

H-89 . 2HCl

Protein kinase inhibitor

Cell permeable potent and selective inhibitor of cAMP- and cGMP-dependent protein kinases (PKA, $K_i=48\text{nM}$, and PKG) and protein kinase $C\mu$ (PKC μ). In contrast, most other protein kinase C (PKC) isotypes are much more weakly inhibited. Also inhibits Ca^{2+} /calmodulin-dependent protein kinase II, casein kinase I and myosin light chain kinase. Induces apoptosis. Concentration range for cultured cells 30-100 μM : osteoblastic cells, renal proximal tubule cells and *Xenopus* oocytes.

Citations: 24

[View Online »](#)

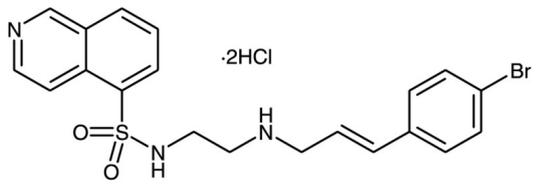
Ordering Information

[Order Online »](#)

BML-EI196-0005	5mg
BML-EI196-0025	25mg

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 4 months in DMSO.
Long Term Storage	+4°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	N-[2-(p-Bromocinnamylamino)ethyl]-5-isoquinolinesulfonamide . 2HCl
Appearance	White to off-white solid.
CAS	127243-85-0
Couple Target	CaM kinase, Casein kinase, MLCK, PKA, PKD, PKG
Couple Type	Inhibitor
Formula	$C_{20}H_{20}BrN_3O_2S \cdot 2HCl$
MW	519.3
MeltingPoint	141-143°C
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (25mg/ml), methanol and 50% ethanol/water (20mg/ml).
Technical Info / Product Notes	Replacement for ADI-HPK-105

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