

Instruction Manual

Catalog Number	PK-AB718-3903P
Quantity	50 µg
Source	14 amino acids near the amino terminus of human SHP2
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 SHP2 peptide at -20°C, stable for one year.
Application	SHP2 peptide is used for blocking activity of SHP2 antibody.
References	<p>Yu Z, Ahmad S, Schwartz JL, et al. Protein-tyrosine phosphatase SHP2 is positively linked to proteinase-activated receptor 2-mediated mitogenic pathway. <i>J. Biol. Chem.</i> 1997; 272:7519-24</p> <p>Chariot A, Leonardi A, Muller J, et al. Association of the adaptor TANK with the IκappaB kinase (IKK) regulator NEMO connects IKK complexes with the IKKepsilon and TBK1 kinases. <i>J. Biol. Chem.</i> 2002; 277:37029-36</p>

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