

## Instruction Manual

<b>Catalog Number</b>	PK-AB718-3119P
<b>Quantity</b>	50 µg
<b>Source</b>	13 amino acids near the C-terminus of human ATR
<b>Formulation</b>	Peptide is supplied as 200 µg/ml solution in PBS pH 7.2 (10 mM NaH <sub>2</sub> PO <sub>4</sub> , 10 mM Na <sub>2</sub> HPO <sub>4</sub> , 130 mM NaCl) containing 0.1% bovine serum albumin and 0.02% sodium azide.
<b>Reconstitution</b>	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.
<b>Storage &amp; Stability</b>	Store ATR / TEM8 peptide at -20°C, stable for one year.
<b>Application</b>	ATR peptide is used for blocking the activity of ATR antibody.
<b>References</b>	<p>Bradley KA, Mogridge J, Mourez M, et al. Identification of the cellular receptor for anthrax toxin. <i>Nature</i> 2001; 414:225-9.</p> <p>Hlaing T, Guo RF, Dilley KA, et al. Molecular cloning and characterization of DEFCAP-L and -S, two isoforms of a novel member of the Ced-4 family of apoptosis proteins. <i>J. Biol. Chem.</i> 2001; 276:9230-8.</p> <p>Sanz C, Calasanz MJ, Andreu E, et al. NALP1 is a transcriptional target for cAMP-response-element-binding protein (CREB) in myeloid leukemia cells. <i>Biochem. J.</i> 2004; 384:281-6.</p>

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