

SARS-CoV-2 PLpro Protease Assay Service

SARS-CoV-2 PLpro [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

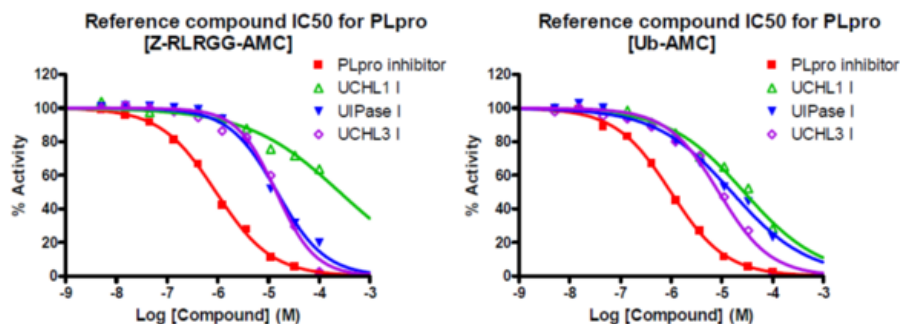
Target Overview

CAT#	PLpro
Enzyme	SARS-CoV-2 PLpro
Synonym	SARS-CoV-2 Papain-like Protease, NS3
Substrates	Fluorogenic substrate [Z-RLRGG-AMC], 10 μ M or [Ubiquitin-AMC], 100 nM

Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	<p>[Z-RLRGG-AMC] Compounds IC50 (μM)</p> <p>PLpro Inhibitor 0.92</p> <p>UCHL1 I 253</p> <p>UIPase I 14.8</p> <p>UCHL3 I 14.2</p> <p>[Ubiquitin-AMC] Compounds IC50 (μM)</p> <p>PLpro Inhibitor 0.98</p> <p>UCHL1 I 27.5</p> <p>UIPase I 15.6</p> <p>UCHL3 I 8.7</p>
Scientific Information	Accession #YP_009725299.1
Screening Location	Malvern, USA
Further information	More information can be found on our website Protease Assay Services .

Reference compound IC50 for SARS-CoV-2 PLpro



Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/sars-cov-2_plpro_protea_malvern

