

Alsterpaullone

Inhibitor of CDK and GSK

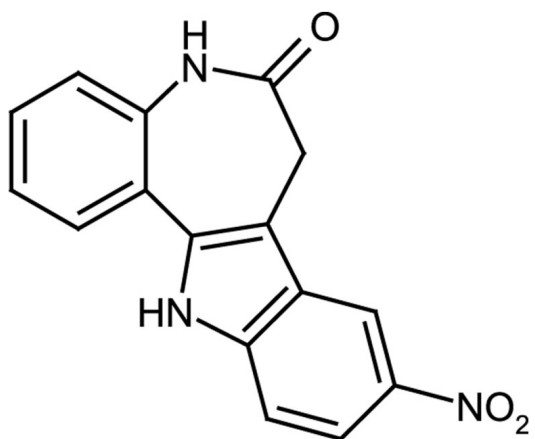
Potent inhibitor of CDK1/cyclin B (IC_{50} =35nM). Potent and selective inhibitor of CDK2/cyclin A, CDK2/cyclin E (IC_{50} =200nM), CDK5/p25 (IC_{50} =40nM), CDK5/p35 (IC_{50} =40nM) and GSK-3 β . Induces apoptosis by activation of caspase-8 and caspase-9 followed by disruption of the mitochondrial potential. *Licensed from NCI.*

Citations: 11

[View Online »](#)

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	+4°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow to brown powder.
CAS	237430-03-4
Couple Target	Caspase, Caspase-8, Caspase-9, CDK, GSK
Couple Type	Activator, Inhibitor
Formula	$C_{16}H_{11}N_3O_3$
Identity	Determined by IR, 1H-NMR and 13C-NMR.
MW	293.3
Purity	≥95%
Solubility	Soluble in DMSO; insoluble in water or 100% ethanol.

Last modified: May 29, 2024



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