

LY 294002

PI3 kinase inhibitor

Potent, cell permeable, inhibitor of PI(3)K (phosphoinositide 3-kinase) that acts on the ATP-binding site of the enzyme. Displays an IC_{50} of 1.4 μ M and inhibits all isoforms equally in in vitro assays using purified PI 3-kinase (p110 α IC_{50} =0.63 μ M). Inhibits Pim-1 kinase with IC_{50} =4.0 μ M. Other kinases such as PKC, PKA, MAP kinase, S6 kinase, EGF receptor tyrosine kinase, c-src kinase, PI 4-kinase, diacylglycerol kinase or rabbit kidney ATPase are not inhibited at 50 μ M.

Induces apoptosis in many cell types by blocking the PI3-K/Akt anti-apoptotic pathway. Useful tool for identifying cellular events which are regulated by the PI3-K/Akt axis. Sensitizes tumor cells to drug-induced apoptosis independent of its PI3K inhibitory activity. Inhibits proliferation and induces apoptosis in human colon cancer cells. Suppresses proliferation of mouse embryonic stem cells. Furthermore, inhibits autophagic sequestration in rat hepatocytes.

Citations: 51

[View Online »](#)

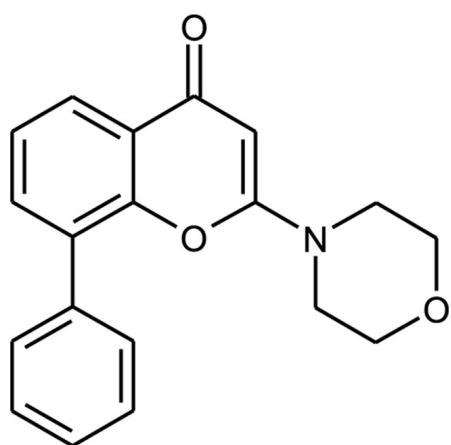
Ordering Information

[Order Online »](#)

BML-ST420-0005	5mg
BML-ST420-0025	25mg

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 at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-(4-Morpholinyl)-8-phenyl-4H-1-benzopyran-4-one
Appearance	White to pale yellow solid.
CAS	154447-36-6
Couple Target	PI3 kinase, Pim
Couple Type	Inhibitor
Formula	$C_{19}H_{17}NO_3$
MW	307.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (25mg/ml, warm), 100% ethanol (25mg/ml, warm) or methanol.

Technical Info / Product Notes Replacement for **ADI-HPK-106**.

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-
fr@enzolifesciences.com](mailto:info-fr@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)