

Indirubin-5-sulfonic acid . sodium salt

Inhibitor of CDK and GSK

Inhibitor of cyclin-dependent kinases (CDKs), with selectivity for CDK1/cyclin B (IC_{50} =55nM), CDK2/cyclin A (IC_{50} =35nM), CDK2/cyclin E (IC_{50} =150nM), CDK4/cyclin D1 (IC_{50} =300nM), CDK5/p35 (IC_{50} =65nM) and GSK-3 β (IC_{50} =280nM).

Citations: 4

[View Online »](#)

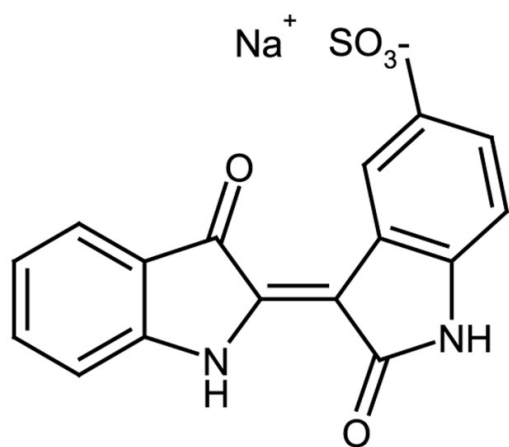
Ordering Information

[Order Online »](#)

ALX-270-296-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name E226, Indirubin-5-sulphonate . Na

Appearance Black powder.

Couple Target CDK, GSK

Couple Type Inhibitor

Formula $C_{16}H_9N_2O_5S \cdot Na$

Identity Determined by 1H -NMR.

MW 341.3 . 23.0

Purity $\geq 95\%$ (1H -NMR)

Solubility Soluble in DMSO or 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)