

Instruction Manual

Catalog Number	PK-AB718-1133P
Quantity	50 µg
Source	14 amino acids near the center of human GFR alpha 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 GFR alpha 1 peptide at -20°C, stable for one year.
Application	GFR alpha 1 peptide is used for blocking the activity of GFR alpha 1 antibody.
References	Jing S; Wen D; Yu Y; Holst PL ; Luo Y; Fang M; Tamir R; Antonio L; Hu Z; Cupples R; Louis JC; Hu S; Altrock BW; Fox GM. GDNF-induced activation of the Ret protein tyrosien kinase is mediated by GDNFR-alpha, a novel receptor for GDNF. Cell, 1996;85:1113-24. Cohen GM. Caspases: the executioners of apoptosis. Biochem J 1997;326:1-16 (RD1299)

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