

SU 9516

CDK/cyclin inhibitor

Cell permeable, potent, selective and ATP-competitive inhibitor of CDK2/cyclin A (IC_{50} =22nM), CDK1/cyclin B (IC_{50} =40nM) and CDK4/cyclin D1 (IC_{50} =200nM). Has no effect on the activity of PKC, p38, PDGFR β and EGFR (IC_{50} >10 μ M). Displays anti-proliferative and pro-apoptotic properties in tumor cells. Inhibits proliferation of growth factor-stimulated colon carcinoma cells by binding to CDK2 and thus preventing the phosphorylation of pRb and its dissociation from E2F.

Citations: 4

[View Online »](#)

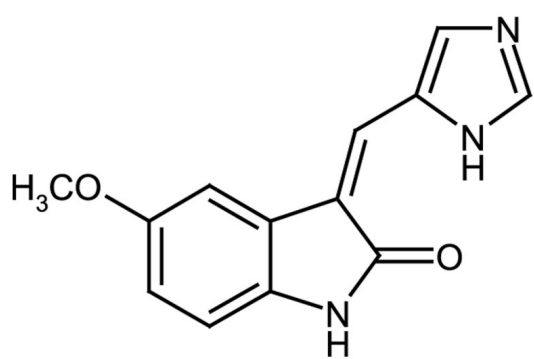
Ordering Information

[Order Online »](#)

ALX-270-400-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-[1-(3H-Imidazol-4-yl)-meth-(Z)-ylidene]-5-methoxy-1,3-dihydro-indol-2-one
Appearance	Yellow to orange solid.
CAS	377090-84-1
Couple Target	CDK, Cyclin
Couple Type	Inhibitor
Formula	$C_{13}H_{11}N_3O_2$
MW	241.3
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (10mg/ml).



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)