

Instruction Manual

Catalog Number	PK-AB718-4031P
Quantity	50 µg
Source	19 amino acid peptide from near the amino terminus of human Emerin
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 Emerin peptide at -20°C, stable for one year.
Application	Emerin peptide is used for blocking the activity of Emerin antibody.
References	Schirmer EC, Florens L, Guan T, et al. Nuclear membrane proteins with potential disease links found by subtractive proteomics. Science 2003; 301:1380-2.

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