## MMP-13 (human) polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ALX-210-833-R100

100µl

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

**Handling** Avoid freeze/thaw cycles.

Long Term Storage -80°C

Shipping Dry Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name Collagenase-3, Matrix metalloproteinase 13

**Application** ELISA, WB

Crossreactivity Does not cross-react with human MMP-2, MMP-9, MMP-

14, MMP-15, MMP-16, MMP-17 or MMP-24.

**Formulation** Liquid. In PBS, pH 7.5, containing 0.01% thimerosal.

**Host** Rabbit

Immunogen Recombinant human MMP-13 (matrix metalloproteinase

13).

Purity Detail Protein A-affinity purified.

**Recommendation Dilutions/Conditions** ELISA (detection, 1:2,000-1:4,000, use Prod. No. ALX-

804-365 for capture)Western Blot (ECL-detection, 1:2,500 – 1:8,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be

determined individually for each application.

Species Reactivity Human

UniProt ID P45452

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee



European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 infofr@enzolifesciences.com Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com