

MMP-13

(Matrix metalloproteinase-13, MMP-13)

- **CAT #:** MMP13
- **Enzyme:** MMP-13, catalytic domain from human recombinant (residues 85-255) 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)	2.57E-09	2.16E-08	1.13E-10	4.40E-10	7.22E-10

- **Scientific Information:** [P45452](#)

* FAM = Carboxyfluorescein

Control Inhibitor IC50s for MMP-13

