

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3417P
<b>Quantity</b>	50 µg
<b>Source</b>	16 amino acids near the carboxy terminus of Anthrax lethal factor
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	Store Anthrax Lethal Factor peptide at -20°C, stable for one year.
<b>Application</b>	Anthrax peptide is used for blocking the activity of Anthrax lethal factor antibody.
<b>References</b>	<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>Rapp G, Freudenstein J, Klaudiny J, et al. Characterization of three abundant mRNAs from human ovarian granulosa cells. <i>DNA Cell Biol.</i> 1990; 9:479-85.</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> <p>Sweet MJ, Leung BP, Kang D, et al. A novel pathway regulating lipopolysaccharide-induced shock by ST2/T1 via inhibition of Toll-like receptor 4 expression. <i>J. Immunol.</i> 2001; 166:6633-9.</p>

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