

2019-nCoV High Sensitivity S1 Protein ELISA Kit

Catalog # C19S1-876

Product Description

The 2019-nCoV High Sensitivity S1 Protein detection Kit is a sandwich Enzyme-Linked ImmunoSorbent Assay (ELISA) kit for the measurement and detection of Coronavirus Spike Protein in complex biological samples. The kit includes one 96-well strip plate pre-coated with a capture anti-S1 protein antibody. A calibration standard is included to allow quantitative determinations of SARS-CoV-2 Spike Protein S1 levels in test samples.

Components – Kit Box

SignalChem Cat#	Component Name	Size
C19S1-HS61DH	Anti-nCoV S1 Detection Antibody, HS-1X	10 mL
CP06-642HL	Anti-2019-nCoV Spike Protein Coated Plate	1 plate
CS02-506	One-step TMB Substrate	10 mL
E01-09	nCoV ELISA Sample Dilution Buffer	2 x 10 mL
SS01-09	Stop Solution	6 mL
C19SD-CAL50	Calibrator: 50 ng/mL	1 mL
WB20-09	Wash Buffer, 20X	2 x 15 mL

Storage and Stability

Store kit and contents at 4°C for up to 6 months. Store unused multi well strips in the provided resealable bag and use within one month after opening.

Other Materials Required

- Micro plate cover
- Milli-Q water or other source of pure water
- Pipettes or multichannel pipettes
- Incubator for 37°C incubation
- Microplate reader capable of measuring absorbance at 450nm

Short Protocol

The following is only a short protocol. For a more detailed protocol, see 2019-nCoV High Sensitivity S1 Protein ELISA Kit Instruction Manual.

2019-nCoV S1 Protein ELISA Assay Protocol:

- Bring all kit reagents to room temperature before use.
- Prepare 1X Wash buffer by adding 30 mL of the 20X concentrate with 570 mL of Milli-Q water.
- Prepare optimal sample dilutions using sample dilution buffer.
- Place the required number of coated strips onto the ELISA plate frame
- Dilute the calibrator with sample dilution buffer to 0 ng/mL (blank), 0.8 ng/mL, 2 ng/mL, and 5 ng/mL.
- Add 100µL each of Sample or Calibrators per well (in duplicate or triplicate).
- Cover plate and incubate 30 minutes at 37°C.
- Decant solution from plate. Wash 4X with 1X wash buffer, 300µL per well. Invert plate and blot dry after last wash.
- Add 100µL 1X Detection Antibody per well.
- Cover plate and incubate 30 minutes at 37°C.
- Repeat decanting/washing step.
- Add 100µL One-Step TMB Substrate per well.
- Cover plate and incubate 5 minutes at RT protected from light.
- Add 50µL stop solution to each well. The color in the wells should change from blue to yellow. Tap the plate to ensure thorough mixing.
- Read OD at 450nm in a microplate reader within 30min of reaction termination.
- Calculate the sample concentration using the standard curve generated by the Calibrators.

Applications Note

This assay kit is for research use only and is NOT intended for use in diagnostic or therapeutic procedures. In native biological samples, there are various interfering substances that affect the accuracy of test results. For more information, visit the 2019-nCoV S1 ELISA Kit page on the SignalChem website.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.