

KN-93 phosphate

CaM kinase II inhibitor

Selective inhibitor of rat brain Ca^{2+} /calmodulin-dependent protein kinase II. Inhibits glucose-induced and forskolin-stimulated insulin release from isolated pancreatic islets of Langerhans. Has no significant effect on protein kinase A (PKA) activity. Induces G_1 cell cycle arrest and apoptosis in NIH 3T3 cells.

Citations: 5

[View Online »](#)

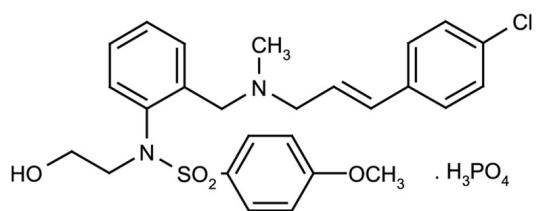
Ordering Information

[Order Online »](#)

ALX-430-127-M001	1mg
ALX-430-127-M005	5mg

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.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	KN-93 (water soluble), 2-[N-(2-Hydroxyethyl)-N-(4-methoxybenzenesulfonyl)]amino-N-(4-chlorocinnamyl)-N-methylbenzylamine . hydrogen phosphate
Appearance	White to off-white solid.
CAS	1188890-41-6
Couple Target	CaM kinase
Couple Type	Inhibitor
Formula	$C_{26}H_{29}ClN_2O_4S \cdot H_3PO_4$
Identity	Determined by NMR.
MW	501.0 . 98.0
Purity	≥98% (1H-NMR)
Solubility	Soluble in DMSO or water.



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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)