
Spike Glycoprotein Polyclonal Antibody

(Catalog # A56214)

Background

S1 attaches the virion to the cell membrane by interacting with sialic acid-containing cell receptors, initiating the infection. S2 is a class I viral fusion protein. Under the current model, the protein has at least 3 conformational states: pre-fusion native state, pre-hairpin intermediate state, and post-fusion hairpin state. During viral and target cell membrane fusion, the coiled coil regions (heptad repeats) assume a trimer-of-hairpins structure, positioning the fusion peptide in close proximity to the C-terminal region of the ectodomain. The formation of this structure appears to drive apposition and subsequent fusion of viral and target cell membranes.

Description

Spike Glycoprotein Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human coronavirus OC43

Isotype

IgG

Uniprot ID

P36334

Purification

>95%, Protein G purified

Immunogen

Recombinant Human coronavirus OC43 Spike glycoprotein (15-344AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

E2, Peplomer protein, S, 3

Application

ELISA
