

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-M001 | 1mg |
| LKT-S7600-M005 | 5mg |
| LKT-S7600-M025 | 25mg |
| LKT-S7600-M100 | 100mg |

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