

AZD7648

DNA-PK inhibitor/CRISPR enhancer

AZD7648 is a very potent and selective DNA-PK inhibitor. It is a sensitizer of radiation- and doxorubicin-induced DNA damage and sustained tumor regression in xenograft and patient-derived xenograft models. AZD7648 significantly enhances precise CRISPR/Cas9 gene editing alone or in combined treatment with a DNA polymerase theta inhibitor (2iHDR cocktail).

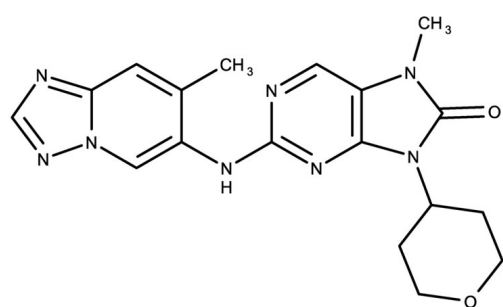
Ordering Information

[Order Online »](#)

ENZ-CHM449-0025	25mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 7-Methyl-2-[(7-methyl-[1,2,4]triazolo[1,5-a]pyridine-6-yl)amino]-9-(oxan-4-yl)purin-8-one

Appearance White solid.

CAS 2230820-11-6

Couple Type Inhibitor

Formula $C_{18}H_{20}N_8O_2$

Identity Determined by NMR.

MW 380.41

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (5 mg/ml with warming).



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)