

## USP34core (Ubiquitin Specific Protease 34 catalytic domain)

Cat. # DB506

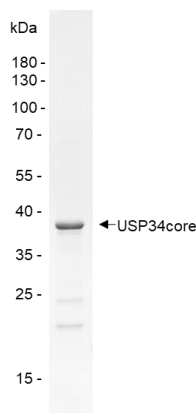
**Background** USP34 (Ubiquitin Specific Protease 34) is a member of the USP family of cysteine proteases. The deubiquitinating action of USP34 leads to the stabilization and nuclear accumulation of AXIN1 and AXIN2 for a positive regulation of beta-catenin (CTNBB1)-mediated transcription. USP34core consists of the catalytic residues 1892-2241.

**Alternate names** FLJ43910, KIAA0570, KIAA0729

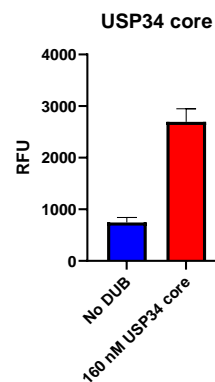
### Product Information

<b>Purity</b>	≥ 85% by SDS-PAGE
<b>Accession No</b>	Q70CQ2
<b>Molecular Weight</b>	42 kDa
<b>Quantity</b>	25 µg
<b>Physical State</b>	Liquid, 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol
<b>Source</b>	Human Recombinant
<b>Tag</b>	His6
<b>Activity</b>	This enzyme is active in the Ub-CHOP assay.
<b>Storage</b>	-80° C. Avoid repeated freeze/thaw cycles

### Product QC



**SDS-Page Analysis of purified USP34core.** Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



**Activity Assay of ATXN3-Like.** Ataxin3-like Protein was tested in a Ub-CHOP assay showing robust DUB activity.

### References

1. Gu, Z., et al., Biol Pharm Bull., 2019. 42(4):573-579.
2. Oh, E., et al., Cell Signal., 2017. 36:230-239.

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