

# Chk2 inhibitor

Cell-permeable, ATP-competitive, potent inhibitor of Chk2.

Displays anti-proliferative and anti-inflammatory properties. Chk2 inhibitor has also been reported to suppress the growth of leukemic T cells.

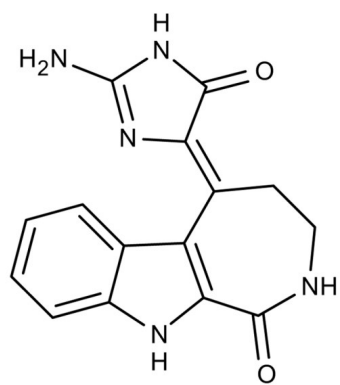
## Ordering Information

[Order Online »](#)

ENZ-CHM115-0001	1mg
ENZ-CHM115-0500	500µg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	5-(2-Amino-5-oxo-1,5-dihydroimidazol-4-ylidene)-3,4,5,10-2H-azepino[3,4-b]indol-1-one, Chk1 inhibitor II, Chk2 inhibitor I
Appearance	Yellow powder solid.
CAS	724708-21-8
Couple Target	Chk
Couple Type	Inhibitor
Formula	$C_{15}H_{13}N_5O_2$
MW	295.3
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO.



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)