

Instruction Manual

Catalog Number	PK-AB718-3529P
Quantity	50 µg
Source	13 amino acids near the carboxy terminus of SARS M 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 Matrix peptide at -20°C, stable for one year.
Application	SARS matrix peptide is used for blocking the activity of the SARS matrix 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>Yonezawa K, Tokunaga C, Oshiro N, et al. Raptor, a binding partner of target of rapamycin. <i>Biochem. Biophys. Res. Commun.</i> 2004; 313:437-441.</p> <p>Hara K, Yonezawa K, Weng QP, et al. Amino acid sufficiency and mTOR regulate p70 S6 kinase and eIF-4E BP1 through a common effector mechanism. <i>J. Biol. Chem.</i> 1998; 273:14484-94.</p>

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