

MMP-9

(MMP-9, 92 kDa type IV collagenase, 92 kDa gelatinase, Gelatinase B, GELB)

- **CAT #:** MMP9
- **Enzyme:** MMP-9, catalytic domain from human recombinant (residues 88-438) 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)	1.43E-09	6.35E-08	1.02E-10	3.63E-10	6.79E-10

- **Scientific Information:** [P14780](#)

* FAM = Carboxyfluorescein

Control Inhibitor IC50s for MMP-9

