

HA-100 . dihydrochloride

Protein kinase inhibitor

Inhibitor of myosin light chain kinase, cAMP-dependent protein kinase (PKA) and protein kinase C (PKC).

Citations: 2

[View Online »](#)

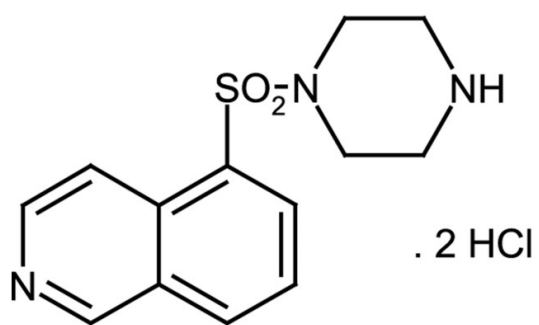
Ordering Information

[Order Online »](#)

ALX-270-069-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 1-(5-Isoquinolinylsulfonyl)piperazine . 2HCl

Appearance Off-white lyophilized solid.

CAS 84468-24-6

Couple Target MLCK, PKA, PKC

Couple Type Inhibitor

Formula $C_{13}H_{15}N_3O_2S \cdot 2HCl$

MW 350.3

Purity ≥98% (HPLC)

Solubility Soluble in water or DMSO.

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-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)