

# CDK1/2 Inhibitor III

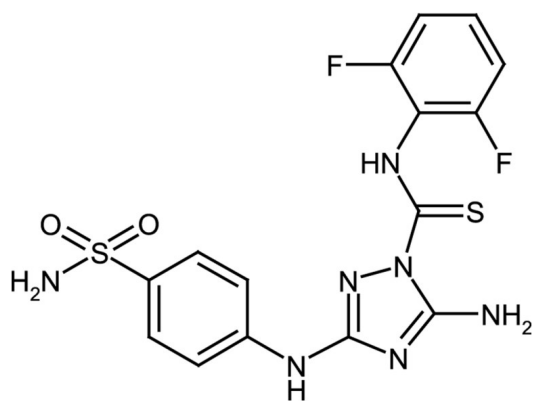
Cell permeable triazolo-diamine compound that displays anti-proliferative properties in various human cancer cells ( $IC_{50}$ =20nM, 35nM and 92nM in HCT-116, HeLa, and A375 cells). Acts as a highly potent, ATP-competitive inhibitor of CDK1/cyclin B and CDK2/cyclin A ( $IC_{50}$ =600pM and 500pM) with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

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

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

**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)