

## **Product Information Sheet**

Product Name:	Human MMP-8, proenzyme (human neutrophils)
Catalog Number:	AS-72008
Size:	1 µg
Concentration:	10 µg/ml
Molecular weight:	85 kDa
Activity (Unit/µg):	Provided on the label
Unit definition:	One unit of protease hydrolyzes 1 picomole of Mca-Pro-Leu-Gly-Leu- Dap(Dnp)-Ala-Arg-NH <sub>2</sub> (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.<sup>1</sup>

Human MMP-8 is a full length pro-enzyme isolated from stimulated human neutrophil granulocytes. Proenzyme can be activated by incubating with 1 mM APMA at 37°C for 1 hr. 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 0.1M Tris, pH 7.5, 0.1M NaCl, 10 mM CaCl2, 0.05% Brij 35, 2mM sodium azide, 1 mg/mL BSA.

## References

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