

# MMP 7 Protease Assay Service

MMP 7 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

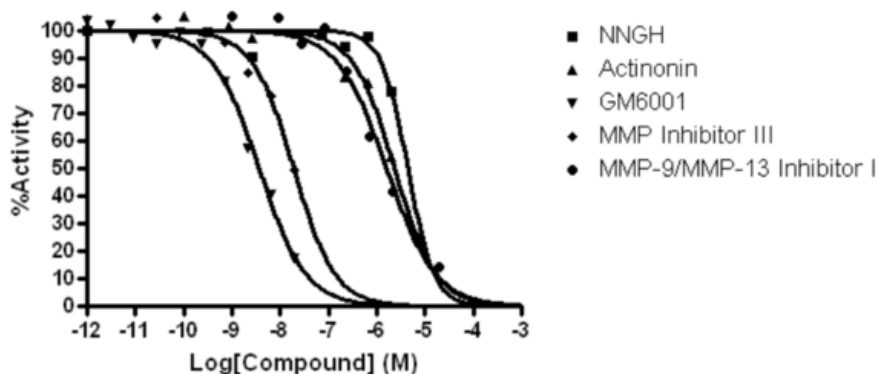
## Target Overview

<b>CAT#</b>	MMP7
<b>Enzyme</b>	MMP-7, catalytic domain from human recombinant (residues 78-250) expressed in E. coli with C-terminal tag
<b>Synonym</b>	Pump-1 protease, Uterine metalloproteinase, Matrix metalloproteinase-7, MMP-7, Matrin
<b>Substrates</b>	(5-FAM/QXL™) FRET peptide

## Assay Properties

<b>Readout</b>	Enzymatic activity	
<b>Reference Compound IC50s</b>	<b>Compounds</b>	<b>IC50 (M)</b>
	NNGH	4.40E-06
	Actinonin	2.33E-06
	GM6001	3.66E-09
	MMP Inhibitor III	1.90E-08
	MMP-9/MMP-13 Inhibitor I	1.61E-06
<b>Scientific Information</b>	<a href="http://www.uniprot.org/uniprot/P09237">P09237</a> ( <a href="http://www.uniprot.org/uniprot/P09237">http://www.uniprot.org/uniprot/P09237</a> )	
<b>Screening Location</b>	Malvern, USA	

## Reference compound IC50 for MMP-7



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/mmp\\_7\\_protea\\_malvern](https://www.reactionbiology.com/datasheet/mmp_7_protea_malvern)

