

USP47 (Ubiquitin specific protease 47)

Cat. # DB510

Background: USP47 deubiquitinates monoubiquitinated DNA polymerase β (PolB). PolB plays a role in DNA proof-reading via base-excision repair[1].

Alternate names: Deubiquitinating enzyme 47, Ubiquitin thioesterase 47, Ubiquitin carboxyl-terminal hydrolase 47

Product Information

Accession No.:	Q96K76
Molecular Weight:	157 kDa
Quantity:	25 μ g
Physical State:	Liquid, 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol
Source:	Human recombinant
Tag:	His ₆
Activity:	Active in the Ub-CHOP assay
Storage:	-80° C. Avoid repeated freeze/thaw cycles

SDS-PAGE analysis of USP47



References

1. Parsons, J.L., Dianova, I.I., Khoronenkova, S.V., Edelman, M.J., Kessler, B.M., and Dianova, G.L., USP47 is a deubiquitylating enzyme that regulates base excision repair by controlling steady-state levels of DNA Polymerase β . *Mol Cell*. **41**:609-615 (2011).

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