

Instruction Manual

Catalog Number	PK-AB718-4247P
Quantity	50 µg
Source	24 amino acids near the center of human IEX-1
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 IEX-1 peptide at -20°C, stable for one year.
Application	IEX-1 peptide is used for blocking the activity of IEX-1 antibody.
References	Kondratyev AD, Chung K-N and Jung MO. Identification and characterization of a radiation-inducible glycosylated human early-response gene. Cancer Res. 1996; 56:1498-502.

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