

KN-62

CaM kinase inhibitor

KN-62 is a potent and selective inhibitor of Ca^{2+} /calmodulin kinase II ($K_i = 0.9 \mu\text{M}$). Causes cells to arrest in S phase. In HEK-293 cells expressing type III adenylyl cyclase (III-AC), $10 \mu\text{M}$ blocked CaMKII phosphorylation of III-AC and inhibited hormone-stimulated Ca^{2+} oscillations.

Citations: 13

[View Online »](#)

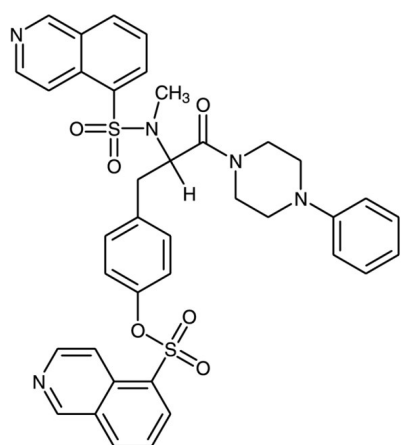
Ordering Information

[Order Online »](#)

BML-EI230-0001	1mg
----------------	-----

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 yellow solid.

CAS 127191-97-3

Couple Target CaM kinase

Couple Type Inhibitor

Formula $C_{38}H_{35}N_5O_6S_2$

Identity Determined by 1H -NMR.

MW 721.8

Purity $\geq 98\%$ (TLC)

Solubility Soluble in methanol and DMSO (50mg/ml)

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)