

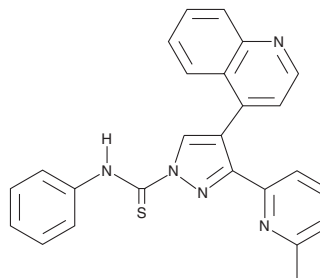
PRODUCT INFORMATION



A 83-01

Item No. 9001799

CAS Registry No.: 909910-43-6
Formal Name: 3-(6-methyl-2-pyridinyl)-N-phenyl-4-(4-quinolinyl)-1H-pyrazole-1-carbothioamide
MF: C₂₅H₁₉N₅S
FW: 421.5
Purity: ≥95%
UV/Vis.: λ_{max}: 227, 307 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

A 83-01 is supplied as a crystalline solid. A stock solution may be made by dissolving the A 83-01 in the solvent of choice. A 83-01 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of A 83-01 in these solvents is approximately 0.3, 14, and 25 mg/ml, respectively.

A 83-01 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, A 83-01 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. A 83-01 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

A 83-01 is an inhibitor of TGF-β type I receptor kinase (ALK5), activin type IB receptor (ALK4), and nodal type I receptor (ALK7) (IC₅₀s = 12, 45, and 7.5 nM, respectively).¹ It blocks the phosphorylation of SMAD2/3 and inhibits TGF-β-induced epithelial-to-mesenchymal transition.¹ It has little effect on bone morphogenic type I receptors, p38 mitogen-activated protein kinase, or ERK.¹ A 83-01 has been used to reprogram fibroblasts into alternative lineages, including neural stem cells and cardiomyocytes.^{2,3} This product is also available as part of Cayman's Hepatocyte Differentiation Reagent Kit (Item No. 44737).

References

1. Tojo, M., Hamashima, Y., Hanyu, A., *et al.* The ALK-5 inhibitor A-83-01 inhibits Smad signaling and epithelial-to-mesenchymal transition by transforming growth factor-β. *Cancer Sci.* **96(11)**, 791-800 (2005).
2. Cao, N., Huang, Y., Zheng, J., *et al.* Conversion of human fibroblasts into functional cardiomyocytes by small molecules. *Science* **Apr**, (2016).
3. Zhang, M., Lin, Y. H., Sun, Y. J., *et al.* Pharmacological reprogramming of fibroblasts into neural stem cells by signaling-directed transcriptional activation. *Cell Stem Cell* **Apr**, (2016).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/27/2025

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM