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

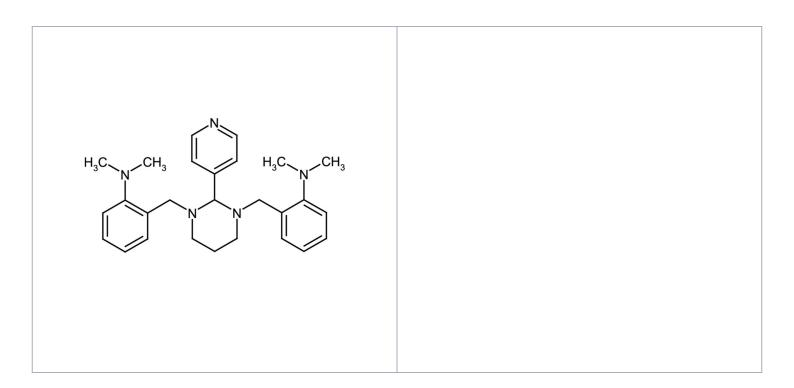
Order Online »

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)

Soluble in DMSO (5mg/ml).

Last modified: May 29, 2024

