

PI-103

Inhibitor of DNA-PK, PI3 kinase and mTOR

Potent, cell permeable and ATP-competitive inhibitor of DNA-PK, phosphoinositide 3-kinase (PI(3)K), the rapamycin-sensitive (mTORC1) and rapamycin-insensitive (mTORC2) complexes of the protein kinase mTOR. Exhibits little activity towards a panel of more than 40 other kinases even at concentrations as high as 10 μ M. Effectively blocks cell proliferation in glioma cell lines both *in vitro* and *in vivo*.

Citations: 7

[View Online »](#)

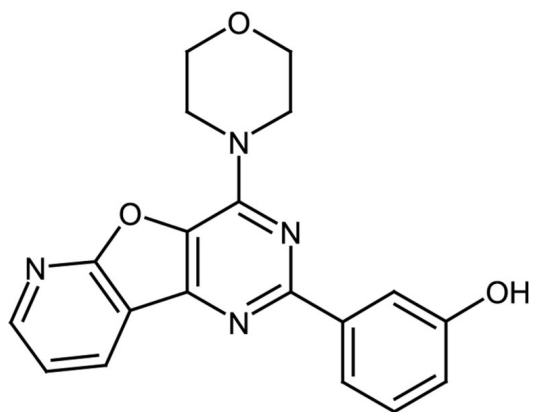
Ordering Information

[Order Online »](#)

ALX-270-460-M001	1mg
ALX-270-460-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from light. After reconstitution, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-(4-(4-Morpholinyl)pyrido[3',2':4,5]furo[3,2-d]pyrimidin-2-yl)phenol
Appearance	Off-white solid.
CAS	371935-74-9
Couple Target	DNA-PK, mTOR, PI3 kinase
Couple Type	Inhibitor
Formula	$C_{19}H_{16}N_4O_3$
Identity	Identity determined by MS and NMR.
MW	348.4
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (5mg/ml; warm).



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-be@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)