

Instruction Manual

Catalog Number	PK-AB718-4641P
Quantity	50 µg
Source	16 amino acids near the amino terminus of HIV-1 gp120
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 gp120 peptide at -20°C, stable for one year.
Application	gp120 peptide is used for blocking the activity of gp120 antibody.
References	Pinter A. Roles of HIV-1 Env variable regions in viral neutralization and vaccine development. Curr. HIV Res. 2007; 5:542-53.

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