

MMP-1

(Matrix metalloproteinase-1, Fibroblast Collagenase)

- **CAT #:** MMP1
- **Enzyme:** MMP-1, catalytic domain from human recombinant (residues 81-249) expressed in *E. coli* with C-terminal tag
- **Substrate:** (5-FAM/QXL™) FRET peptide
- **References Compound IC50s:**

<u>Compounds</u>	<u>NNGH</u>	<u>Actinonin</u>	<u>GM6001</u>	<u>MMP Inhibitor III</u>	<u>MMP-9/MMP-13 Inhibitor I</u>
IC50(M)	7.51E-08	5.19E-08	5.19E-10	2.10E-10	4.75E-08

- **Scientific Information:** [P03956](#)

* FAM = Carboxyfluorescein

Control Inhibitor IC50s for MMP-1

