

ALK5 Inhibitor I

Cell permeable, potent, selective and ATP-competitive inhibitor of TGF- β RI kinase (IC_{50} =51nM). Displays ~15-fold greater selectivity over p38 α MAPK (IC_{50} =740nM). Inhibits TGF- β -dependent cellular growth (IC_{50} =89nM in NIH 3T3 mouse fibroblasts) and transcription activation (IC_{50} =47nM in mink lung cells). Suppresses invasion of MDA-MB-231 breast cancer cells in a matrigel invasion assay.

Citations: 5

[View Online »](#)

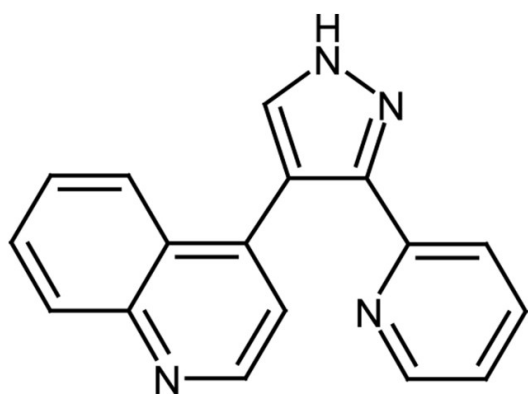
Ordering Information

[Order Online »](#)

ALX-270-448-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 2 years after receipt when stored at -20°C. Stock solutions are stable for up to 3 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	TGF-β RI Kinase Inhibitor I, 3-(Pyridin-2-yl)-4-(4-quinonyl)]-1H-pyrazole
Appearance	Yellow to brown solid.
CAS	396129-53-6
Couple Target	TGF-beta receptor
Couple Type	Inhibitor
Formula	$C_{17}H_{12}N_4$
MW	272.3
Purity	≥97% (TLC)
Solubility	Soluble in DMSO or methanol.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com