

# 1-Azakenpauillone

GSK-3 $\beta$  inhibitor

Analog of kenpauillone (Prod. No. BML-EI310) acting as a potent and ATP-competitive inhibitor of GSK-3 $\beta$  ( $IC_{50}$ =18nM). Displays ~100-200 fold greater selectivity on GSK-3 $\beta$  than over CDK1/cyclin B and CDK5/p25 ( $IC_{50}$ =2.0 $\mu$ M and 4.2 $\mu$ M, respectively).

Citations: 4

[View Online »](#)

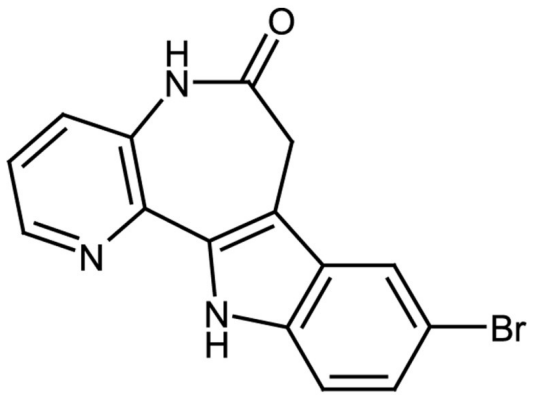
Ordering Information

[Order Online »](#)

ALX-270-430-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended. Stock solutions are stable for up to 6 months when stored at -20°C.
<b>Handling</b>	Protect from moisture. Protect from light.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	9-Bromo-7,12-dihydropyrido[3',2':2,3]azepino[4,5-b]indol-6(5H)-one, 1-AKP
<b>Appearance</b>	Off-white to brown powder.
<b>CAS</b>	676596-65-9
<b>Couple Target</b>	GSK
<b>Couple Type</b>	Inhibitor
<b>Formula</b>	$C_{15}H_{10}BrN_3O$
<b>MW</b>	328.2
<b>Purity</b>	≥95% (HPLC, TLC)
<b>Solubility</b>	Soluble in DMSO; insoluble in water.

Last modified: May 29, 2024



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)