



Product Information Sheet

Product Name:	Human MMP-8, Recombinant
Catalog Number:	72008
Amount:	10 µg/ml, 100µl
Molecular weight:	75 kD
Activity (Unit/µg):	>250 Unit/µg.
Unit definition:	One unit of protease hydrolyzes 1 picomole of Mca-Pro-Leu-Gly-Leu-Dap(Dnp)-Ala-Arg-NH ₂ (AnaSpec Cat#27076) per minute at pH 7.5 at 25° C.
Storage:	Store at -80°C. Avoid multiple thaw-freeze cycles.

Instruction:

Matrix metalloproteinases (MMPs) belong to a family of secreted or membrane-associated zinc endopeptidases capable of digesting extracellular matrix components. MMP-8 (neutrophil collagenase) is synthesized by neutrophils and stored in the specific granules until it is secreted.¹

Human MMP-8 was expressed as a full length pro-enzyme in *E. coli*. Proenzyme can be activated by incubating with 1 mM APMA at 37°C for 1 hr. and generate 64-kDa active form. Its activity can be measured in FRET-based enzymatic assays (AnaSpec Cat#71133, Cat#71154). 10-20 ng of enzyme is sufficient for FRET-based assay.

MMP-8 is stored in 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM CaCl₂, 1 mg/mL BSA.

References

1. Bainton, D. et al. *J. Exp. Med.* 134, 907 (1971).