

# MMP 10 Protease Assay Service

MMP 10 (Catalytic domain) [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

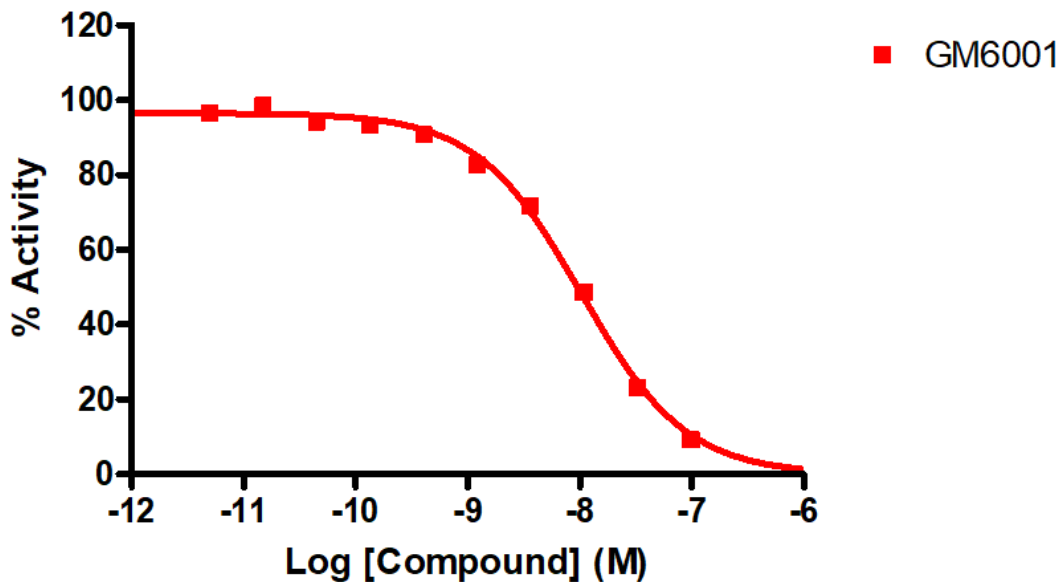
## Target Overview

CAT#	MMP-10
Enzyme	Human MMP-10
Synonym	Matrix metalloproteinase-10, Transin-2
Substrates	5 $\mu$ M Fluorogenic peptide 5-FAM/QXLTM

## Assay Properties

Readout	Enzymatic activity
Reference Compound IC50s	<b>Compounds IC50 (nM)</b> GM6001 13.4
Scientific Information	<a href="https://www.uniprot.org/uniprotkb/P09238/entry">P09238</a> ( <a href="https://www.uniprot.org/uniprotkb/P09238/entry">https://www.uniprot.org/uniprotkb/P09238/entry</a> )
Screening Location	Malvern, PA, USA
Further Information	More information can be found on our website <a href="#">Protease Assay Services</a>

## Reference compound IC50 for MMP-10 (Catalytic domain)



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/mmp\\_10\\_protea\\_malvern](https://www.reactionbiology.com/datasheet/mmp_10_protea_malvern)

