

SARS 3CL Protease Antibody (Chymotrypsin-like Protease)

Cat. # CV4002

Background

Within the last two decades, SARS and MERS coronaviruses emerged as global health concerns causing severe acute respiratory syndromes. In December 2019, a novel coronavirus (SARS-CoV-2) was identified in Wuhan, Hubei province in China. The SARS-CoV genome encodes several non-structural proteins that are essential for the life cycle of the virus. The viral 3-chymotrypsin-like cysteine protease (3CL-PRO) enzyme controls coronavirus replication. 3CL-PRO is a proven drug discovery target in the case of SARS-CoV and MERS-CoV. Since the genome sequence of SARS-CoV-2 is very similar to that of SARS-CoV, inhibition of 3CL-PRO activity represents a viable strategy at the basis of a COVID-19 therapy. The SARS 3CL-PRO Protein Antibody recognizes and binds to both SARS-CoV and SARS-CoV-2 3CL-PRO enzymes.

Target Molecular Weight: 34 kDa

Product Information

Description: Rabbit, polyclonal antibody to SARS 3CL Protease

Species Cross Reactivity: SARS-CoV, SARS-CoV-2

Source: Rabbit

Applications: WB, ELISA

Recommended Antibody Dilutions:

Western Blotting: Robust detection of 10 ng of recombinant protein was possible when antibody was used at a final concentration of 1 µg/mL

Storage/Purification

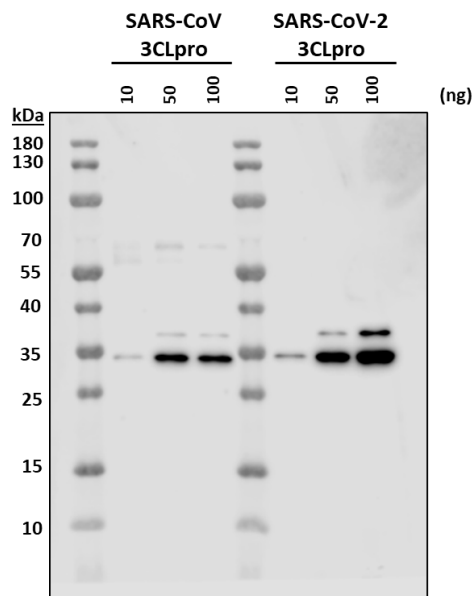
Polyclonal antibodies are produced by repeatedly immunizing rabbits with purified recombinant full-length protein. Antibodies are purified from monospecific antiserum by protein A affinity purification.

Storage: Supplied in 0.02 M Potassium Phosphate, 0.15 M Sodium Chloride pH 7.2.

Store at -20°C.

Avoid Freeze/Thaw Cycles.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
Copyright © 2007 LifeSensors, Inc. All Rights Reserved



References

- 1) Wang D, Hu B, Hu C, et al. Clinical Characteristics of 138 Hospitalized Patients With 2019 Novel Coronavirus–Infected Pneumonia in Wuhan, China. *JAMA*. 2020;323(11):1061.
- 2) Zhou P, Yang X-L, Wang X-G, et al. A pneumonia outbreak associated with a new coronavirus of probable bat origin. *Nature*. 2020;579(7798):270–273.
- 3) Zhu N, Zhang D, Wang W, et al. A Novel Coronavirus from Patients with Pneumonia in China, 2019. *N. Engl. J. Med.* 2020;382(8):727–733.

All products are for research use only • Not intended for human or animal diagnostic or therapeutic uses
 Copyright © 2007 LifeSensors, Inc. All Rights Reserved