

MMP-2 (mouse fibroblasts), (purified)

Ordering Information

[Order Online »](#)

BML-SE505-0005	5µg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Matrix metalloproteinase 2, Gelatinase A, 72 kDa Type IV collagenase

Application Notes Useful for the study of enzyme regulation and kinetics, cleavage of substrates *in vitro* or in tissue culture, zymography or immunoblot positive control.

Formulation Liquid. In 50mM Tris-HCl, pH 7.0, containing 200mM NaCl, 5mM CaCl₂, 1µM ZnCl₂, 0.05% sodium azide and 0.05% Brij 35.

Purity Detail Free of TIMPs and other MMPs.

Source Isolated from mouse fibroblasts. Primarily latent form.

UniProt ID P08253



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

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

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

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

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

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