

Instruction Manual

Catalog Number	PK-AB718-3415P
Quantity	50 µg
Source	16 amino acids in the middle of Anthrax lethal factor
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 Lethal Factor peptide at -20°C, stable for one year.
Application	Anthrax peptide is used for blocking the activity of Anthrax lethal factor antibody.
References	<p>Bragg TS and Robertson DL. Nucleotide sequence and analysis of the lethal factor gene (lef) from <i>Bacillus anthracis</i>. <i>Gene</i> 1989; 81:45-54.</p> <p>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.</p> <p>Wang K, Yin XM, Chao DT, et al. BID: a novel BH3 domain-only death agonist. <i>Genes Dev.</i> 1996; 10:2859-69.</p> <p>YI J-S, Lee, S-K, Sato T-A, et al. Co-induction of p75NTR and the associated death executor NADE in degenerating hippocampal neurons after kainate-induced seizures in the rat. <i>Neurosci. Lett.</i> 2003; 347:126-30.</p>

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