

## USP47 (Ubiquitin specific protease 47)

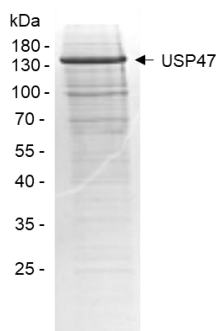
Cat. # DB510

|                        |                                                                                                                                                                             |
|------------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Background</b>      | USP47 (Ubiquitin specific protease 47) deubiquitinates monoubiquitinated DNA polymerase beta (POLB), stabilizing POLB thereby playing a role in base-excision repair (BER). |
| <b>Alternate names</b> | Deubiquitinating enzyme 47, Ubiquitin thioesterase 47, Ubiquitin carboxyl-terminal hydrolase 47                                                                             |

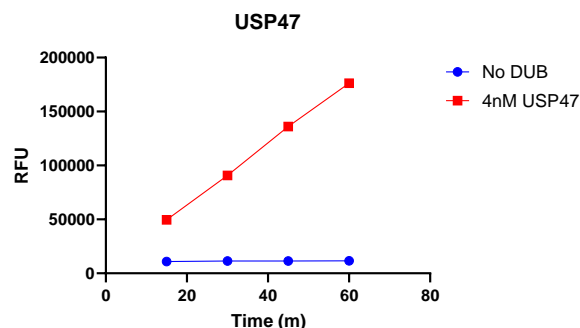
### Product Information

|                         |                                                                     |
|-------------------------|---------------------------------------------------------------------|
| <b>Purity</b>           | ≥ 90% by SDS-PAGE                                                   |
| <b>Accession No.</b>    | Q96K76                                                              |
| <b>Molecular Weight</b> | 157 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>         | Active in the Ub-CHOP assay                                         |
| <b>Storage</b>          | -80° C. Avoid repeated freeze/thaw cycles                           |

### Product QC



**SDS-Page Analysis of purified USP47.** Two µg of the enzyme was loaded on a 10-20% SDS-PAGE gel and stained with Coomassie brilliant blue.



**Activity Assay of USP47.** 4 nM USP47 was tested in a Ub-CHOP assay showing robust DUB activity.

### References

1. Cho, J., et al., Cancers (Basel), 2020. 12(5):1137.
2. Lei, H., et al., Nat Commun., 2021. 12(1):51.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses  
Copyright © 2012 LifeSensors, Inc. All Rights Reserved