

KN-93

CaM kinase inhibitor

KN-93 is a potent, selective and cell-permeable inhibitor of CaM kinase II (IC_{50} =370 nM). Does not significantly inhibit PKA. Causes cell-cycle arrest in NIH 3T3 fibroblasts and HeLa cells. Inhibits glucose-induced and forskolin-stimulated insulin release from isolated pancreatic islets of Langerhans. Induces G_1 cell cycle arrest and apoptosis in NIH 3T3 cells.

Citations: 9

[View Online »](#)

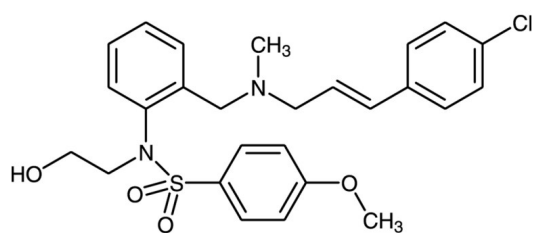
Ordering Information

[Order Online »](#)

BML-EI268-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 2 years after receipt. Stock solutions are stable for up to 3 months when stored at -20°C.
Handling	Protect from light. Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	white to off-white solid
CAS	139298-40-1
Couple Target	CaM kinase
Couple Type	Inhibitor
Formula	$C_{26}H_{29}ClN_2O_4S$
MW	501
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (25mg/ml).



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)