

Aminopurvalanol A

CDK/cyclin inhibitor

Cell permeable, reversible and competitive inhibitor of CDK1/cyclin B (IC_{50} =33nM), CDK2/cyclin A (IC_{50} =33nM), CDK2/cyclin E (IC_{50} =28nM) and CDK5/p35 (IC_{50} =20nM).

Citations: 3

[View Online »](#)

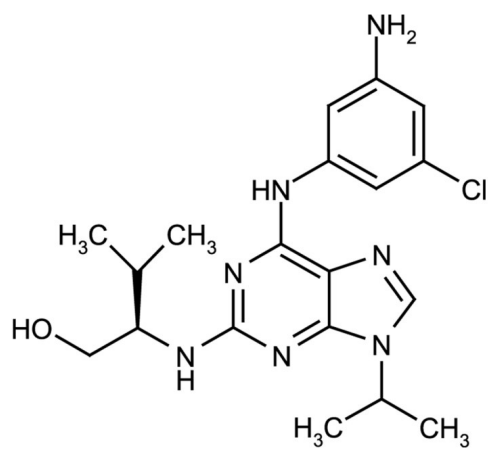
Ordering Information

[Order Online »](#)

ALX-270-249-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. Packaged under inert gas.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	(2R)-2-((6-((3-Amino-5-chlorophenyl)amino)-9-(1-methylethyl)-9H-purin-2-yl)amino)-3-methyl-1-butanol, NG-97
Appearance	White to brownish solid.
CAS	220792-57-4
Couple Target	CDK, Cyclin
Couple Type	Inhibitor
Formula	$C_{19}H_{26}N_7OCl$
Identity	Determined by NMR.
MW	403.9
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO or methanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com