

USP10 (Ubiquitin Specific Protease 10)

Cat. # DB514

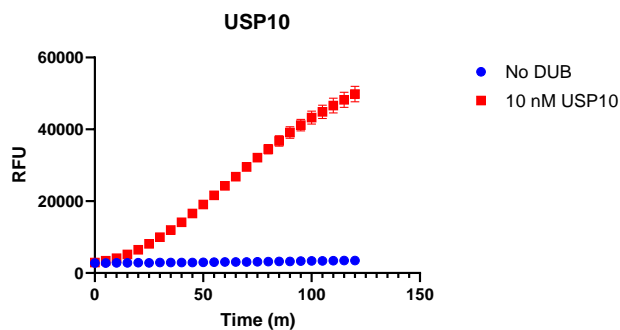
Background USP10 (Ubiquitin Specific Protease 10, UBPO) is a deubiquitinase that can remove conjugated ubiquitin from target proteins such as p53/TP53, BECN1, SNX3 and CFTR. In response to DNA damage, USP10 is translocated to the nucleus where it is involved in the deubiquitination of p53.

Alternate names Deubiquitinating enzyme 10, KIAA0190, MGC2621, Ubiquitin carboxyl-terminal hydrolase 10, Ubiquitin-specific-processing protease 10, Ubiquitin thioesterase 10, UBPO

Product Information

Accession No.	Q14694
Molecular Weight	99 kDa
Quantity	25 µg
Physical State	Liquid, 25 mM Tris-HCl, pH 7.5, 150 mM NaCl, 5 mM DTT, 10% glycerol
Source	Human Recombinant
Tag	His6-SUMO
Suggested Substrate	Ub-CHOP Reporter
Storage	-80° C. Avoid repeated freeze/thaw cycles

Product QC



Activity Assay of USP10. 10 nM USP10 was tested in a CHOP assay showing robust DUB activity.

References

1. Li, H., et al., Cell Rep., 2022. 41(1):111435.
2. Tao, L., et al., Genes (Basel), 2022. 13(5):831.

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