

# ACE2 Protease Assay Service

ACE2 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

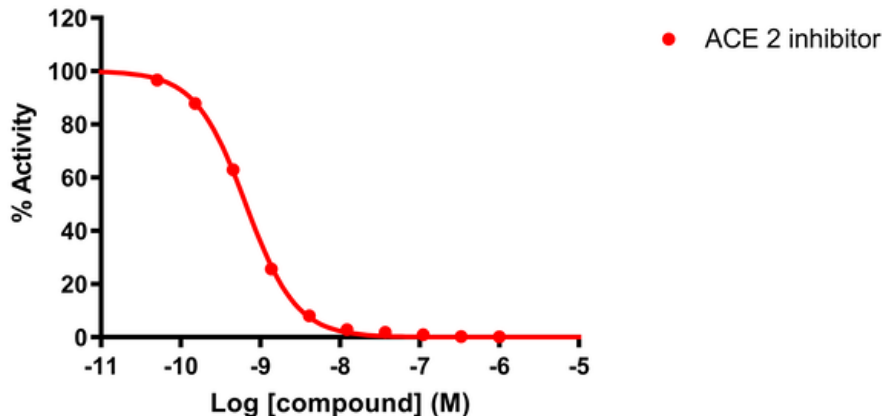
## Target Overview

CAT#	ACE2
Enzyme	Human ACE2
Synonym	Angiotensin-converting enzyme 2, ACE-related carboxypeptidase, Angiotensin-converting enzyme homolog, Metalloprotease MPROT15
Substrates	Fluorogenic peptide [MCA-YVADAPK(Dnp)], 10 $\mu$ M

## Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	<b>Compounds</b> IC50 (nM) ACE2 Inhibitor 0.5
Scientific Information	<a href="https://www.uniprot.org/uniprot/Q9BYF1">Q9BYF1</a> ( <a href="https://www.uniprot.org/uniprot/Q9BYF1">https://www.uniprot.org/uniprot/Q9BYF1</a> )
Screening Location	Malvern, USA
Further information	More information can be found on our website <a href="#">Protease Assay Services</a> .

## Reference Compound IC50 for ACE2



Let's Discover Together.

© Jun 2023 **Reaction Biology**. All rights reserved.

[https://www.reactionbiology.com/datasheet/ace2\\_protea\\_malvern](https://www.reactionbiology.com/datasheet/ace2_protea_malvern)

