

SARS-CoV-2 Mpro Protease Assay Service

SARS-CoV-2 Mpro [Protease Assay](#) (3CL Protease Assay) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

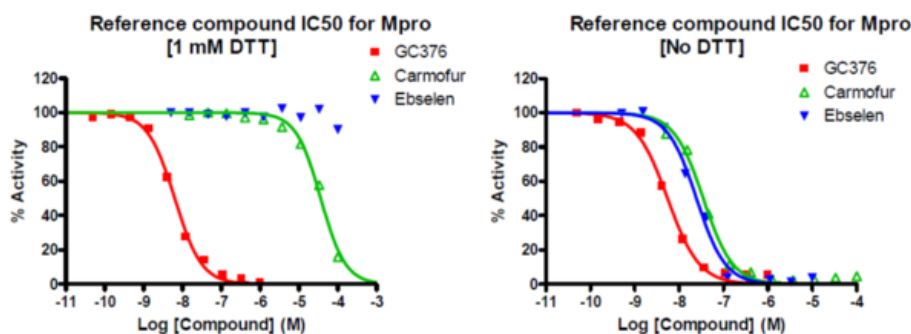
Target Overview

CAT#	Mpro
Enzyme	SARS-CoV-2 Mpro
Synonym	3CL Protease, SARS-CoV-2 Main Protease, SARS-CoV-2 3C-like Proteinase, 3CLpro
Substrates	Fluorogenic substrate [NH ₂ -C(EDANS)VNSTQSLRK(DABCYL)M-COOH], 5 μM

Assay Properties

Readout	Enzymatic activity												
Reference Compound IC50s	<p>Presence of 1 mM DTT Compounds IC50 (nM)</p> <table border="1"> <tr> <td>GC376</td> <td>6.5</td> </tr> <tr> <td>Carmofur</td> <td>36,000</td> </tr> <tr> <td>Ebselen</td> <td>>100,000</td> </tr> </table> <p>Absence of DTT Compounds IC50 (nM)</p> <table border="1"> <tr> <td>GC376</td> <td>5.5</td> </tr> <tr> <td>Carmofur</td> <td>35</td> </tr> <tr> <td>Ebselen</td> <td>25</td> </tr> </table>	GC376	6.5	Carmofur	36,000	Ebselen	>100,000	GC376	5.5	Carmofur	35	Ebselen	25
GC376	6.5												
Carmofur	36,000												
Ebselen	>100,000												
GC376	5.5												
Carmofur	35												
Ebselen	25												
Scientific Information	Accession #QHD43415.1												
Screening Location	Malvern, USA												
Further information	More information can be found on our website Protease Assay Services .												

Reference compound IC50 for SARS-CoV-2 Mpro



Let's Discover Together.

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

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

