

Purvalanol A

CDK/cyclin inhibitor

Inhibitor of human CDK1 ($IC_{50}=4nM$), CDK2/cyclin A ($IC_{50}=70nM$), Cdc2/cyclin B, CDK2/cyclin E ($IC_{50}=35nM$), CDK4/cyclin D1 ($IC_{50}=850nM$) as well as CDK5/p35 ($IC_{50}=75nM$). More membrane permeable than purvalanol B.

Citations: 5

[View Online »](#)

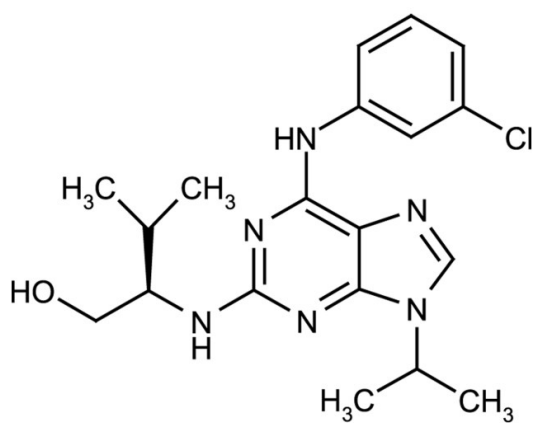
Ordering Information

[Order Online »](#)

ALX-270-246-M001	1mg
ALX-270-246-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Keep under inert gas. Keep cool and dry.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 6-[(3-Chloro)anilino]-2(1R)-(isopropyl-2-hydroxyethylamino)-9-isopropylpurine

Appearance Off-white to yellow powder.

CAS 212844-53-6

Couple Target CDK, Cyclin

Couple Type Inhibitor

Formula $C_{19}H_{25}N_6OCl$

MW 388.9

Purity ≥99% (NMR)

Solubility Soluble in DMSO, methanol or dichloromethane.

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-
be@enzolifesciences.com](mailto:info-be@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)