

Instruction Manual

Catalog Number	PK-AB718-1135P
Quantity	50 µg
Source	15 amino acids near the center of human GFR- 2
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 2 peptide at -20°C, stable for one year.
Application	GFR alpha 2 peptide is used for blocking the activity of GFR alpha 2 antibody.
References	<p>Klein RD, Sherman D, Ho WH, et al. A GPI-linked protein that interacts with Ret to form a candidate neurturin receptor. <i>Nature</i> 1997; 387:717-21.</p> <p>Wolf BB and Green DR. Suicidal tendencies: apoptotic cell death by caspase family proteinases. <i>J. Biol. Chem.</i> 1999; 274:20049-52.</p> <p>Muzio M, Chinnaiyan AM, Kischkel FC, et al. FLICE, a novel FADD-homologous ICE/CED-3-like protease, is recruited to the CD95 (Fas/APO-1) death-inducing signaling complex. <i>Cell</i> 1996; 85:817-27.</p>

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