

Indirubin-3'-monoxime

Inhibitor of CDK and GSK

A potent inhibitor of glycogen synthase kinase-3 β (IC_{50} =22 nM). Also inhibits CDK1 (p34^{cdc2}; IC_{50} =180 nM) and CDK5 (IC_{50} =100 nM). It reversibly arrests asynchronous HBL-100 cells at G2. Induces apoptosis in the mammary carcinoma cell line MCF-7 (10 μ M).

Citations: 6

[View Online »](#)

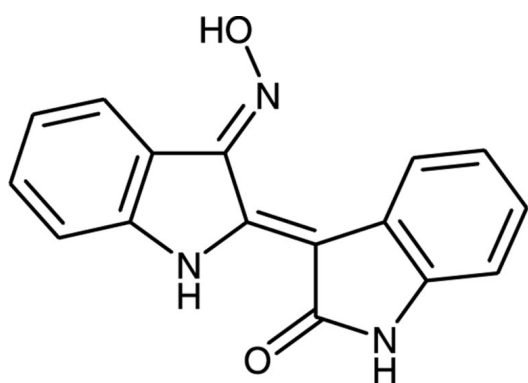
Ordering Information

[Order Online »](#)

BML-CC207-0005	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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Dark red solid.
CAS	160807-49-8
Couple Target	CDK, GSK
Couple Type	Inhibitor
Formula	$C_{16}H_{11}N_3O_2$
MW	277.3
Purity	≥98%
Solubility	Soluble in DMSO (6mg/ml) and ethanol (4mg/ml).

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-
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)