

# MMP 1 Protease Assay Service

MMP 1 [Protease Assay](#) for compound screening and profiling via fluorescence-based quantification of enzyme activity.

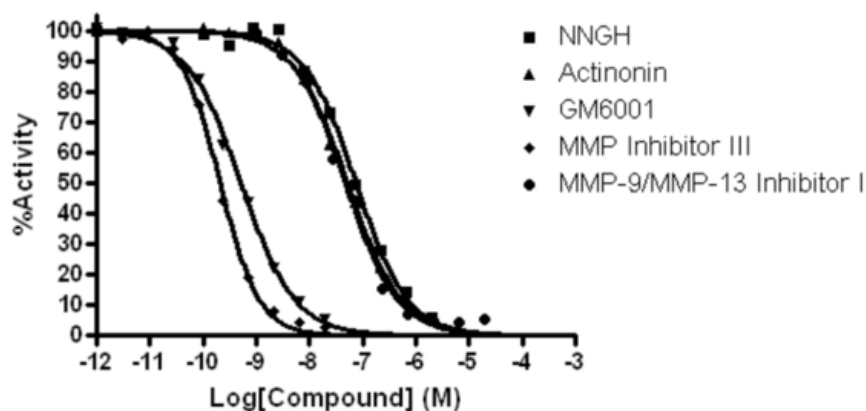
## Target Overview

|                   |   |
|-------------------|---|
| <b>CAT#</b>       | MMP1  |
| <b>Enzyme</b>     | MMP-1, catalytic domain from human recombinant (residues 81-249) expressed in E. coli with C-terminal tag |
| <b>Synonym</b>    | Matrix metalloproteinase-1, Fibroblast Collagenase  |
| <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   | 7.51E-08        |
|                                 | Actinonin  | 5.19E-08        |
|                                 | GM6001   | 5.19E-10        |
|                                 | MMP Inhibitor III  | 2.10E-10        |
|                                 | MMP-9/MMP-13 Inhibitor I   | 4.75E-08        |
| <b>Scientific Information</b>   | <a href="http://www.uniprot.org/uniprot/P03956">P03956</a> ( <a href="http://www.uniprot.org/uniprot/P03956">http://www.uniprot.org/uniprot/P03956</a> ) |                 |
| <b>Screening Location</b>       | Malvern, USA   |                 |

## Reference compound IC50 for MMP-1



Let's Discover Together.

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

[https://www.reactionbiology.com/datasheet/mmp\\_1\\_protea\\_malvern](https://www.reactionbiology.com/datasheet/mmp_1_protea_malvern)

