

SARS E (NT) antibody (pAb)

Rabbit Anti-SARS Virus Envelope Protein

Instruction Manual

| | |
|---------------------|--|
| Catalog Number | PK-AB718-3531 |
| Synonyms | SARS Envelope Antibody: SARS |
| Description | A novel coronavirus has recently been identified as the causative agent of SARS (Severe Acute Respiratory Syndrome). Coronaviruses are a major cause of upper respiratory diseases in humans. The genomes of these viruses are positive-stranded RNA approximately 27-31kb in length. SARS infection can be mediated by the binding of the viral spike protein, a glycosylated 139 kDa protein and the major surface antigen of the virus, to the angiotensin-converting enzyme 2 (ACE2) on target cells. This binding can be blocked by a soluble form of ACE2. |
| Quantity | 100 µg |
| Source / Host | Rabbit |
| Immunogen | Rabbit polyclonal SARS E antibody was raised against a synthetic peptide corresponding to amino acids at the amino-terminus of the SARS E protein (Genbank accession no. P59637). |
| Purification Method | Affinity chromatography purified via peptide column. |
| Clone / IgG Subtype | Polyclonal antibody |
| Species Reactivity | Virus |
| Specificity | |
| Formulation | Antibody is supplied in PBS containing 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 | Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures. |
| Applications | E Note: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. SARS envelope antibody can be used for the detection of SARS envelope protein in ELISA. It will detect 10 ng of free peptide at 1 µg/mL. |
| Images | NA |
| References | Marra MA, Jones SJ, Astell CR, et al. The Genome sequence of the SARS-associated corona virus. Science 2003;300:1399-404. Rota PA, Oberste MS, Monroe SS, et al. Characterization of a novel coronavirus associated with severe acute respiratory syndrome. Science 2003;300:1394-9. Navas-Nartin SR and Weiss S. Coronavirus replication and pathogenesis: Implications for the recent outbreak of severe acute respiratory syndrome (SARS), and the challenge for vaccine development. J Neurovirol. 2004;10:75-85. Arbely E, Khattari Z, Brotons G, et al. A highly unusual palindromic transmembrane helical hairpin formed by SARS coronavirus E protein. J Mol. Biol. 2004;3414:769-79. |
| Images | NA |
| Related Products | Blocking Peptide, Cat. No. PK-AB718-3531P SARS E protein Antibody (CT), Cat. No. PK-AB718-3533; SARS Spike Antibody (NT), Cat. No. PK-AB718-3219 SARS Spike Antibody (IN1), Cat. No. PK-AB718-3221; SARS M protein Antibody (NT), Catalog No. PK-AB718-3529 SARS Spike Antibody (IN2), Cat. No. PK-AB718-3223; SARS Spike Antibody (IN3), Cat. No. PK-AB718-3225 |
| FOR IN VITRO RES | SARS Spike Antibody (CT), Cat. No. PK-AB718-3525, SARS M protein Antibody (CT), Cat. No. PK-AB718-3527 |