

ALK5 Inhibitor II

TGF- β type I receptor kinase (ALK5) inhibitor

ALK5 inhibitor II is a cell permeable, potent, selective and ATP-competitive inhibitor of TGF- β RI kinase (IC_{50} =23nM, 4nM and 18nM for binding, auto-phosphorylation, and cellular assay in HepG2 cells of TGF- β RI kinase, respectively). It minimally affects a panel of nine closely related kinases including p38 MAPK and GSK3 (IC_{50} >16 μ M). It can be used for the induction of the pluripotency of stem cells by replacing the transcription factor SOX2 via the inhibition of TGF- β signaling and inducing the expression of the transcription factor Nanog.

Citations: 50

[View Online »](#)

Ordering Information

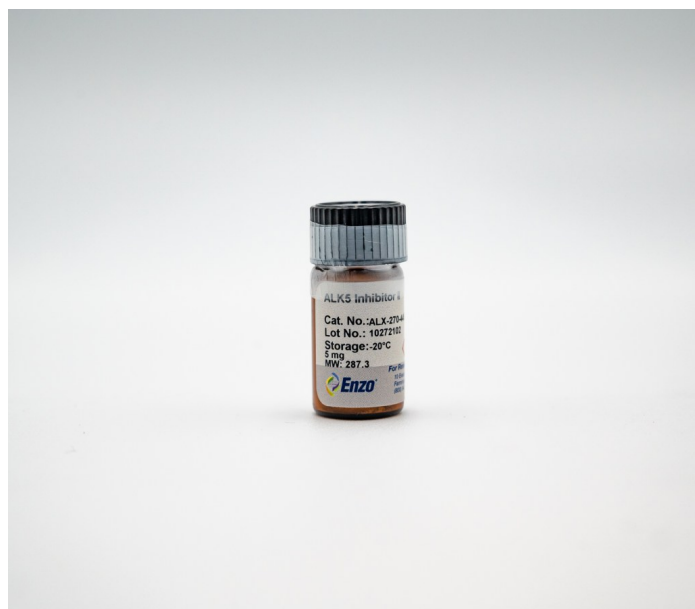
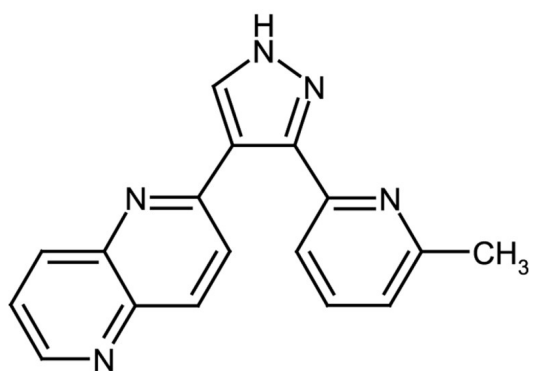
[Order Online »](#)

ALX-270-445-M001	1mg
ALX-270-445-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

- Potent ATP-competitive inhibitor of the TGF- β type I receptor ALK5
- Useful for cellular reprogramming and the culture of human pluripotent stem cells
- Active *in vitro* and *in vivo*
- GMP format available



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 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	TGF- β RI Kinase Inhibitor II, 2-(3-(6-Methylpyridin-2-yl)-1H-pyrazol-4-yl)-1,5-naphthyridine, RepSox, E-616452, SJN 2511
Appearance	White to off-white or yellow solid.
CAS	446859-33-2
Couple Target	TGF-beta receptor
Couple Type	Inhibitor
Formula	C ₁₇ H ₁₃ N ₅
Identity	Identity determined by NMR.
MW	287.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (5 mg/ml) or methanol (5 mg/ml).



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-be@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