Olomoucine II

CDK/cyclin inhibitor

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

Citations: 2

View Online »

Ordering Information

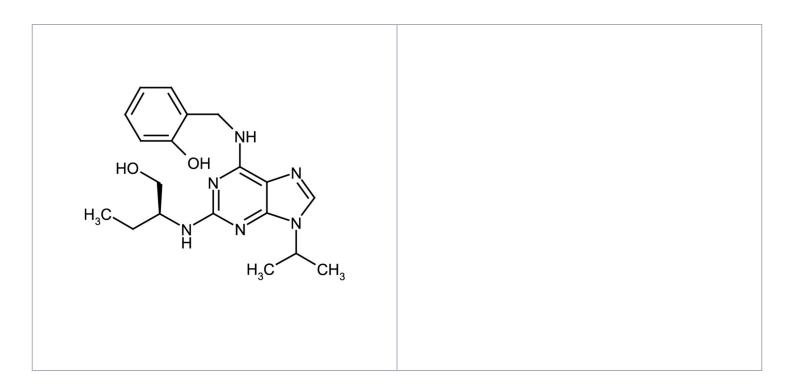
Order Online »

ALX-270-396-M005

5mg

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