

Instruction Manual

Catalog Number	PK-AB718-3043P
Quantity	50 µg
Source	15 amino amino 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	Nakano K, Vousden KH. PUMA, a novel proapoptotic gene, is induced by p53. <i>Mol Cell.</i> 2001;7(3):683-94. Robinson DN and Cooley L. Drosophila kelch is an oligomeric ring canal actin organizer. <i>J. Cell Biol.</i> 1997; 138:799-810.

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