

Instruction Manual

Catalog Number	PK-AB718-3521P
Quantity	50 µg
Source	16 amino acids near the center of human IRS-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 IRS-1 peptide at -20°C, stable for one year.
Application	IRS-1peptide is used for blocking the activity of IRS-1 antibody.
References	<p>Giovannone B, Scaldaferrri ML, Federici M, et al. Insulin receptor substrate (IRS) transduction system: distinct and overlapping signaling potential. <i>Diabetes Metab. Res. Rev.</i> 2000; 16:434-41.</p> <p>McGettrick AF and O'Neill LAJ. The expanding family of MyD88-like adaptors in Toll-like receptor signal transduction. <i>Mol. Imm.</i> 2004; 41:577-82.</p> <p>Zhang G and Ghosh S. Negative regulation of Toll-like receptor-mediated signaling by Tollip. <i>J. Biol. Chem.</i> 2002; 277:7059-65.</p> <p>Kozopas KM, Yang T, Buchan HL, et al. MCL1, a gene expressed in programmed myeloid cell differentiation, has sequence similarity to BCL2. <i>Proc. Natl. Acad. Sci. USA</i> 1993; 90:3516-20.</p>

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