

Matriptase 2 Protease Assay Service

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

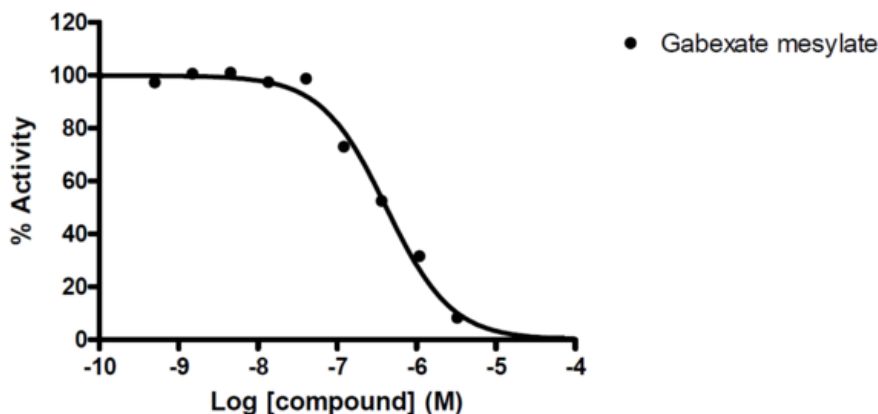
Target Overview

CAT#	Matriptase 2
Enzyme	Human Matriptase 2
Synonym	Transmembrane protease serine 6, TMPRSS6
Substrates	Fluorogenic peptide [Boc-QAR-AMC], 10 μ M

Assay Properties

Readout	Enzymatic activity				
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (μM)</th> </tr> </thead> <tbody> <tr> <td>Gabexate mesylate</td> <td>0.42</td> </tr> </tbody> </table>	Compounds	IC50 (μ M)	Gabexate mesylate	0.42
Compounds	IC50 (μ M)				
Gabexate mesylate	0.42				
Scientific Information	Q8IU80 (http://www.uniprot.org/uniprot/Q8IU80)				
Screening Location	Malvern, USA				

Reference compound IC50 for Matriptase 2



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/matriptase_2_protea_malvern

