

W-7

Calmodulin inhibitor

Binds to calmodulin, inhibiting Ca^{2+} -calmodulin-regulated enzyme activity. It inhibits the Ca^{2+} /calmodulin-induced activation of myosin light chain kinase in a concentration dependent manner ($\text{IC}_{50}=25\text{ }\mu\text{M}$). Cell permeable.

Citations: 6

[View Online »](#)

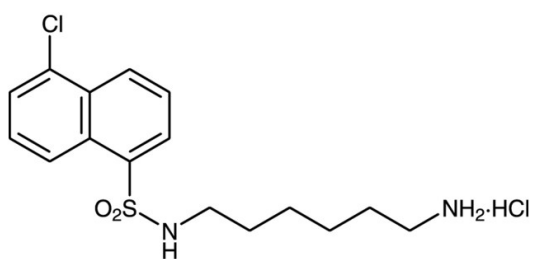
Ordering Information

[Order Online »](#)

BML-CA320-0050	50mg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name N-(6-Aminohexyl)-5-chloro-1-naphtalene-sulfonamide . HCl

Appearance White to off-white solid.

CAS 61714-27-0

Couple Target Calmodulin, MLCK

Couple Type Inhibitor

Formula $C_{16}H_{21}ClN_2O_2S \cdot HCl$

MW 377.3

Purity ≥98% (TLC)

Solubility Soluble in water (warm), methanol, 100% ethanol (10mg/ml) (warm), DMSO (25mg/ml), dimethyl formamide or ethylene glycol.

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-
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)