

Instruction Manual

Catalog Number	PK-AB718-4205P
Quantity	50 µg
Source	19 amino acids near the center of human HAP1
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 HAP1 peptide at -20°C, stable for one year.
Application	HAP1 peptide is used for blocking the activity of HAP1 antibody.
References	Li X-J, L S-H, Sharp AH, et al. A huntingtin-associated protein enriched in brain with implications for pathology. Nature 1995; 378:398-402.

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