

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 deubiquitinating 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](#).

Ordering Information

[Order Online »](#)

BML-PW0585-0025	25µl
BML-PW0585-0100	100µl

Manuals, SDS & CofA

[View Online »](#)

250kDa -

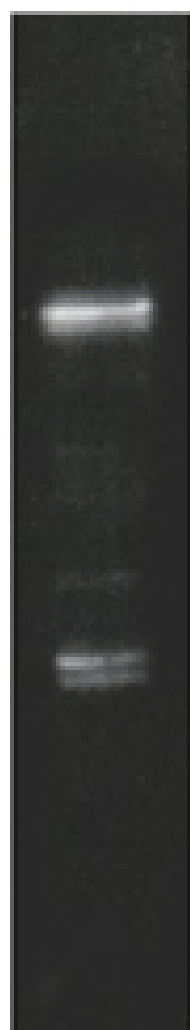
148kDa -

98kDa -

64kDa -

50kDa -

36kDa -



- USP21

Handling & Storage

Use/Stability	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.
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

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



Worry-Free Guarantee

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

This antibody is covered by our Worry-Free Guarantee

Belgium, The Netherlands
& Luxembourg

Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France

Phone: +33 1 72 440 1335
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