

Instruction Manual

| | |
|--------------------------------|---|
| Catalog Number | PK-AB718-3553P |
| Quantity | 50 µg |
| Source | 15 amino acids near the amino terminus of human adiponectin |
| 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 Adiponectin peptide at -20°C, stable for one year. |
| Application | Adiponectin peptide is used for blocking the activity of the adiponectin antibody. |
| References | <p>Tsao T-S, Lodish HF, and Fruebis J. ACRP30, a new hormone controlling fat and glucose metabolism. <i>Euro. J. Pharmacol.</i> 2002; 440:213-21.</p> <p>Inoki K, Li Y, Xu T, et al. Rheb GTPase is a direct target of TSC2 GAP activity and regulates mTOR signaling. <i>Genes Dev.</i> 2003; 17:1829-34.</p> <p>Stocker H, Radimerski T, Schindelholz B, et al. Rheb is an essential regulator of S6K in controlling cell growth in <i>Drosophila</i>. <i>Nat. Cell Biol.</i> 2003; 5:559-65.</p> |

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