

XL413 . HCl

Cdc7 inhibitor/CRISPR enhancer

XL413 is a potent ATP-competitive inhibitor of the important DNA replication initiation kinase Cdc7 (DDK). It increases the efficiency of homology directed DNA repair in CRISPR-Cas9 gene editing. XL413 acted synergistically with other chemotherapy agents in various cancer models.

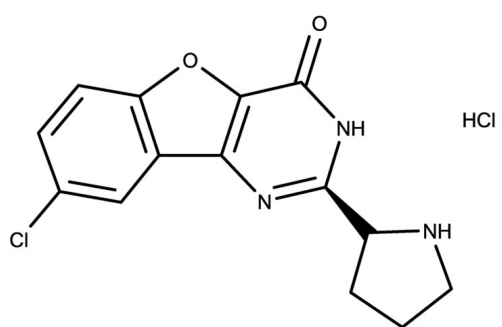
Ordering Information

[Order Online »](#)

ENZ-CHM453-0025	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 8-Chloro-2-[(2S)-pyrrolidin-2-yl]-3H-[1]benzofuro[3,2-d]pyrimidin-4-one hydrochloride

Appearance White solid.

CAS 2062200-97-7

Couple Type Inhibitor

Formula $C_{14}H_{12}ClN_3O_2 \cdot HCl$

Identity Determined by NMR.

MW 326.18

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (2 mg/ml with warming) or water (7 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)