

**SARS-ACE (residues 1-76)**

SARS-Associated Coronavirus Envelope
recombinant, *E. coli*

Cat. No.	Amount
PR-1100	100 µg

For in vitro use only!

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Molecular Weight: 34 kDa

Purity: > 95 % (SDS-PAGE)

Form: liquid (Supplied in 50 mM Tris-HCl, 60 mM NaCl and 50% glycerol)

Applications:

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

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.

Specificity: Immunoreactive with sera of SARS infected individuals.

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.