

# MMP-7

(Pump-1 protease, Uterine metalloproteinase, Matrix metalloproteinase-7, MMP-7, Matrin)

- **CAT #:** MMP7
- **Enzyme:** MMP-7, catalytic domain from human recombinant (residues 78-250) expressed in *E. coli* with C-terminal tag
- **Substrate:** (5-FAM/QXL™) FRET peptide
- **References Compound IC50s:**

Compounds	NINGH	Actinonin	GM6001	MMP Inhibitor III	MMP-9/MMP-13 Inhibitor I
IC50(M)	4.40E-06	2.33E-06	3.66E-09	1.90E-08	1.61E-06

- **Scientific Information:** [P09237](#)

\* FAM = Carboxyfluorescein

**Control Inhibitor IC50s for MMP-7**

