

(S)-Roscovitine

CDK1 inhibitor

(S)-Enantiomer of the CDK inhibitor roscovitine (Prod. No. BML-CC205).
Potent inhibitor of CDK1/cyclin B kinase (IC_{50} =950nM).

Citations: 6

[View Online »](#)

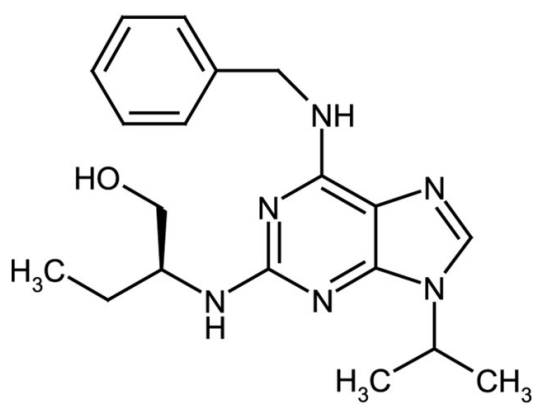
Ordering Information

[Order Online »](#)

ALX-350-293-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 6-Benzylamino-2-(S)-[(1-Ethyl)-2-hydroxyethylamino]-9-isopropylpurine

Appearance White to off-white solid.

CAS 186692-45-5

Couple Target CDK

Couple Type Inhibitor

Formula $C_{19}H_{26}N_6O$

MW 354.5

Purity ≥98% (HPLC)

Solubility Soluble in DMSO, dimethyl formamide, 100% ethanol 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-be@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