

Instruction Manual

Catalog Number	PK-AB718-2383P
Quantity	50 µg
Source	16 amino amino terminus of human p53R2
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 p53R2 peptide at -20°C, stable for one year.
Application	p53R2 peptide is used for blocking the activity of p53R2 antibody.
References	<p>Matsuoka S, Huang M, Elledge SJ. Linkage of ATM to cell cycle regulation by the Chk2 protein kinase. <i>Science</i>. 1998;282:1893-7.</p> <p>Li Y, Kang J, Friedman J, Tarassishin L, Ye J, Kovalenko A, Wallach D, Horwitz MS. Identification of a cell protein (FIP-3) as a modulator of NF-kappaB activity and as a target of an adenovirus inhibitor of tumor necrosis factor alpha-induced apoptosis. <i>Proc Natl Acad Sci U S A</i> 1999;96(3):1042-7</p> <p>Mercurio F, Murray BW, Shevchenko A, Bennett BL, Young DB, Li JW, Pascual G, Motiwala A, Zhu H, Mann M, Manning AM. IkappaB kinase (IKK)-associated protein 1, a common component of the heterogeneous IKK complex. <i>Mol Cell Biol</i>. 1999;19(2):1526-38.</p>

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