

2019-nCoV N-S1 Human IgA ELISA Kit

Catalog # C19NS1-A877

Product Description

The 2019-nCoV N-S1 Human IgA Kit is an Enzyme-Linked ImmunoSorbent Assay (ELISA) kit for the detection of Human anti-nCoV N & Spike protein S1 IgA in serum or complex biological samples. The kit includes one 96-well strip plate pre-coated with a recombinant 2019-nCoV N & Spike protein RBD Chimera protein. Positive and Negative Controls are included to ensure reliable interpretation of anti-CoV N & S1 Human IgA levels in test samples.

Components – Kit Box

SignalChem Cat#	Component Name	Size
H38A-63DG	Anti-Human IgA Fc Detection Antibody, 1X	10 mL
CP04-C19NSD	2019-nCoV RBD-N Protein Chimera Coated Plate	1 plate
CS02-506	One-step TMB Substrate	10 mL
E02-09	nCoV Ig Sample Dilution Buffer	2 x 10 mL
SS01-09	Stop Solution	6 mL
C19NS1A-NEG C19NS1A-POS	IgA Negative Control IgA Positive Control	1 ml each
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 the 2019-nCoV N-S1 Human IgA ELISA Kit Instruction Manual.

2019-nCoV N-S1 Human IgA 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.
- Add 100µL each of Sample or Control per well (in duplicate or triplicate).
- Cover plate and incubate 1hr at 37°C.
- Decant solution from plate. Wash 4X with 300µL per well of 1X Wash buffer. Invert plate and blot dry after last wash.
- Add 100µL Detection Antibody per well.
- Cover plate and Incubate 1hr at 37°C.
- Repeat decanting/washing step.
- Add 100µL One-Step TMB Substrate per well.
- Cover plate and incubate 10min 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.
- Classify samples as positive or negative based on the O.D readings relative to controls.

Applications Note

This assay kit is for research use only and is NOT intended for use in diagnostic or therapeutic procedures. This kit is limited to the qualitative detection of antibodies specific to the 2019-nCoV virus. A negative or non-reactive result can occur if the quantity of antibodies for the 2019-nCoV virus is below the detection limit. This kit is not validated for use in donor serum screening. For more information, visit the 2019-nCoV N-S1 Human IgA 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.