

Instruction Manual

Catalog Number	PK-AB718-3679P
Quantity	50 µg
Source	16 amino acids near the amino terminus of human ASK1
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 ASK1 peptide at -20°C, stable for one year.
Application	ASK1 peptide is used for blocking the activity of ASK1 antibody.
References	<p>Wang XS, Diener K, Jannuzzi D, et al. Molecular cloning and characterization of a novel protein kinase with a catalytic domain homologous to mitogen-activated protein kinase kinase kinase. <i>J. Biol. Chem.</i> 1996; 271:31607-11.</p> <p>Grech AP, Amesbury M, Chan T, et al. TRAF2 differentially regulates the canonical and noncanonical pathways of NF-kappaB activation in mature B cells. <i>Immunity</i> 2004; 21:629-42.</p> <p>Davies CC, Mak TW, Young LS, et al. TRAF6 is required for TRAF2-dependent CD40 signal transduction in nonhemopoietic cells. <i>Mol. Cell. Biol.</i> 2005; 25:9806-19. 5. Kuai J, Nickbarg E, Wooters J, et al. Endogenous association of TRAF2, TRAF3, cIAP1, and Smac with lymphotoxin-beta receptor reveals a novel mechanism of apoptosis. <i>J. Biol. Chem.</i> 2003; 278:14363-9.</p>

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