

# 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 $\mu$ M. Effectively blocks cell proliferation in glioma cell lines both *in vitro* and *in vivo*.

Citations: 7

[View Online »](#)

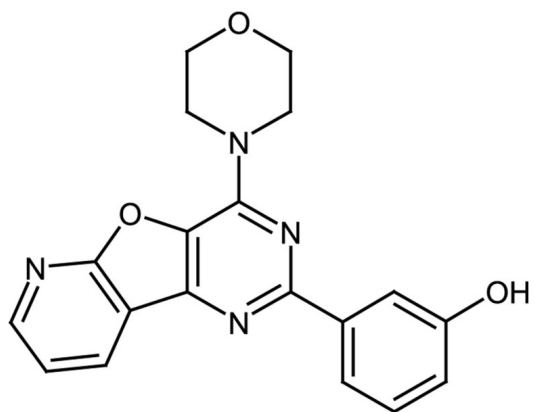
### Ordering Information

[Order Online »](#)

ALX-270-460-M001	1mg
------------------	-----

### Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	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.
<b>Handling</b>	Protect from light. After reconstitution, prepare aliquots and store at -20°C.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

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

**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)