

# Purvalanol B

Cyclin-dependent kinase inhibitor.

Purvalanol B inhibits cyclin-dependent kinases with  $IC_{50}$  values of 6, 6, 9, > 10,000, and 6 nM for cdc2/cyclin B, cdk2/cyclin A, cdk2/cyclin E, cdk4/cyclin D1 and cdk5-p35 respectively. Has also been shown to have anti-proliferative properties, mediated by p42/p44 MAPK.

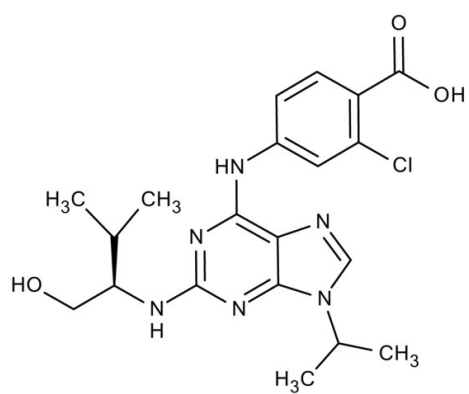
## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-270-247-M001 | 1mg |
|------------------|-----|

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Short Term Storage** +4°C

**Long Term Storage** -20°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Appearance** White powder.

**CAS** 212844-54-7

**Couple Target** CDK

**Couple Type** Inhibitor

**Formula**  $C_{20}H_{25}ClN_6O_3$

**Identity** Determined by EM-MS and NMR.

**MW** 432.9

**Purity** ≥98% (HPLC)

**Solubility** Soluble in DMSO (112mg/ml) or ethanol (4mg/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-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@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)