

Matriptase (human), (recombinant) (His- tag)

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-246-U250	250U
------------------	------

Manuals, SDS & CofA

[View Online »](#)

kDa

75—

50—

37—

25—

20—



← Activated matriptase

Handling & Storage

Use/Stability	Stable for at least 6 months when stored at -20°C. 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	Trypsin-like serine protease.
Alternative Name	MT-SP1, Membrane-type serine protease 1, ST14, Suppression of tumorigenicity protein 14, TADG-15, Tumor-associated differentially expressed gene 15 protein, SNC19
Concentration	100nM
Formulation	Liquid. In 50mM TRIS, pH 8.0, containing 200mM sodium chloride, 1mM β -mercaptoethanol and 10% glycerol.
MW	~25kDa
Purity	≥98% (SDS-PAGE)
Source	Recombinant human matriptase (aa 596-855) fused at the N-terminus to a His-tag.
Specific Activity	~2.5U/μl. One unit is defined as the amount of enzyme that releases 10pmol of AMC per min. at 37°C in the presence of 50μM of Boc-Gln-Ala-Arg-AMC.
Technical Info / Product Notes	Recommended Assay Buffer is HEPES 50mM pH 7.4, 150 mM NaCl and 5mg/ml BSA
UniProt ID	Q9Y5Y6



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)