

# ALK5 Inhibitor II

## TGF- $\beta$ type I receptor kinase (ALK5) inhibitor

ALK5 inhibitor II is a cell permeable, potent, selective and ATP-competitive inhibitor of TGF- $\beta$  RI kinase ( $IC_{50}$ =23nM, 4nM and 18nM for binding, auto-phosphorylation, and cellular assay in HepG2 cells of TGF- $\beta$  RI kinase, respectively). It minimally affects a panel of nine closely related kinases including p38 MAPK and GSK3 ( $IC_{50}$ >16 $\mu$ 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- $\beta$  signaling and inducing the expression of the transcription factor Nanog.

Citations: 52

[View Online »](#)

## Ordering Information

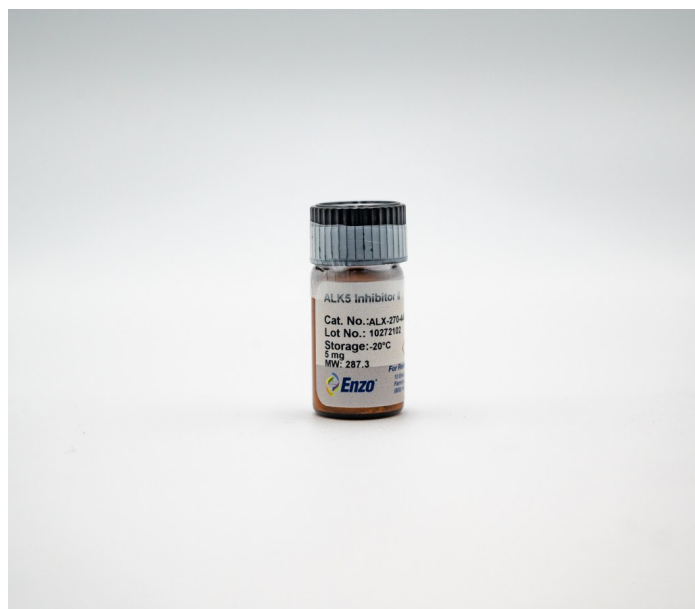
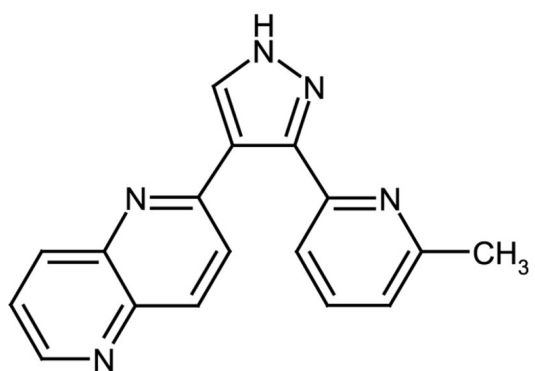
[Order Online »](#)

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

## Manuals, SDS & CofA

[View Online »](#)

- Potent ATP-competitive inhibitor of the TGF- $\beta$  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 <sub>17</sub> H <sub>13</sub> N <sub>5</sub>
Identity	Identity determined by NMR.
MW	287.3
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (5 mg/ml) or methanol (5 mg/ml).

Last modified: February 18, 2025



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-  
be@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)