

CDK1/2 Inhibitor III

Cell permeable triazolo-diamine compound that displays anti-proliferative properties in various human cancer cells (IC_{50} =20nM, 35nM and 92nM in HCT-116, HeLa, and A375 cells). Acts as a highly potent, ATP-competitive inhibitor of CDK1/cyclin B and CDK2/cyclin A (IC_{50} =600pM and 500pM) with selectivity over VEGF-R2 and GSK-3 β (IC_{50} =32nM and 140nM).

Citations: 4

[View Online »](#)

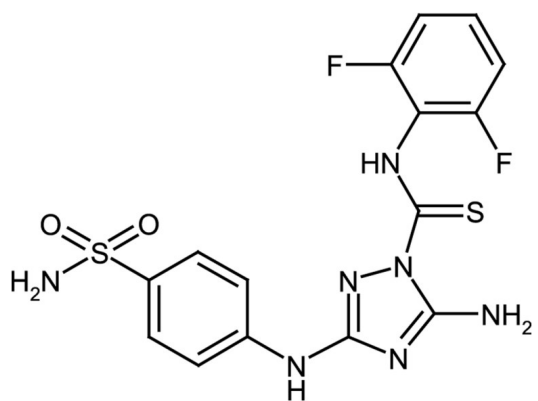
Ordering Information

[Order Online »](#)

ALX-270-442-M001	1mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	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.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	5-Amino-3-((4-(aminosulfonyl)phenyl)amino)-N-(2,6-difluorophenyl)-1H-1,2,4-triazole-1-carbothioamide
Appearance	White to off-white solid.
CAS	443798-55-8
Couple Target	CDK, GSK, VEGFR
Couple Type	Inhibitor
Formula	$C_{15}H_{13}F_2N_7O_2S_2$
Identity	Identity determined by NMR.
MW	425.4
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO (10mg/ml), tetrahydrofuran or acetone.



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)