

SB-525334

ALK5 inhibitor

SB-525334 is a potent inhibitor of the transforming growth factor β 1 receptor (ALK5). It has an IC50 of 14.3 nM for ALK5.

Key features and applications include:

- **High Selectivity:** Specifically inhibits ALK5 with minimal activity against other ALK receptors.
- **Potency:** Demonstrates strong inhibitory effects with an IC50 of 14.3 nM.
- **Block Phosphorylation and Nuclear Translocation of Smad2/3:** This inhibition reduces PAI-1 and procollagen α 1 (I) mRNA expression in renal epithelial carcinoma cells.
- **Reduce Fibrosis:** Effective in models of pulmonary fibrosis and renal disease by decreasing TGF- β -mediated responses.
- **Attenuate Pulmonary Arterial Hypertension:** Lowers pulmonary arterial pressure and right ventricular hypertrophy in relevant models.

Relevant disease states include:

- **Pulmonary Fibrosis:** Reduces myofibroblast proliferation and fibrosis.
- **Renal Disease:** Decreases markers of fibrosis in renal epithelial carcinoma cells.
- **Pulmonary Arterial Hypertension:** Mitigates symptoms by reducing TGF- β sensitivity.

- GMP format available

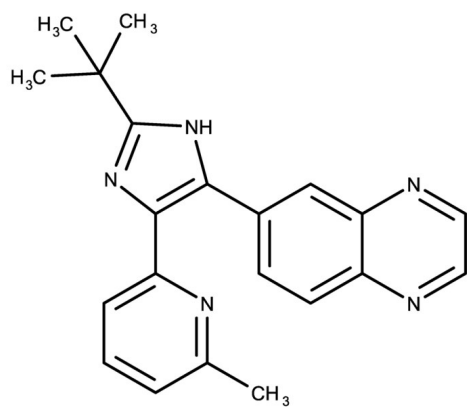
Ordering Information

[Order Online »](#)

ENZ-CHM492-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20°C for up to 3 months.
Handling	Tightly closed. Dry. Keep in a well-ventilated place. Keep locked up or in an area accessible only to qualified or authorized persons
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	6-[2-(1,1-Dimethylethyl)-5-(6-methyl-2-pyridinyl)-1H-imidazol-4-yl]quinoxaline
Appearance	Yellow solid.
CAS	356559-20-1
Couple Target	Activin receptor, TGF-beta receptor
Couple Type	Inhibitor
Formula	$C_{21}H_{21}N_5$
Identity	Determined by NMR.
MW	343.42
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (up to 30 mg/ml).



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)