

Instruction Manual

Catalog Number	PK-AB718-3525P
Quantity	50 µg
Source	20 amino acids near the carboxy terminus of SARS Spike protein
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 SARS Spike peptide at -20°C, stable for one year.
Application	SARS Spike peptide is used for blocking the activity of the SARS Spike antibody.
References	<p>Marra MA, Jones SJ, Astell CR, et al. The Genome sequence of the SARS-associated corona virus. <i>Science</i> 2003;300:1399-404.</p> <p>Opferman JT, Letai A, Beard C, et al. Development and maintenance of B and T lymphocytes require antiapoptotic MCL-1. <i>Nature</i> 2003; 426:671-6.</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> <p>Fingar DC and Blenis J. Target of rapamycin (TOR): an integrator of nutrient and growth factor signals and coordinator of cell growth and cell cycle progression. <i>Oncogene</i> 2004; 23:3151-71.</p>

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