

Instruction Manual

Catalog Number	PK-AB718-3513P
Quantity	50 µg
Source	14 amino acids near the carboxy terminus of human 4E-BP1
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 4E-BP1 peptide at -20°C, stable for one year.
Application	4E-BP1 peptide is used for blocking the activity of 4E-BP1 antibody.
References	<p>Shah OJ, Anthony JC, Kimball SR, et al. 4E-BP1 and S6K1: translational integration sites for nutritional and hormonal information in muscle. <i>Am. J. Physiol. Endocrinol. Metab.</i> 2000; 279:E715-29.</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>Fernandes-Alnemri T, Litwack G, and Alnemri ES. Mch2, a new member of the apoptotic Ced-3/Ice cysteine protease gene family. <i>Cancer Res.</i> 1995; 55:2737-42.</p> <p>Nishiya T and DeFranco AL. Ligand-regulated chimeric receptor approach reveals distinctive subcellular localization and signaling properties of the Toll-like receptors. <i>J. Biol. Chem.</i> 2004; 279:19008-17.</p>

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