

Instruction Manual

Catalog Number	PK-AB718-3399P
Quantity	50 µg
Source	14 amino acids near the center of human IRF3
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 IRF3 peptide at -20°C, stable for one year.
Application	IRF3 peptide is used for blocking the activity of IRF3 antibody.
References	<p>Au WC, Moore PA, Lowther W, et al. Identification of a member of the interferon regulatory factor family that binds to the interferon-stimulated response element and activates expression of interferon-induced genes. <i>Proc. Natl. Acad. Sci. USA</i> 1995; 92:11657-61.</p> <p>Heiser D, Labi V, Erlacher M, et al. The Bcl-2 protein family and its role in the development of neoplastic disease. <i>Exp. Gerontol.</i> 2004; 39:1125-35.</p> <p>Hockenbery D, Nunez G, Milliman C, et al. Bcl-2 is an inner mitochondrial membrane protein that blocks programmed cell death. <i>Nature</i> 1990; 348:334-6.</p> <p>Gonzalez-Garcia M, Perez-Ballestro R, Ding L et al. bcl-xL is the major bcl-x mRNA form expressed during murine development and its product localizes to mitochondria. <i>Development</i> 1994; 120:3033-42.</p>

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