

H-9 . HCl

PKA inhibitor

An excellent affinity ligand for chromatographic purification of protein kinases. Inhibits protein kinases (K_i 's: PKA=1.9 μ M; PKC=18 μ M; PKG=0.87 μ M; MLCK=70 μ M).

Citations: 5

[View Online »](#)

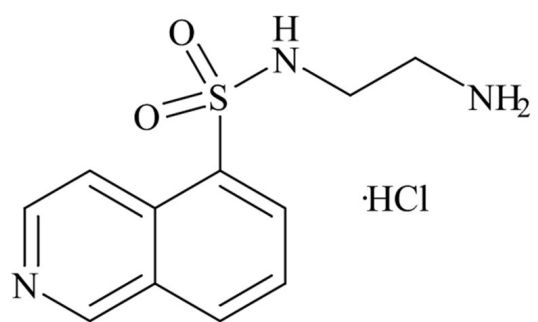
Ordering Information

[Order Online »](#)

BML-EI195-0010	10mg
BML-EI195-0050	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	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(2-Aminoethyl)-5-isoquinolinesulfonamide . HCl
Appearance	White solid.
CAS	116970-50-4
Couple Target	PKA, PKG
Couple Type	Inhibitor
Formula	$C_{11}H_{13}N_3O_2S \cdot HCl$
Identity	Determined by ¹ H-NMR.
MW	287.8
Purity	≥98% (TLC)
Solubility	Soluble in water.

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)