

ULP1 Antibody (Ubiquitin-like-specific protease 1)

Cat. # AB702

Background

Protease that specifically cleaves Smt3p protein conjugates; required for cell cycle progression; associates with nucleoporins and may interact with septin rings during telophase; sequestered to the nucleolus under stress conditions^{1,2,3,4}

Alternate Names: NIB1

Molecular Weight: 72kDa

Specificity/Applications

Species Cross Reactivity: Human

Source: Chicken

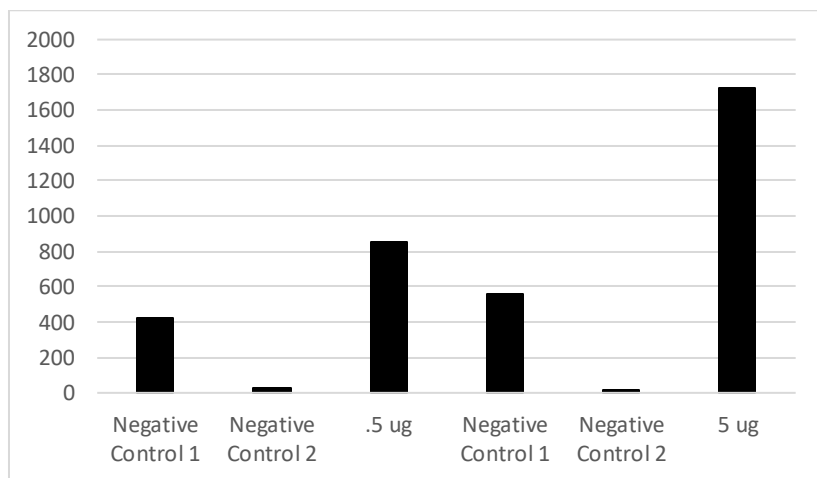
Applications: WB

Recommended Antibody Dilutions:

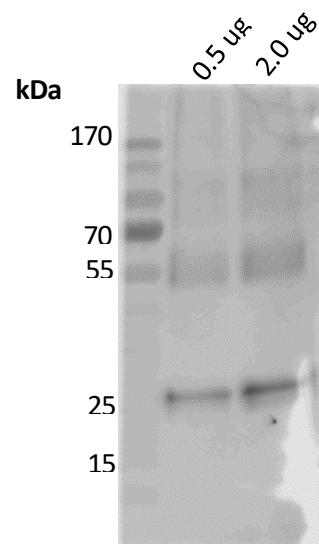
Western Blotting: Robust detection of 500ng of recombinant protein was possible when antibody was used at a final concentration of 5µg/mL

Storage: Store at -20°C; Avoid repeated freeze thaw. Supplied in phosphate buffered saline pH 7.4.

Detection of ULP1 by AB4010 using ELISA and Western Blot



0.5 or 5.0 ug of ULP1 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. ULP1 was detected by 1.54 ug/mL of AB703 using traditional ELISA detection reagent. 2° Antibody: α-Chicken HRP (1:5000). Negative Control1: No AB4010. Negative Control2: no detection reagent. 2° Antibody: α-Chicken HRP (1:5000).



Indicated amounts of ULP1 were loaded on SDS-PAGE gel followed by Western Blot. The blot was detected by 7.71 ug/mL of AB4010.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2007 LifeSensors, Inc. All Rights Reserved

References

- 1) Hochstrasser M (2001) SP-RING for SUMO: new functions bloom for a ubiquitin-like protein. Cell 107(1):5-8 PMID: 11595179
- 2) Li SJ and Hochstrasser M (2003) The Ulp1 SUMO isopeptidase: distinct domains required for viability, nuclear envelope localization, and substrate specificity. J Cell Biol 160(7):1069-81 PMID: 12654900
- 3) Soustelle C, et al. (2004) A new *Saccharomyces cerevisiae* strain with a mutant Smt3-deconjugating Ulp1 protein is affected in DNA replication and requires Srs2 and homologous recombination for its viability. Mol Cell Biol 24(12):5130-43 PMID: 15169880
- 4) Sydorsky Y, et al. (2010) A novel mechanism for SUMO system control: regulated Ulp1 nucleolar sequestration. Mol Cell Biol 30(18):4452-62 PMID: 20647537

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2007 LifeSensors, Inc. All Rights Reserved