

Instruction Manual

Catalog Number	PK-AB718-2143P
Quantity	50 µg
Source	15 amino amino terminus of human SODD
Formulation	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH ₂ PO ₄ , 10 mM Na ₂ HPO ₄ , 130 mM NaCl) containing 0.1% bovine serum albumin and 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	Store SODD peptide at -20°C, stable for one year.
Application	SODD peptide is used for blocking the activity of SODD antibody.
References	<p>Jiang Y, Woronicz JD, Liu W, Goeddel DY. Prevention of constitutive TNF receptor 1 signaling by silencer of death domains. <i>Science</i> 1999;283:543-6.</p> <p>Wesche H, Korherr C, Kracht M, Falk W, Resch K, Martin MU. The interleukin-1 receptor accessory protein (IL-1RAcP) is essential for IL-1-induced activation of interleukin-1 receptor-associated kinase (IRAK) and stress-activated protein kinases (SAP kinases). <i>J Biol Chem</i> 1997;272:7727-31</p> <p>Hofmeister R, Wiegmann K, Korherr C, Bernardo K, Kronke M, Falk W. Activation of acid sphingomyelinase by interleukin-1 (IL-1) requires the IL-1 receptor accessory protein. <i>J Biol Chem</i> 1997;272:27730-6 (RD1299)</p>

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