

MMP 2 Protease Assay Service

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

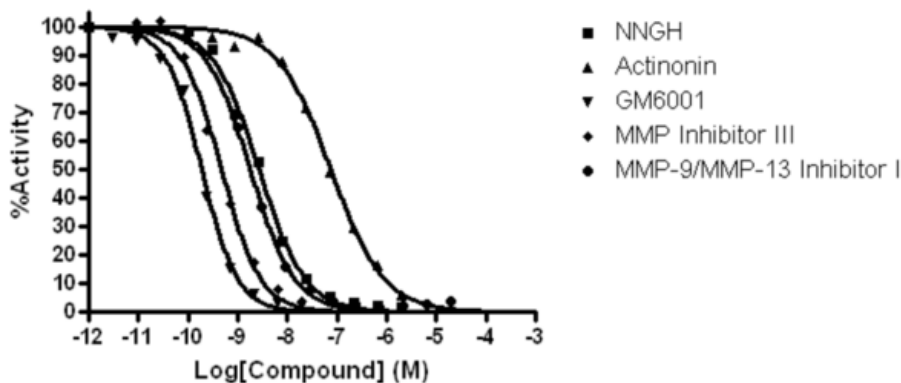
Target Overview

CAT#	MMP2
Enzyme	MMP-2, catalytic domain from human recombinant (residues 81-423) in E. coli with C-terminal tag
Synonym	Matrix metalloproteinase-2, MMP-2, Gelatinase A, TBE-1
Substrates	(5-FAM/QXL™) FRET peptide

Assay Properties

Readout	Enzymatic activity	
Reference Compound IC50s	Compounds	IC50 (M)
	NNGH	2.71E09
	Actinonin	7.50E-08
	GM6001	1.88E-10
	MMP Inhibitor III	4.70E-10
	MMP-9/MMP-13 Inhibitor I	1.79E-09
Scientific Information	P08253 (http://www.uniprot.org/uniprot/P08253)	
Screening Location	Malvern, USA	

Reference compound IC50 for MMP-2



Let's Discover Together.

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

https://www.reactionbiology.com/datasheet/mmp_2_protea_malvern

