

MMP-3

(Matrix metalloproteinase-3, MMP-3, Transin-1, SL-1)

- **CAT #:** MMP3
- **Enzyme:** MMP-3, catalytic domain from human recombinant (residues 83-255) 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-9/MMP-13 Inhibitor I</u>	<u>MMP-2/MMP-3 Inhibitor I</u>
IC50(M)	1.65E-08	7.23E-07	1.80E-8	1.41E-08	5.16E-06

- **Scientific Information:** [P08254](#)

* FAM = Carboxyfluorescein

Control Inhibitor IC50s for MMP-3

