

Instruction Manual

Catalog Number	PK-AB718-3041P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human PUMA
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 PUMA peptide at -20°C, stable for one year.
Application	PUMA peptide is used for blocking the activity of PUMA antibody.
References	<p>Nakano K, Vousden KH. PUMA, a novel proapoptotic gene, is induced by p53. <i>Mol Cell.</i> 2001;7(3):683-94</p> <p>Fischer KD, Tedford K, Wirth T. New roles for Bcl10 in B-cell development and LPS response. <i>Trends Immunol.</i> 2004; 25:113-6.</p> <p>Stilo R, Liguoro D, Di Jeso B, et al. Physical and functional interaction of CARMA1 and CARMA3 with Ikappa kinase kappa-NF-kappaB essential modulator. <i>J. Biol. Chem.</i> 2004; 279:34323-31.</p>

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