

Anti-huSUMO3 antibody

Cat. # AB7030

Description

Our huSUMO3 antibody is affinity-purified, polyclonal chicken IgY. The antibody has not been tested for immunohistochemistry or immunoprecipitation applications.

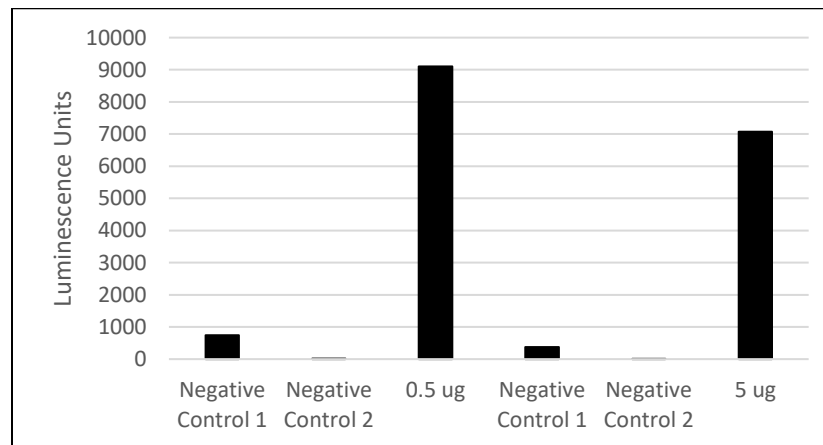
Product Information

Specificity:	HuSUMO3
Source:	Chicken, affinity purified IgY

Activity & Storage

Physical State:	Liquid
Concentration:	100 µL at 0.5 mg/mL
Buffer:	25 mM Tris-HCl, pH 7.3, 150 mM NaCl
Recommended Dilution:	1:1000 to 1:10000(Western)
Storage/Stability:	Store vial at -20°C. For extended storage, aliquot contents and freeze at -20°C or below. Avoid cycles of freezing and thawing. This product is stable for several weeks at 4°C. Make final dilution immediately prior to use.

Detection of Sumo3 by AB7030 using ELISA and Western Blot



0.5 or 5 ug of Sumo3 was coated on ELISA plate. Subsequently, unbound proteins were washed away and blocked with BSA. Sumo3 was detected by 2.51 µg/mL of AB7030 using traditional ELISA detection reagent. 2° Antibody: α-Chicken HRP (1:5000). Control 1: No AB7030
Negative Control 2: no detection reagent.

