

DESC1 (human), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-201-751-1	125U
---------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Stable for at least 6 months when stored at -20°C. Recommended Assay Buffer: HEPES 50mM, pH 7.4, 150mM NaCl, 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

Alternative Name	Transmembrane protease serine 11E, TMPRSS11E
Biological Activity	Trypsin-like serine protease.
Concentration	100nM
Formulation	Liquid. In 100 mM Tris-HCl, pH 8.0, containing 150 mM sodium chloride and 20% glycerol.
MW	~45kDa; a-subunit: ~13kDa; b-subunit: ~32kDa
Source	Produced in <i>Drosophila</i> Schneider 2 (S2) cells. Recombinant human DESC1 (aa 44-423) is fused at the C-terminus to a V5-His-tag. Produced by plasmid transfection expression.
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 Boc-Gln-Ala-Arg-AMC.
UniProt ID	Q9UL52

Last modified: May 29, 2024



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