

Matriptase-2 (human), (recombinant)

Integral player in iron homeostasis, acts as a potent negative regulator of hepcidin.

Matriptase-2 (TMPRSS6) is a type II transmembrane serine protease mainly produced by the liver. It acts as a potent negative regulator of hepcidin, a hepatic peptide hormone and integral player in regulating iron homeostasis.

Matriptase-2 plays an important role in understanding iron disorders; including various anemias and iron overload conditions, affecting millions of people worldwide.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-752-1	125U
ALX-201-752-0500	500U

Manuals, SDS & CofA

[View Online »](#)

- Highly active serine protease providing superior results.
- Potent negative regulator of hepcidin.
- Eukaryotic cell-derived protein guarantees proper PTM.

Handling & Storage

Use/Stability	Stable for at least 6 months when stored at -20°C. Recommended Assay Buffer: 100mM TRIS, pH 9.0, containing 0.5mg/ml BSA. Do not store diluted solutions.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Activity	≥2.5 U/μl Please see Certificate of Analysis for detailed activity information.
Alternative Name	Transmembrane protease serine 6, TMPRSS6
Biological Activity	Trypsin-like serine protease.
Formulation	Liquid. In 100mM TRIS-HCl, pH 9.0, containing 150 mM sodium chloride and 20% glycerol.
MW	~84kDa; a-subunit: ~56kDa; b-subunit: ~28kDa
Purity Detail	Purified by multi-step chromatography.
Source	Produced in <i>Drosophila</i> Schneider 2 (S2) cells. Recombinant human matriptase-2 (aa 78-811) is fused at the C-terminus to a V5-His-tag. Produced by plasmid transfection expression.
Specific Activity	One U = 10 pmol/min, using Boc-Gln-Ala-Arg-AMC (50μM, Prod. No. BML-P237) as substrate at 37°C.

UniProt ID	Q8IU80
------------	--------



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024