

Olomoucine II

CDK/cyclin inhibitor

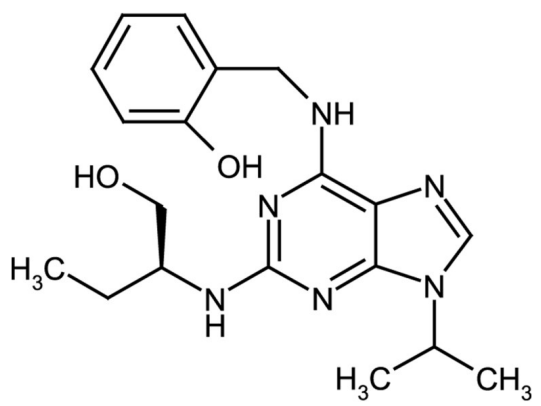
Selective inhibitor of CDK1/cyclin B (IC_{50} =20nM). Potent cytotoxin in tumor cells: IC_{50} =5.3 μ M against MCF7 cells, IC_{50} =11.1 μ M against K562 cells, IC_{50} =3.0 μ M against CEM cells, IC_{50} =6.3 μ M against HOS cells, and IC_{50} =6.3 μ M against G361 cells.

Citations: 2

[View Online »](#)

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-(2-Hydroxybenzylamino)-2(R)-[[1-hydroxymethyl)propyl]amino]-9-isopropylpurine

Appearance White to off-white crystalline solid.

CAS 500735-47-7

Couple Target CDK, Cyclin

Couple Type Inhibitor

Formula $C_{19}H_{26}N_6O_2$

MW 370.5

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

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-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-uk@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com