

H-1152 . dihydrochloride

ROCK inhibitor

A cell permeable, highly specific, potent and ATP-competitive inhibitor of Rho kinase (ROCK) ($K_i=1.6\text{nM}$). Exhibits a much weaker affinity for other serine/threonine kinases ($K_i=630\text{nM}$ for PKA, 9.27mM for PKC and 10.1mM for MLCK). More potent and selective than Y-27632 (Prod. No. ALX-270-333). Inhibits the phosphorylation of MARCKS in cells stimulated by lysophosphatidic acid. Inhibits EP3-stimulated NO formation. Prevents fragmentation of apoptotic cells. Relieves neuropathic pain.

Citations: 18

[View Online »](#)

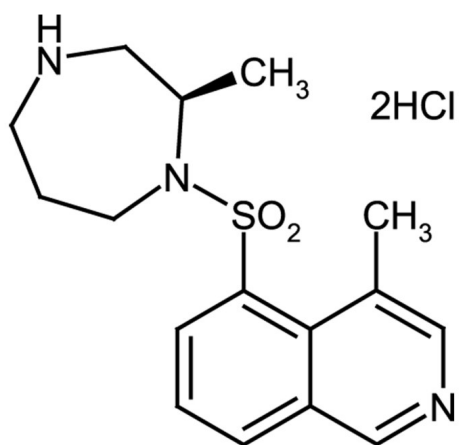
Ordering Information

[Order Online »](#)

ALX-270-423-M001	1mg
ALX-270-423-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	H-1152P . 2HCl, (S)-(+)-2-Methyl-1-[(4-methyl-5-isoquinolinyl)sulfonyl]homopiperazine . 2HCl
Appearance	White to off-white solid.
CAS	451462-58-1
Couple Target	ROCK
Couple Type	Inhibitor
Formula	$C_{16}H_{21}N_3O_2S \cdot 2HCl$
Identity	Identity determined by 1H-NMR and 13C-NMR.
MW	319.4 . 73.0
Purity	≥98% (TLC)
Solubility	Soluble in water.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com