

(R)-Roscovitine

CDK inhibitor

Potent and selective inhibitor of CDKs. Selective for CDK1/cyclin B kinase (IC_{50} =450nM) [2], CDK2 (IC_{50} =700nM) [4] and CDK5/p35 (IC_{50} =160nM) [4,5]. More potent than olomoucine (Prod. No. BML-CC200). Inhibits M phase promoting factor (MPF) kinase activity. Has effects on calcium channel gating.

Citations: 22

[View Online »](#)

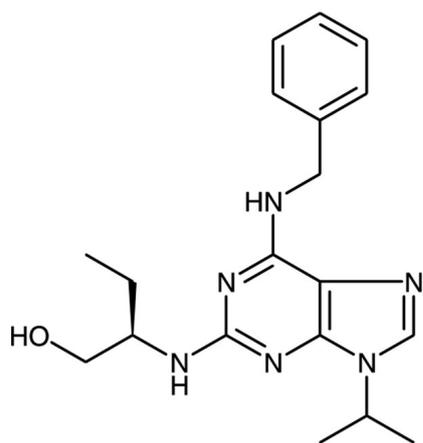
Ordering Information

[Order Online »](#)

BML-CC205-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	6-Benzylamino-2-(R)-[(1-ethyl)-2-hydroxyethylamino]-9-isopropylpurine, 2(R)-(1-Ethyl-2-hydroxyethylamino)-6-benzylamino-9-isopropylpurine, Seliciclib
Appearance	White to off-white solid.
CAS	186692-46-6
Couple Target	CDK
Couple Type	Inhibitor
Formula	$C_{19}H_{26}N_6O$
MW	354.4
Purity	≥95% (HPLC, TLC)
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-be@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