

MMP-7 (catalytic domain) (human), (recombinant)

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE181-0010	10µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Activity	Preincubation of MMP-7 catalytic domain at 19nM with the broad-spectrum inhibitor GM6001 (Prod. No. BML-EI300) at 100nM for 1 hour completely inhibits enzymatic activity.
Alternative Name	Matrix metalloproteinase 7, Matrilysin, Pump
Application Notes	Useful tool to study enzyme kinetics, cleave target substrates, and screen for inhibitors.
Formulation	Liquid. In 50 mM Tris-HCl pH 7.5, 5 mM CaCl ₂ , 300 mM NaCl, 20 µM ZnCl ₂ , 0.5% Brij-35, and 30% glycerol.
MW	20.4 kDa
Purity	≥95% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>E. coli</i> . Active Matrix Metalloproteinase-7 (MMP-7, matrilysin, pump) catalytic domain from human cDNA. The enzyme consists of the catalytic domain of human MMP-7 (Tyr ⁹⁵ -Lys ²⁶⁷ , NM_002423) with a C-terminal purification tag.
Specific Activity	≥200 pmol/min/µg at 37°C using the colorimetric thiopeptolide Ac-Pro-Leu-Gly-S-Leu-Leu-GlyOEt (100 µM; Prod. No. BML-P125) as substrate.
UniProt ID	P09237