

CGP 74514A

CDK/cyclin inhibitor

Cell permeable, potent and selective inhibitor of CDK1/cyclin B (IC_{50} =25nM). Reported to affect the activities of other kinases only at much higher concentrations: PKC α (IC_{50} =6.1 μ M), PKA (IC_{50} =125 μ M) and EGFR (IC_{50} >10 μ M). Induces mitochondrial damage and apoptosis in several human leukemia cell lines (>3 μ M).

Citations: 2

[View Online »](#)

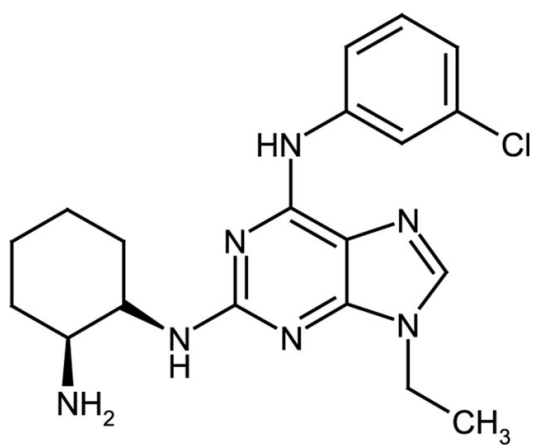
Ordering Information

[Order Online »](#)

ALX-270-391-M005	5mg
------------------	-----

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	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(cis-2-Aminocyclohexyl)-N-(3-chlorophenyl)-9-ethyl-9H-purine-2,6-diamine
Appearance	White to yellow solid.
Couple Target	CDK, EGFR, PKA, PKC
Couple Type	Inhibitor
Formula	$C_{19}H_{24}ClN_7$
MW	385.9
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

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-fr@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