

7-Oxostaurosporine

Apoptosis inducer. PKC inhibitor.

Antibiotic. Potent inhibitor of protein kinase C (PKC). Inhibits the cell cycle at the G2 stage.

Citations: 2

[View Online »](#)

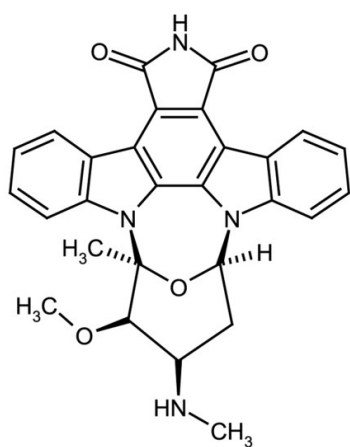
Ordering Information

[Order Online »](#)

ALX-380-210-M001	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

Alternative Name Antibiotic BMY 41950, Antibiotic RK 1409, LCM76-L

Appearance Yellow solid.

CAS 125035-83-8

Couple Target PKC

Couple Type Inhibitor

Formula $C_{28}H_{24}N_4O_4$

MW 480.5

Purity ≥98% (HPLC)

Solubility Soluble in DMSO, dimethyl formamide, methanol or 100% ethanol; slightly soluble in water.

Source Isolated from *Streptomyces* sp. MST-AS5345.

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-
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)