

RO-3306

CDK/cyclin inhibitor

Inhibitor of CDK1/cyclin B1 and CDK1/cyclin A. Reversibly arrests human cells at the G₂/M border of the cell cycle. Allows for effective cell synchronisation in early mitosis. Downregulates expression of the antiapoptotic proteins Bcl-2 and survivin and blocks p53-mediated induction of p21 and MDM2.

Citations: 31

[View Online »](#)

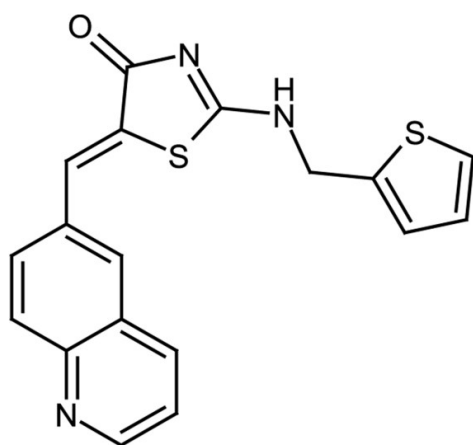
Ordering Information

[Order Online »](#)

ALX-270-463-M001	1mg
ALX-270-463-M005	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 6 months after receipt when stored at -20°C in the dark.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(5Z)-2-((Thiophen-2-yl)methylamino)-5-((quinolin-6-yl)methylene)thiazol-4(5H)-one
Appearance	Brown solid.
CAS	872573-93-8
Couple Target	CDK, Cyclin
Couple Type	Inhibitor
Formula	$C_{18}H_{13}N_3OS_2$
Identity	Determined by ¹ H-NMR and MS.
MW	351.4
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (50mM).

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-
be@enzolifesciences.com](mailto:info-be@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)