

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.

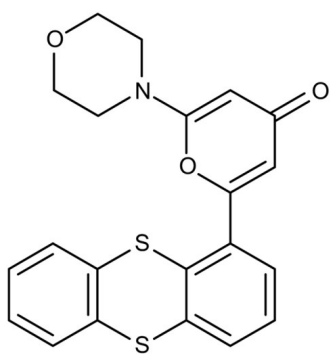
Ordering Information

[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 (



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com