

CVT-313

CDK/cyclin inhibitor

Specific inhibitor of CDK2/cyclin A (IC_{50} =500nM) and CDK2/cyclin E (IC_{50} =500nM). Also inhibits CDK1/cyclin B (IC_{50} =4.2 μ M) and CDK4/cyclin D1 (IC_{50} =215 μ M).

Citations: 3

[View Online »](#)

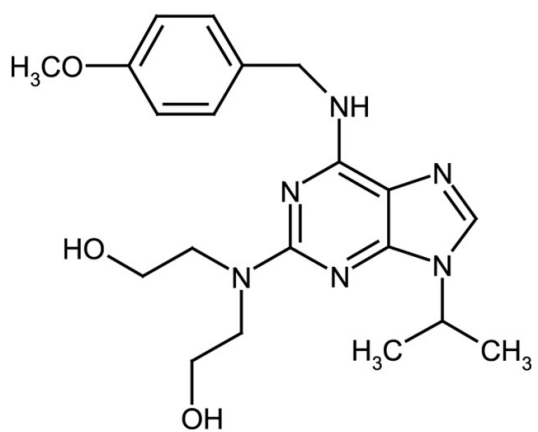
Ordering Information

[Order Online »](#)

ALX-270-393-M001	1mg
ALX-270-393-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 moisture.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 6-(4-Methoxybenzylamino)-2-[bis(2-hydroxyethylamino)]-9-isopropylpurine

Appearance White crystalline solid.

CAS 199986-75-9

Couple Target CDK, Cyclin

Couple Type Inhibitor

Formula $C_{20}H_{28}N_6O_3$

MW 400.5

Purity ≥98% (HPLC)

Solubility Soluble in DMSO, dimethyl formamide or 100% ethanol.

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)