

Staurosporine

Alkaloid produced by *Streptomyces*, precursor in the synthesis of K252c; PKC and mammalian RNA splicing inhibitor.

A protein kinase C inhibitor that induces apoptosis in many cell types. Experimental results obtained in a cell-free assay suggest that cytoplasmic proteins directly modulated by staurosporine may be involved in signaling the induction of DNA fragmentation and apoptosis.

Ordering Information

[Order Online »](#)

LKT-S7600-M005	5mg
LKT-S7600-M025	25mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	AM-2282
CAS	62996-74-1
Couple Type	Inhibitor
Formula	$C_{28}H_{26}N_4O_3$
MW	466.5
Purity	≥98%
Solubility	Soluble in DMSO (50 mM), ethanol, or methanol (2 mg/mL). Insoluble in water.



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-be@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