



Product Information Sheet

Revision#1.3	Last Updated: October 2024
Catalog #	AS-72007

Product Name:	MMP-7, Human, Catalytic domain
Catalog Number:	AS-72007
Size:	1 µg
Concentration:	10 µg/mL
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 ₂ 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.^{1,2} The importance of MMPs in tumor development and invasion as well as other diseases is well known. MMP-7,³ (matrilysin, PUMP-1) is proposed as a potential anti-cancer drug target.

Recombinant human MMP-7 was expressed in E.coli as catalytic domain. The molecular mass of the catalytic domain is 19-kDa. It's activity can be measured by FRET peptides (AnaSpec Cat#71153). 5-10 ng of enzyme is sufficient for FRET-based assay.

MMP-7 is stored in 0.1M Tris, pH 7.5, 0.1M NaCl, 10mM CaCl₂, 0.05% Brij 35, 2mM sodium azide, and 1mg/mL BSA.

References

1. Woessner, J. et al. *J. Biol. Chem.* 263, 16918 (1988).
2. Woessner, J. et al. *FASEB. J.* 5, 2145 (1991).