

Instruction Manual

Catalog Number	PK-AB718-1137P
Quantity	50 µg
Source	14 amino acids near the center of human GFR alpha 3
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 3 peptide at -20°C, stable for one year.
Application	GFR alpha 3 peptide is used for blocking the activity of GFR alpha 3 antibody.
References	<p>Naveilhan P, Baudet C, Mikaels A, Shen L, Westphal H, Ernfors P. Expression and regulation of GFRalpha3, a glial cell line-derived neurotrophic factor family receptor. Proc Natl Acad Sci USA 1998;95:1295-1300.</p> <p>Woronicz JD, Gao X, Cao Z, Rothe M, Goeddel DV. IkappaB kinase-beta: NF-kappaB activation and complex formation with IkappaB kinase-alpha and NIK. Science 1997;278:866-869</p> <p>Ling L, Cao Z, Goeddel D. NF-kappaB-inducing kinase activates IKK-alpha by phosphorylation of Ser-176. Proc Natl Acad Sci USA 1998;95:3792-3797</p>

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