

Instruction Manual

Catalog Number	PK-AB718-3659P
Quantity	50 µg
Source	15 amino acids near the center of human IRE1p
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 IRE1p peptide at -20°C, stable for one year.
Application	IRE1p peptide is used for blocking the activity of IRE1p antibody.
References	<p>Miyoshi K, Katayama T, Imaizumi K, et al. Characterization of mouse Ire1 alpha: cloning, mRNA localization in the brain and functional analysis in a neural cell line. <i>Brain Res. Mol. Brain Res.</i> 2000; 85:68-76.</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>Hayashi F, Smith KD, Ozinsky A, et al. The innate immune response to bacterial flagellin is mediated by Toll-like receptor 5. <i>Nature</i> 2001; 410:1099-103.</p> <p>Hawn TR, Verbon A, Lettinga KD, et al. A common dominant TLR5 stop codon polymorphism abolishes flagellin signaling and is associated with susceptibility to legionnaires' disease. <i>J. Exp. Med.</i> 2003; 198:1563-72.</p>

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