

## **Product Information Sheet**

Product Name: MMP-3, Human, Catalytic domain

Catalog Number: AS-72006

Size:  $1 \mu g$ 

Concentration: 10 µg/mL

Activity (Unit/µg): Provided on the label

Unit definition: One unit of protease hydrolyzes 1 picomole of Mca-Arg-Pro-Lys-Pro-Val-Glu

Nva-Trp-Arg-Lys(Dnp)-NH2 (AnaSpec Cat#AS-27114) per minute at pH 7.5 at

25°C.

Storage: Store at -80°C. Avoid multiple freeze/thaw cycles.

## **Instruction:**

Matrix metalloproteinases (MMPs) belong to a family of secreted or membrane-associated zinc endopeptidases capable of digesting extracellular matrix components.  $^{1,2}$  The importance of MMPs in tumor development and invasion as well as other diseases is well known. MMP-3 (stromelysin-1, transin-1) has been shown to involved in tumor metastasis and rheumatoid arthritis  $^{4,5}$ . Therefore it is proposed as a therapeutic target for these diseases. The native pro-MMP-3 is Mr 59/57-kDa doublet, which can be autocatalyzed to an active form of 45-kDa, and is then processed partially to a second active form of 28-kDa.  $^{6}$ 

Recombinant human MMP-3 was expressed as catalytic domain in *E. coli*. The molecular mass is 19.5 kDa. Purity is > 95% by SDS-PAGE.

Recombinant human MMP-3 enzyme catalytic domain does not need APMA activation before enzyme assay. Its activity can be measured in FRET-based enzymatic assays (AnaSpec Cat#71130, Cat#71152). 10-20 ng of enzyme is sufficient for FRET-based assay. MMP-3 is stored in 50 mM HEPES, pH 7.5, 0.05% Brij 35, 10 mM CaCl<sub>2</sub>, 1 mg/mL BSA, 2mM sodium azide.

## References

- 1. Woessner, J. et al. J. Biol. Chem. 263, 16918 (1988).
- 2. Woessner, J. et al. FASEB. J. 5, 2145 (1991).
- 3. Matrisian, L. et al. *Proc. Natl. Acad. Sci. U.S.A* 83, 9413 (1986).
- 4. Chin, J. et al. J. Biol. Chem. 260, 12367 (1985).
- 5. Okada, Y. et al. J Biol. Chem. 261, 14245 (1986).
- 6. Okada, Y. et al. J. Biol. Chem. 261, 14245 (1986)