

H-7

PKC inhibitor

Inhibits PKA ($K_i=3.0\mu\text{M}$), PKG ($K_i=5.8\mu\text{M}$), MLCK ($K_i=97\mu\text{M}$) and PKC ($K_i=6\mu\text{M}$). Also inhibits myosin light chain kinase. Induces apoptotic DNA fragmentation and cell death. Inhibits telomerase activity in treated NPC-076 cells. Cell permeable.

Citations: 11

[View Online »](#)

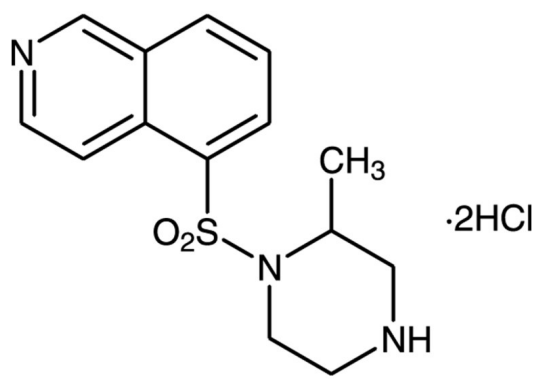
Ordering Information

[Order Online »](#)

BML-EI148-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at 0-4°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from light.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(±)-1-(5-Isoquinolinesulfonyl)-2-methylpiperazine . 2HCl
Appearance	White to light yellow solid.
CAS	108930-17-2
Couple Target	MLCK, PKA, PKC, PKG, Telomerase
Couple Type	Inhibitor
Formula	$C_{14}H_{17}N_3O_2S \cdot 2 HCl$
MW	364.3
Purity	≥98% (TLC)
Solubility	Soluble in water (50mg/ml), 100% ethanol 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-
be@enzolifesciences.com](mailto:info-be@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)