

LY-364947

ALK5 Inhibitor

Catalog # T07-901

Lot # D3370-8

CAS # 396129-53-6

Product Description

Molecular Formula: C₁₇H₁₂N₄

Appearance: Brown Powder

Molecular Weight: 272.31

Purity: >98% (TLC); NMR (Conforms)

Solubilization: May be dissolved in DMSO (25 mg/ml)

Alias

4-(3-Pyridin-2-yl-1H-pyrazol-4-yl)quinolone; HTS-466284

Specific Activity

ALK5 kinase inhibitor

Storage and Stability

Store desiccated as supplied at ambient temperature for up to 2 year. Store solutions at -20°C for up to 3 months.

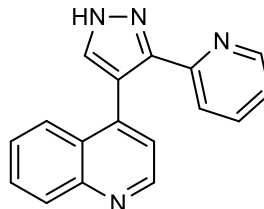
Scientific Background

Selective ALK5 inhibitor, IC₅₀=59, 400 and 1400 nM for TGF-β RI, TGF-β RII and MLK-7K, respectively¹. Inhibits Smad2 phosphorylation induced by TGF-β as well as fibronectin expression and MDA-MB-231 cell invasion².

References

1. JS Sawyer *et al.* *J. Med. Chem.* 2003 46:3953
2. SR Shiou *et al.* *J. Biol. Chem.* 2006 281:33971
3. ME Hardee *et al.* *Cancer Res.* 2012 72:4119

Molecular Structure



LY-364947

ALK5 kinase inhibitor

Catalog Number	T07-901
CAS Number	396129-53-6
Specific Lot Number	D3370-8
Purity	>98% (TLC); NMR (Conforms)
Format	Lyophilized brown powder
Solubilization	May be dissolved in DMSO (25 mg/ml)
Storage & Shipping	Store desiccated as supplied at ambient temperature for up to 2 year. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.