

SODD (NT) antibody (pAb)

Rabbit Anti-Human/Mouse/Rat Silencer of Death Domains

Instruction Manual

Catalog Number	PK-AB718-2143
Synonyms	SODD Antibody:
Description	Apoptosis is induced by certain cytokines including TNF and Fas ligand of the TNF family through their death domain containing receptors, TNF-R1 and Fas. Several novel death receptors including DR3, DR4, DR5, and DR6 were recently identified. Cell death signal is transduced by death domain containing adapter molecules through the interaction with death domain of these death receptors. A novel TNF-R1 interacting protein was recently identified and designated SODD for silencer of death domains. SODD associates with the death domain of TNF-R1 and prevents constitutive activation of TNF-R1 signaling. TNF treatment releases SODD and permits adapter molecules such as TRADD recruiting to the active TNF-R1 complex, which activates TNF signaling pathways. SODD also interacts with DR3. SODD is ubiquitously expressed in human tissues and cell lines.
Quantity	100 µg
Source / Host	Rabbit
Immunogen	Rabbit polyclonal SODD antibody was raised against a peptide corresponding to amino acids 2 to 16 of human SODD.
Purification Method	Antibody is purified via DEAE-column chromatography.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human, Mouse, Rat
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, ICC, IF INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. SODD antibody can be used for detection of SODD by Western blot at 0.5 µg/mL. An approximately 60 kDa band can be detected. Antibody can also be used for immunocytochemistry starting at 5 µg/mL. For immunofluorescence start at 20 µg/mL.
Images	Available upon request.
References	Jiang Y, Woronicz JD, Liu W, Goeddel DY. Prevention of constitutive TNF receptor 1 signaling by silencer of death domains. Science 1999;283:543-6 (RD1299)
Images	Available upon request.
Related Products	DR3 Antibody, Cat. No. PK-AB718-1158 DR4 Antibody, Cat. No. PK-AB718-1139 DR5 Antibody, Cat. No. PK-AB718-2019 HeLa Lysate, Cat. No. PK-AB718-1201 SODD Peptide, Cat. No. PK-AB718-2143P

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.