

Instruction Manual

Catalog Number	PK-AB718-3257P
Quantity	50 µg
Source	15 amino acids near the amino terminus of human APOBEC3G
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 APOBEC3G peptide at -20°C, stable for one year.
Application	The peptide is used for blocking the activity of APOBEC3G antibody.
References	Jarmuz A, Chester A, Bayliss J, et al. An anthropoid-specific locus of Orphan C to U RNA-editing enzymes on chromosome 22. <i>Genomics</i> 2002; 79:285-96. Nishigaki K, Thompson D, Yugawa T, et al. (2003) Identification and characterization of a novel Ste20/germinal center kinase-related kinase, polyploidy-associated protein kinase. <i>J. Biol. Chem.</i> 278:13520-30.

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