

A-3

PKA inhibitor

Inhibitor of cAMP- and cGMP-dependent protein kinase (PKA, $K_i=4.3\mu\text{M}$ and PKG, $K_i=3.8\mu\text{M}$), protein kinase C (PKC, $K_i=47\mu\text{M}$), casein kinase I and II, and myosin light chain kinase (MLCK, $K_i=7.4\mu\text{M}$).

Citations: 2

[View Online »](#)

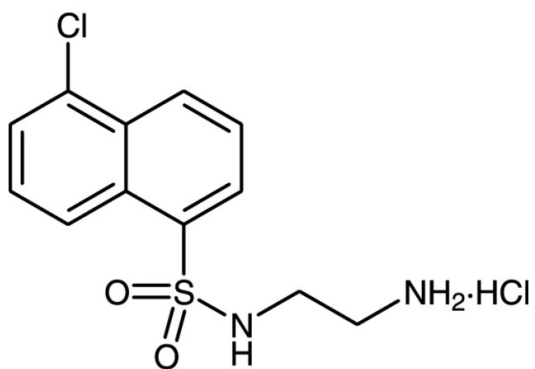
Ordering Information

[Order Online »](#)

BML-EI145-0010	10mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white solid.

CAS 78957-85-4

Couple Target Casein kinase, MLCK, PKA, PKC, PKG

Couple Type Inhibitor

Formula $C_{12}H_{13}ClN_2O_2S \cdot HCl$

MW 284.8 . 36.5

Purity ≥98% (TLC)

Solubility Soluble in water (50mg/ml), DMSO (50mg/ml) or 100% ethanol.

Technical Info / Product Notes Note: Product is not sterile.

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)