## MMP-13 (catalytic domain) (human), (recombinant)

Citations: 8

View Online »

**Ordering Information** 

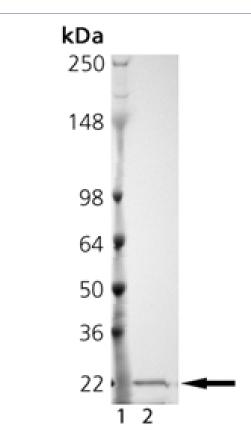
Order Online »

BML-SE246-0010

10µg

Manuals, SDS & CofA

**View Online »** 



SDS-PAGE Analysis: Lane1: MW Marker; Lane 2: 2µg of purified MMP-13 (catalytic domain) (human), (recombinant).

## **Handling & Storage**

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

## **Product Details**

**Activity** Preincubation of MMP-13 catalytic domain at 13nM with

the broad-spectrum inhibitor GM6001 (Prod. No. BML-El300) at 20nM for 1 hour completely inhibits enzymatic

activity.

Alternative Name Matrix metalloproteinase 13, Collagenase-3

**Application Notes**Useful tool to study enzyme kinetics, cleave target

substrates, and screen for inhibitors.

Formulation Liquid. In 50mM TRIS, 5mM CaCl<sub>2</sub>, 300mM NaCl, 20µM

ZnCl<sub>2</sub>, 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 *E. coli*. Active Matrix Metalloproteinase-13

(MMP-13, collagenase-3) catalytic domain from human cDNA. The enzyme consists of the catalytic domain of human MMP-13 (Tyr<sup>104</sup>-Asn<sup>274</sup>, NM\_002427) with a Cterminal purification tag. This represents a naturally-occurring active form of MMP-13 which lacks the Cterminal hemopexin domain. MMPs lacking this domain cannot cleave native collagens; however, activity toward other targets such as gelatin, casein, or peptide substrates

is unaffected.

Specific Activity ≥2000 pmol/min/µg at 37°C using the colorimetric

thiopeptolide Ac-Pro-Leu-Gly-S-Leu-Leu-Gly-OEt (100 μM;

Prod. No. BML-P125) as substrate.

UniProt ID P45452

Last modified: May 29, 2024



info-

eu@enzolifesciences.com