

W-5

Calmodulin inhibitor

Binds to calmodulin, inhibiting Ca^{2+} -calmodulin-regulated enzyme activity.
Cell permeable.

Citations: 3

[View Online »](#)

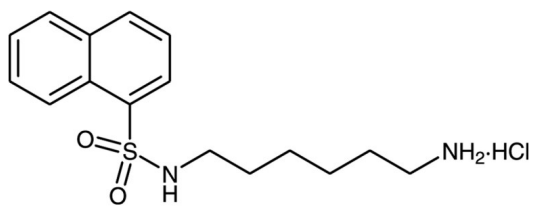
Ordering Information

[Order Online »](#)

BML-CA315-0020	20mg
----------------	------

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 at room temperature. Stock solutions are stable for up to 6 months at -20°C.
Handling	Protect from light.
Long Term Storage	Ambient
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(6-Aminohexyl)-1-naphthalenesulfonamide . HCl
Appearance	White to off-white solid.
CAS	61714-25-8
Couple Target	Calmodulin
Couple Type	Inhibitor
Formula	$C_{16}H_{22}N_2O_2S \cdot HCl$
MW	342.9
Purity	≥98% (HPLC)
Solubility	Soluble in water, DMSO (20mg/ml), 100% ethanol (hot) or methanol (hot).

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-fr@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