

3-(4-Pyridyl)-1H-indole

ROCK inhibitor

A cell permeable, selective, and ATP-competitive inhibitor of Rho kinase (ROCK) activity ($IC_{50}=25\mu M$). Does not inhibit the activation of Rho nor does it affect the *in vitro* activities of MLCK, PKC α , and SAPK2a/p38 α . Shown to be less potent than Y-27632 (Prod. No. ALX-270-333).

Citations: 2

[View Online »](#)

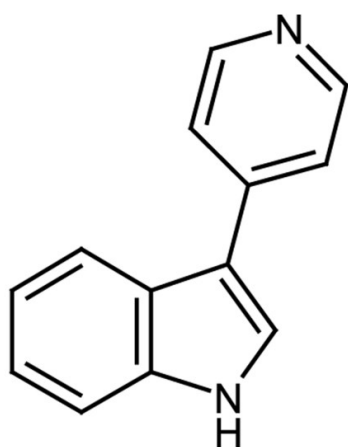
Ordering Information

[Order Online »](#)

ALX-270-429-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. 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

Appearance	Off-white to beige solid.
CAS	7272-84-6
Couple Target	ROCK
Couple Type	Inhibitor
Formula	$C_{13}H_{10}N_2$
Identity	Determined by NMR.
MW	194.2
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO or dimethyl formamide.

Last modified: May 29, 2024



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)