

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

Cat. No.	Amount
PR-1100	100 μ 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 SARSinfected individuals.

Application

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

Purity

>95% by SDS-PAGE (coomassie staining) and RP-HPLC.

Description

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