

Purmorphamine

Smoothened activator

Cell permeable purine compound. Induces osteoblast differentiation. Agonist of smoothened by activation of the Hedgehog (Hh) pathway. Also shown to induce transdifferentiation in pre-adipocytes and myoblasts. Binds to and activates 7-transmembrane receptor of Hh.

Citations: 14

[View Online »](#)

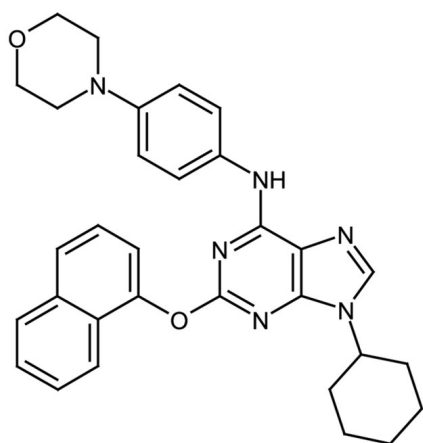
Ordering Information

[Order Online »](#)

ALX-420-045-M001	1mg
ALX-420-045-M005	5mg
ALX-420-045-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from light. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-(1-Naphthoxy)-6-(4-morpholinoanilino)-9-cyclohexylpurine
Appearance	Off-white to beige solid.
CAS	483367-10-8
Couple Target	Smoothened
Couple Type	Activator
Formula	$C_{31}H_{32}N_6O_2$
MW	520.6
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (10mM).

Last modified: April 24, 2025



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-be@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)