

MMP-12

(HME, Matrix metalloproteinase-12, MMP-12, Macrophage elastase, ME)

- **CAT #:** MMP12
- **Enzyme:** MMP-12, catalytic domain from human recombinant (residues 84-255) expressed in *E. coli* with C-terminal tag
- **Substrate:** (5-FAM/QXL™) FRET peptide
- **References Compound IC50s:**

Compounds	NNGH	Actinonin	GM6001	MMP Inhibitor III	MMP-9/MMP-13 Inhibitor I
IC50(M)	2.04E-09	1.36E-07	3.07E-10	2.35E-09	1.41E-09

- **Scientific Information:** [P39900](#)

* FAM = Carboxyfluorescein

Control Inhibitor IC50s for MMP-12

