

MMP-2

(Matrix metalloproteinase-2, MMP-2, Gelatinase A, TBE-1)

- **CAT #:** MMP2
- **Enzyme:** MMP-2, catalytic domain from human recombinant (residues 81-423) 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)	2.71E-09	7.50E-08	1.88E-10	4.70E-10	1.79E-09

- **Scientific Information:** [P08253](#)
- FAM = Carboxyfluorescein

Control Inhibitor IC50s for MMP-2

