

## SARS-ACE (residues 1-76) SARS-Associated Coronavirus Envelope recombinant, *E. coli*

Cat. No.	Amount
PR-1100	100 $\mu$ g

For *in vitro* use only  
Quality guaranteed for 12 months  
Store at -20°C

### Avoid freeze / thaw cycles

### Form

Liquid. Supplied in 50 mM Tris-HCl, 60 mM NaCl and 50% glycerol.

### Specificity

Immunoreactive with sera of SARS infected individuals.

### Application

Recombinant SARS-ACE Antigen may be used in ELISA and Western blots, excellent for detection of SARS with minimal specificity problems.

### Molecular Weight

43 kDa

### Purity

>95% by SDS-PAGE

### Description

SARS-ACE contains the N-terminus Envelope protein immunodominant regions, amino acids: 1-76. SARS-ACE is purified by proprietary chromatographic techniques.

### Background

SARS (Severe Acute Respiratory Syndrome) Coronavirus is an enveloped virus containing three outer structural proteins, namely the membrane (M), envelope (E), and spike (S) proteins.

Spike (S)-glycoprotein of the virus interacts with a cellular receptor and mediates membrane fusion to allow viral entry into susceptible target cells.

Accordingly, S-protein plays an important role in virus infection cycle and is the primary target of neutralizing antibodies.

### Selected References:

- Liao *et al.* (2004) Expression of SARS-coronavirus envelope protein in *Escherichia coli* cells alters membrane permeability. *Biochem. Biophys. Res Commun.* **325**:374.  
Shen *et al.* (2003) Small envelope protein E of SARS: cloning, expression, purification, CD determination, and bioinformatics analysis. *Acta Pharmacol. Sin.* **24**:505.