

Data Sheet

3CL Protease Antibody (SARS-CoV)

Rabbit, Polyclonal
Catalog #: 100865
Lot #: 200820

Host Species | Isotype:

Rabbit | IgG

Formulated in: 0.02 M Potassium Phosphate,
 0.15 M NaCl pH 7.2.

Stability: At least 6 months at -20°C. *Avoid freeze/thaw cycles.*

References:

1. Wang, D., *et al.*, *JAMA*. 2020; **323(11)**: 1061-1069.
2. Zhou, P., *et al.*, *Nature*. 2020; **579(7798)**: 270-273.

Description: Rabbit polyclonal antibody recognizing the SARS-CoV 3CL Protease (MPro) as a 34 kDa protein. This antibody recognizes 3CL protease in both SARS-CoV and SARS-CoV-2.

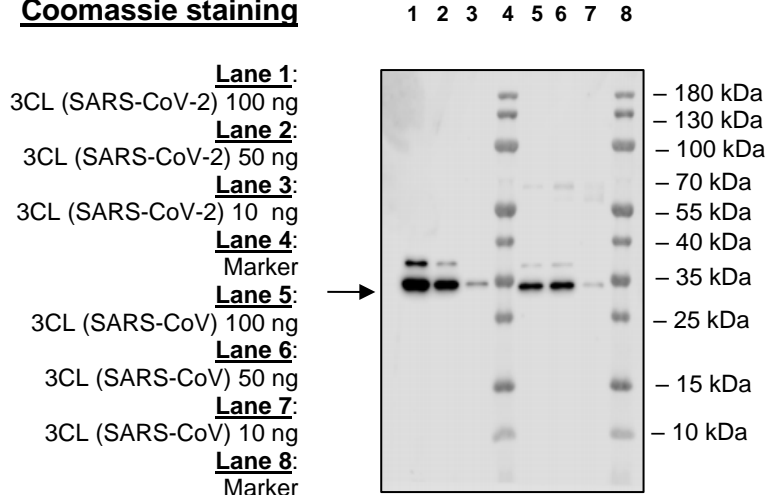
Immunogen: Purified, full-length 3CL protease.

Purification: Purified using protein A resin.

Applications: Useful in ELISA and Western blotting assays (1 µg/ml) for detecting SARS 3CL protease.

Quality Assurance

4-20% SDS-PAGE
Coomassie staining



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com