

Instruction Manual

Catalog Number	PK-AB718-3411P
Quantity	50 µg
Source	16 amino acids near the middle of Anthrax protective antigen
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 Anthrax PA peptide at -20°C, stable for one year.
Application	Anthrax peptide is used for blocking the activity of Anthrax PA antibody.
References	Welkos SL, Lowe JR, Eden-McCutchan F, et al. Sequence and analysis of the DNA encoding protective antigen of Bacillus anthracis. <i>Gene</i> 1988; 69:287-300. Heiser D, Labi V, Erlacher M, et al. The Bcl-2 protein family and its role in the development of neoplastic disease. <i>Exp. Gerontol.</i> 2004; 39:1125-35. Oltvai ZN, Millman CL, and Korsmeyer SJ. Bcl-2 heterodimerizes in vivo with a conserved homolog, Bax, that accelerates programmed cell death. <i>Cell</i> 1993; 74:609-19. Wang K, Yin XM, Chao DT, et al. BID: a novel BH3 domain-only death agonist. <i>Genes Dev.</i> 1996; 10:2859-69.

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