

Asunaprevir

HCV NS3 protease inhibitor/CRISPR
enhancer

Asunaprevir is a potent hepatitis C virus NS3 protease inhibitor. Asunaprevir was used as part of a repressible Cas9 system (small molecule-assisted shut off - SMASh) capable of degrading newly synthesized Cas9 protein rapidly to precisely control gene editing. The Asunaprevir/SMASh system has also been using to target endogenous PD-1 protein for proteasomal degradation.

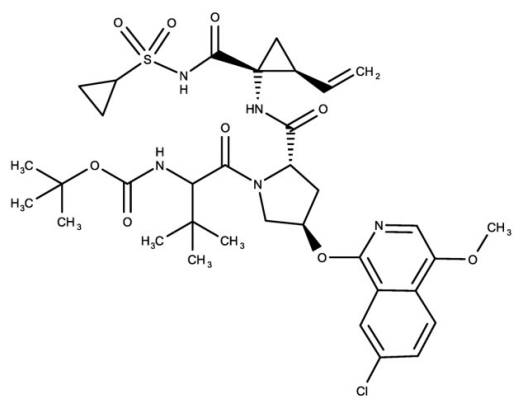
Ordering Information

[Order Online »](#)

ENZ-CHM460-0025	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 tert-Butyl N-[(2S)-1-[(2S,4R)-4-(7-chloro-4-methoxyisoquinolin-1-yl)oxy-2-[[[(1R,2S)-1-(cyclopropylsulfonylcarbamoyl)-2-ethenylcyclopropyl]carbamoyl]pyrrolidine-1-yl]-3,3-dimethyl-1-oxobutan-2-yl]carbamate

Appearance White solid.

CAS 630420-16-5

Couple Type Inhibitor

Formula $C_{35}H_{46}ClN_5O_9S$

Identity Determined by NMR.

MW 748.29

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (>50 mg/ml).

Last modified: January 24, 2025



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)