

USP19 polyclonal antibody

The deubiquitylating enzyme USP19 is induced in skeletal muscle atrophy in response to numerous catabolic stimuli, including cancer, fasting, diabetes, and glucocorticoids. Loss of USP19 interferes with the growth of L6 myoblasts through its modulation of the transcription of major myofibrillar proteins.

USP19 regulates cell growth, partly through indirect control of p27Kip1 degradation in the G1 phase of the cell cycle. Since a number of cancers are associated with enhanced degradation of p27Kip1, the overexpression of USP19 may also be a pathogenetic mechanism of the disordered cell growth in cancer and a potential novel therapeutic target for its treatment.

In addition USP19 has been shown to be a tail anchored DUB, localised to the ER, suggesting a role in the turnover of ERAD substrates and to be targeted by the unfolded protein response.

USP19 is expressed in a wide range of tissues and is localised to the cytoplasm and endoplasmic reticulum.

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

Ordering Information

[Order Online »](#)

BML-PW1025-0025	25µl
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

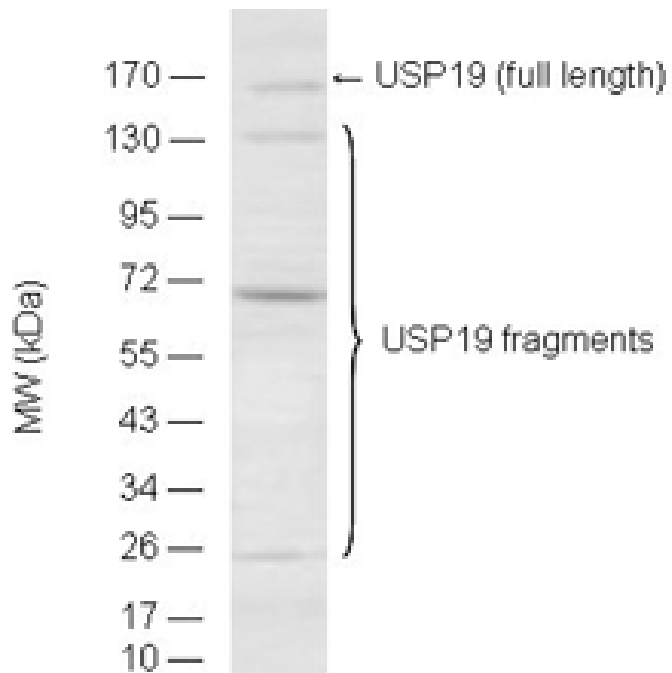


Figure: Detection of full length USP19 and its degradation fragments in mouse testis lysate (50 μg) by Western blot using USP19, pAb (Prod. No. BML-PW1025) (1:100).

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
Handling	Do not freeze.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	ZMYND9, Zinc finger MYND domain-containing protein 9
Application	WB
Formulation	Liquid. In 100mM TRIS, pH 7.0 containing 100mM glycine and 0.02% sodium azide.
Host	Rabbit
Immunogen	Recombinant fragment of human USP19 (aa 1-129).
Purity Detail	Immunoaffinity purified.
Recommendation Dilutions/Conditions	Western Blot (1:100)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, Mouse, Rat
UniProt ID	O94966

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
Last modified: **May 29, 2024**
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com
uk@enzolifesciences.com