

# USP21 (human) polyclonal antibody

USP21 is a ubiquitin-specific protease that catalyzes the hydrolysis of ubH2A thereby acting to relieve ubH2A-specific repression. USP21 is also thought to be involved with hepatocyte regeneration. Furthermore, in addition to its deubiquitylating activity, USP21 is capable of removing NEDD8 from NEDD8 conjugates but has no effect on SUMO conjugates.

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-PW0585-0025	25µl
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

250kDa -

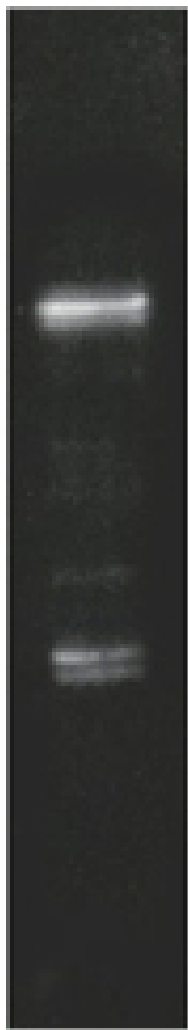
148kDa -

98kDa -

64kDa -

50kDa -

36kDa -



- USP21

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 1 year after receipt when stored at -20°C. Dilute with PBS, pH 7.2-7.4, containing 1% normal goat serum (if a goat anti-rabbit IgG linker antibody is to be used). Store diluted antibody at +4°C, do not freeze and use within 1 month.
<b>Handling</b>	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Ubiquitin thioesterase 21, Ubiquitin-specific-processing protease 21, Deubiquitinating enzyme 21, NEDD8-specific protease
<b>Application</b>	WB
<b>Formulation</b>	Liquid. In PBS containing 0.01% sodium azide.
<b>Host</b>	Rabbit
<b>Immunogen</b>	Synthetic peptide corresponding to aa 60-74 of human USP21.
<b>Purity Detail</b>	Purified by salt precipitation.
<b>Recommendation Dilutions/Conditions</b>	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Source</b>	Purified from rabbit serum.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	Q9UK80

Last modified: May 29, 2024



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)