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-
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)