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 deubiquitinylating 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.

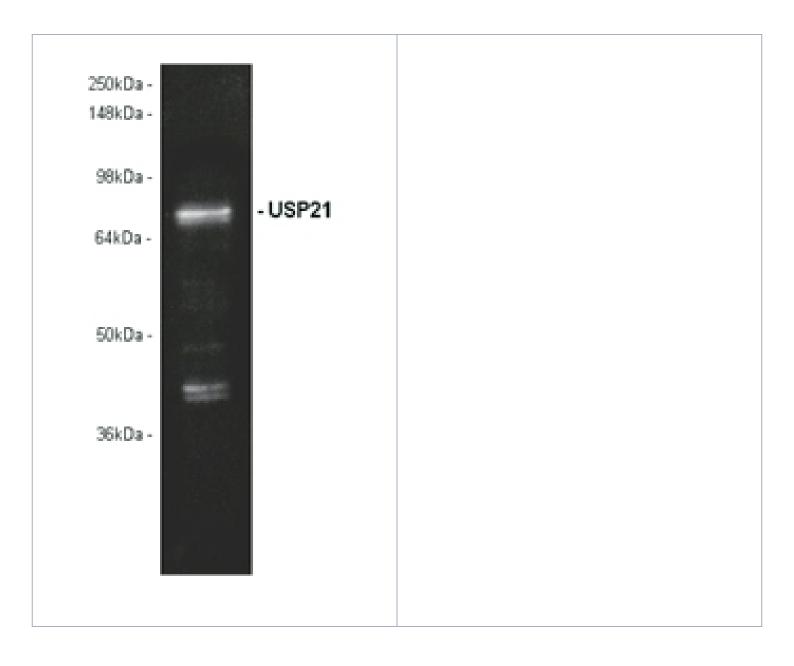
Ordering Information

Order Online »

BML-PW0585-0025	25µI
BML-PW0585-0100	100μΙ

Manuals, SDS & CofA

View Online »



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/ConditionsWestern 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com