

Kenpaullone

Inhibitor of CDK and GSK

Potent inhibitor of CDK1/cyclin B ($IC_{50}=400nM$). Also inhibits CDK2/cyclin A ($IC_{50}=680nM$), CDK5 ($IC_{50}=850nM$) and with much less effect other kinases. More recently, kenpaullone has been found to be a useful GSK-3 β inhibitor ($IC_{50}=23nM$). Induces pluripotency in somatic cells when used in combination with reprogramming factors. Increases neurogenesis of human neural progenitor cells through stimulation of Wnt/ β -catenin signaling pathway.

Citations: 10

[View Online »](#)

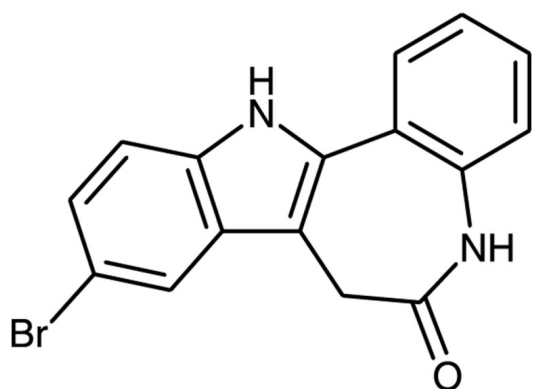
Ordering Information

[Order Online »](#)

BML-EI310-0005	5mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	9-Bromo-7,12-dihydroindolo-[3,2-d][1]benzazepin-6(5H)-one
Appearance	Tan solid.
CAS	142273-20-9
Couple Target	CDK, GSK
Couple Type	Inhibitor
Formula	$C_{16}H_{11}BrN_2O$
Identity	Determined by 1H -NMR.
MW	327.2
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (>25mg/ml); insoluble in water or ethanol.



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)