USP19 polyclonal antibody

The deubiquitinylating 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	25µI
BML-PW1025-0100	100µl

Manuals, SDS & CofA

View Online »

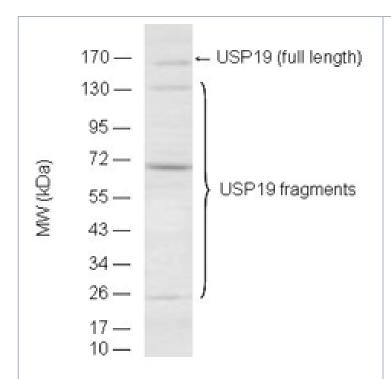


Figure: Detection of full length USP19 and its degradation fragments in mouse testis lysate (50μg) by Western blot using USP19, pAb (Prod. No. BML-PW1025) (1:100).

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/ConditionsWestern 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 094966

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

