

# TMPRSS2 Protease Assay Service

TMPRSS2 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

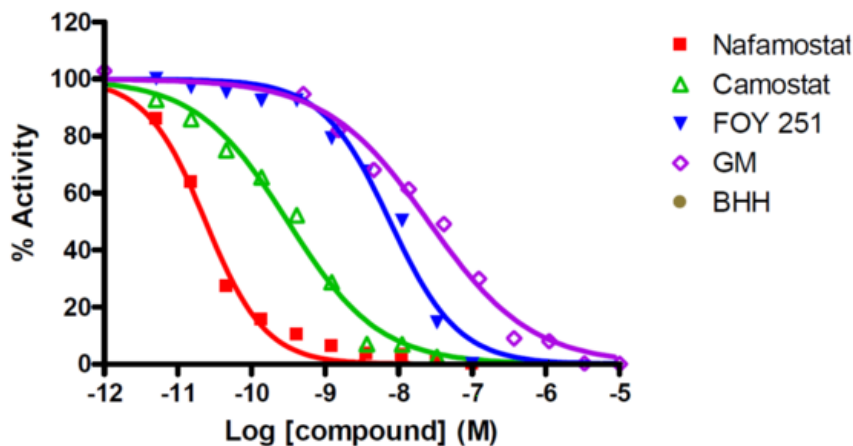
## Target Overview

<b>CAT#</b>	TMPRSS2
<b>Enzyme</b>	Human TMPRSS2
<b>Synonym</b>	Transmembrane protease serine 2, Serine protease 10, PRSS10, TMPS2
<b>Substrates</b>	Fluorogenic peptide [Boc-Gln-Ala-Arg-AMC], 25 $\mu$ M

## Assay Properties

<b>Readout</b>	Enzymatic activity	
<b>Reference Compound IC50s</b>	<b>Compounds</b>	<b>IC50 (nM)</b>
	Nafamostat Mesylate	0.024
	Camostat Mesylate	0.31
	FOY 251	7.6
	Gabexate Mesylate (GM)	25
	Bromhexine hydrochloride (BHH)	>100,000
<b>Scientific Information</b>	O15393 ( <a href="http://www.uniprot.org/uniprot/O15393">http://www.uniprot.org/uniprot/O15393</a> )	
<b>Screening Location</b>	Malvern, PA, USA	
<b>Further information</b>	More information can be found on our website <a href="#">Protease Assay Services</a> .	

## Reference compound IC50 for TMPRSS2



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/tmprrs2\\_protea\\_malvern](https://www.reactionbiology.com/datasheet/tmprrs2_protea_malvern)

