

Instruction Manual

Catalog Number	PK-AB718-1151P
Quantity	50 µg
Source	20 amino acids near the carboxy 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>Tobiume K, Inage T, Takeda K, et al. Molecular cloning and characterization of the mouse apoptosis signal-regulating kinase 1. <i>Biochem. Biophys. Res. Commun.</i> 1997; 239:905-10.</p> <p>Wolf BB and Green DR. Suicidal tendencies: apoptotic cell death by caspase family proteinases. <i>J. Biol. Chem.</i> 1999; 274:20049-52.</p> <p>Muzio M, Chinnaiyan AM, Kischkel FC, et al. FLICE, a novel FADD-homologous ICE/CED-3-like protease, is recruited to the CD95 (Fas/APO-1) death-inducing signaling complex. <i>Cell</i> 1996; 85:817-27.</p>

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