

ZSTK474

Inhibits all four PI3K isoforms.

Cell permeable, reversible, potent and specific inhibitor of phosphoinositide 3-kinase (IC_{50} =17nM for PI(3)K p110 β , IC_{50} =53nM for PI(3)K p110 γ and IC_{50} =6nM for PI(3)K p110 δ). Blocks cellular PI(3)K/Akt (PKB) signalling *in vitro* and inhibits Akt(PKB)-dependent tumor growth in mice *in vivo* without significant toxic effect. More effective than LY 294,002 (Prod. No. BML-ST420) *in vitro* and as *in vivo*.

Citations: 4

[View Online »](#)

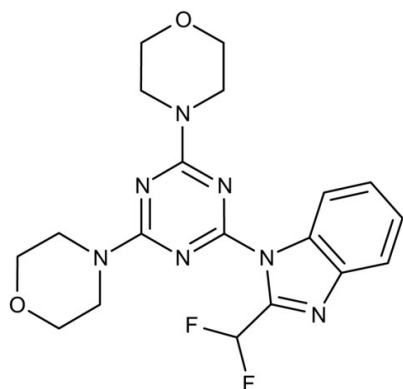
Ordering Information

[Order Online »](#)

ALX-270-454-M010	10mg
------------------	------

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 6 months when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2-(2-Difluoromethylbenzimidazol-1-yl)-4,6-dimorpholino-1,3,5-triazine
Appearance	White powder.
CAS	475110-96-4
Couple Target	PI3 kinase
Couple Type	Inhibitor
Formula	$C_{19}H_{21}F_2N_7O_2$
Identity	Determined by EM-MS and NMR.
MW	417.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (21mg/ml). Sparingly soluble in ethanol (<1mg/ml).

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)