

# HA-1077

## Rho kinase (ROCK) inhibitor

Inhibits Rho kinase (ROCK;  $K_i=0.33\mu\text{M}$ ). Selective protein kinase A and G inhibitor (both  $K_i$ 's =  $1.6\mu\text{M}$ ). Inhibits agonist-induced activation of NADPH oxidase in human neutrophils ( $\text{IC}_{50}=15\mu\text{M}$ ). Inhibitor of myosin light chain kinase and  $\text{Ca}^{2+}$ /calmodulin-dependent protein kinase II. Inhibits translocation of  $\text{PKC}\beta\text{I}$ ,  $\text{PKC}\beta\text{II}$  and  $\text{PKC}\zeta$ . Cell permeable  $\text{Ca}^{2+}$  antagonist with antivasospastic properties.

Citations: 14

[View Online »](#)

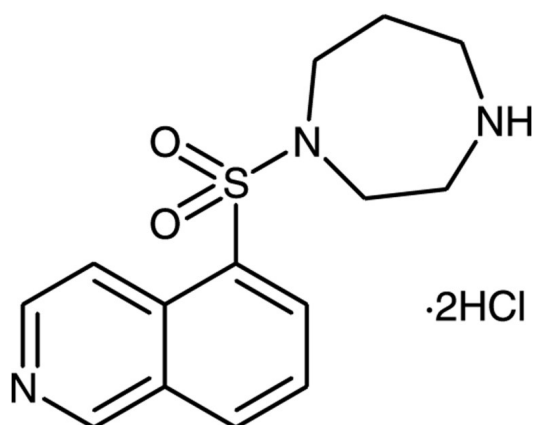
## Ordering Information

[Order Online »](#)

BML-EI233-0010	10mg
BML-EI233-0050	50mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at 0-4°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	1-(5-Isoquinolinesulfonyl)homopiperazine . 2HCl, Fasudil
Appearance	White to off-white solid.
CAS	103745-39-7
Couple Target	CaM kinase, MLCK, PKA, PKG, ROCK
Couple Type	Inhibitor
Formula	$C_{14}H_{17}N_3O_2S \cdot 2HCl$
Identity	Determined by <sup>1</sup> H-NMR and MS.
MI	14: 3942
MW	364.3
MeltingPoint	238-242°C
Purity	≥98% (TLC)
Solubility	Soluble in water (>20mg/ml), methanol or DMSO.



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-  
be@enzolifesciences.com](mailto:info-be@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)