

GRL-0667

Noncovalent inhibitor of SARS-CoV PLpro

Potent noncovalent inhibitor of severe acute respiratory syndrome-coronavirus papain-like protease (SARS-CoV PLpro). GRL-0667 inhibited SARS-CoV viral replication in Vero E6 cells with no cytotoxicity.

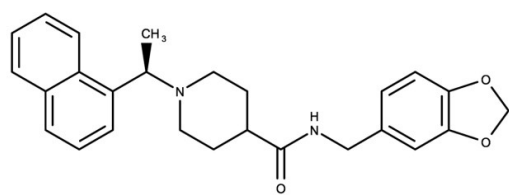
Ordering Information

[Order Online »](#)

ENZ-CHM478-0025	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 1-[(R)-1-(1-Naphthyl)ethyl]-4-[3,4-(methylenedioxy)benzylamino]carbonylpiperidine

Appearance Off-white solid.

CAS 1234195-91-5

Couple Target Protease

Couple Type Inhibitor

Formula $C_{26}H_{28}N_2O_3$

Identity Determined by NMR.

MW 416.52

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (>25 mg/ml) or ethanol (15 mg/ml).

Last modified: February 27, 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-
be@enzolifesciences.com](mailto:info-be@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)