

Chicken anti-SUMO/SUMOstar, affinity purified

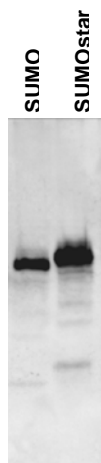
Cat. # AB7002

Background: Our Anti-SUMO antibody is affinity- purified, polyclonal chicken IgY. It can be used to localize yeast SUMO- and SUMOstar-tagged polypeptides by Western blot and ELISA. The antibody has not been tested for immunohistochemistry and immunoprecipitation applications.

Application: Western blot, recommended dilution 1:1000 to 1:5000

Product Information

Species:	Chicken
Isotype:	IgY
Purification:	Affinity purified
Physical State:	liquid
Applications:	WB (1:1000 to 1:5000)
Quantity:	50 µg
Storage:	For extended storage aliquot contents and freeze at -20°C or below. Avoid repeated freeze/thaw cycles



1:100 dilution of crude *E. coli* lysate expressing either SUMO-GFP or SUMOstar-GFP

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