

MMP-14

(MMP-14, Membrane-type matrix metalloproteinase 1, MT-MMP 1, MTMMP1, Membrane-type-1 matrix metalloproteinase, MT1-MMP, MT1MMP, MMP-X1)

- **CAT #:** MMP14
- **Enzyme:** MMP-14, catalytic domain from human recombinant (residues Tyr92-Arg278) 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)	2.53E-09	7.93E-07	4.71E-10	3.70E-09	4.19E-09

- **Scientific Information:** [P50281](#)
- FAM = Carboxyfluorescein

Control Inhibitor IC50s for MMP-14

