

# NU6102

## CDK/cyclin inhibitor

1,000-fold more potent inhibitor of CDK1/cyclin B ( $IC_{50}=9.5nM$ ) and CDK2/cyclin A3 ( $IC_{50}=5.4nM$ ) than NU2058 (Prod. No. ALX-270-394). Selective for CDK1 and CDK2 compared to CDK4/D1 ( $IC_{50}=1.6\mu M$ ), DYRK1A ( $IC_{50}=0.9\mu M$ ), PDK1 ( $IC_{50}=0.8\mu M$ ) and ROCKII ( $IC_{50}=0.6\mu M$ ).

Citations: 4

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-270-419-M001	1mg
------------------	-----

## 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	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Appearance	White to off-white solid.
CAS	444722-95-6
Couple Target	CDK, Cyclin
Couple Type	Inhibitor
Formula	$C_{18}H_{22}N_6O_3S$
MW	402.5
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO.

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)