

## USP18 (Ubiquitin specific protease 18)

Cat. # DB519

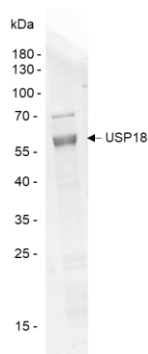
**Background** USP18 (Ubiquitin specific protease 18, ISG15-specific protease) is a member of the USP family of deubiquitinases and specifically removes the ubiquitin-like molecule ISG15 from conjugated proteins. It is involved in the regulation of the inflammatory response to interferon type 1.

**Alternate names** 43 kDa ISG15-specific protease, hUBP43, ISG15-specific-processing protease, ISG43, Ubl carboxyl-terminal hydrolase 18, Ubl thioesterase 18, UBP43

### Product Information

<b>Purity</b>	≥ 85% by SDS-PAGE
<b>Accession No.</b>	Q9UMW8
<b>Molecular Weight</b>	55 kDa
<b>Quantity</b>	25 µg
<b>Physical State</b>	Liquid, 50 mM Hepes, pH 7.5, 200 mM NaCl, 2 mM DTT, 0.5 mM EDTA, 10% glycerol
<b>Source</b>	Human Recombinant
<b>Tag</b>	His6-SUMO
<b>Activity</b>	This enzyme is active in the ISG15-CHOP assay.
<b>Storage</b>	-80° C. Avoid repeated freeze/thaw cycles

### Product QC



**SDS-Page Analysis of purified USP18.** Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.

### References

1. Arimoto, KI., et al., Nat Commun., 2023. 14(1):251.
2. Fernández, D., et al., Front Chem., 2020. 7:923.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2009 LifeSensors, Inc. All Rights Reserved