

MMP-13 (human) polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

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](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-](mailto:info-usa@enzolifesciences.com)
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-](mailto:info-eu@enzolifesciences.com)
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-](mailto:info-be@enzolifesciences.com)
be@enzolifesciences.com

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
[info-](mailto:info-de@enzolifesciences.com)
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-](mailto:info-uk@enzolifesciences.com)
uk@enzolifesciences.com

Last modified: May 29, 2024