

# UCN-01

## PKC and CDK inhibitor

Inhibitor of protein kinase C (PKC) and cyclin-dependent kinases (CDKs) resulting in accumulation of cells in the G1 phase and induction of apoptosis. Enhances the cytotoxicity of other anticancer drugs, such as DNA-damaging agents and anti-metabolite drugs, through putative abrogation of G2 and/or S phase accumulation induced by these agents.

Citations: 8

[View Online »](#)

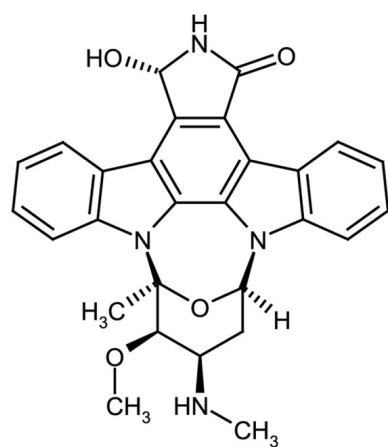
## Ordering Information

[Order Online »](#)

ALX-380-222-MC25	0.25mg
------------------	--------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	7-Hydroxystaurosporine
Appearance	White film.
CAS	112953-11-4
Couple Target	PKC
Couple Type	Inhibitor
Formula	$C_{28}H_{26}N_4O_4$
MW	482.5
Purity	≥99% (HPLC)
RTECS	KC6600010
Solubility	Soluble in methanol, 100% ethanol, dimethyl formamide or DMSO.
Source	Isolated from <i>Streptomyces</i> sp. MST-AS5345.
Technical Info / Product Notes	<b>Note:</b> Some interconversion of UCN-01 to UCN-02 may occur under acidic HPLC conditions).



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-  
be@enzolifesciences.com](mailto:info-be@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)