

# KU55933

Potent and selective ATP-competitive inhibitor of ATM

KU55933 is a potent and selective ATP-competitive inhibitor of ATM, with an  $IC_{50}$  of 13 nmol/l and a  $K_i$  of 2.2 nmol/l. KU55933 has potential for use as a new radio- and chemo-sensitizing agent for the treatment of cancer.

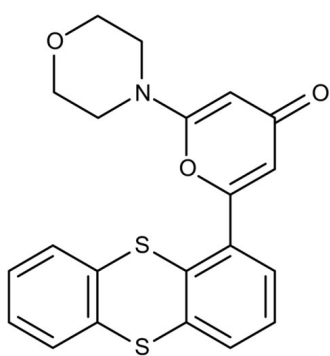
## Ordering Information

[Order Online »](#)

ENZ-CHM150-0010	10mg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	ATM kinase inhibitor, 2-Morpholino-6-(thianthren-1-yl)-4H-pyran-4-one
Appearance	Beige powder.
CAS	587871-26-9
Couple Target	ATM
Couple Type	Inhibitor
Formula	$C_{21}H_{17}NO_3S_2$
Identity	Determined by EM-MS, NMR
MW	395.5
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (50 mg/ml) or ethanol (

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-  
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)