

Product datasheet

COVID-19 N-Protein Human IgG ELISA Kit ab274339

1 Image

Overview

---

<b>Product name</b>	COVID-19 N-Protein Human IgG ELISA Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Serum
<b>Assay type</b>	Quantitative
<b>Range</b>	1.37 U/ml - 1000 U/ml
<b>Assay duration</b>	Multiple steps standard assay
<b>Product overview</b>	<p>COVID-19 N-Protein Human IgG ELISA Kit (ab274339) is an <i>in vitro</i> indirect ELISA for the quantitative measurement of human IgG antibody against SARS-CoV-2 N protein in human serum.</p>

Standard 96-well plates (12 strips with 8 wells/strip) are coated with the SARS-CoV-2 N protein, which combines with the corresponding antibody present in a sample and Positive Control which used as calibration curve for interpretation purposes. The wells are washed, and biotinylated anti-human IgG antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of COVID19 N protein human IgG antibody bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

The Positive Controls are from an inactivated serum sample which contains SARS-COV-2 N protein human IgG antibody. We do not know the exact amount of SARS-COV-2 N protein human IgG antibody in the Positive Control sample. The Positive Control can be used as calibration curve for interpretation purposes in different assays.

**Notes**

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

**Platform** Pre-coated microplate (12 x 8 well strips)

Properties

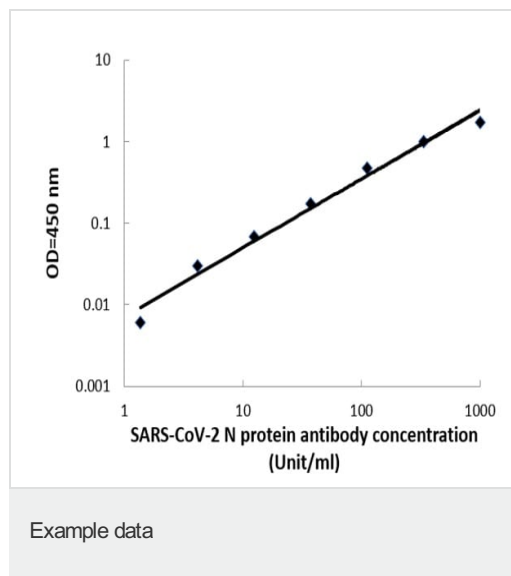
---

## Storage instructions

Store at -80°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
5X Sample Diluent	1 x 25ml
Biotinylated anti-Human IgG Antibody	1 vial
HRP-Streptavidin Concentrate	1 vial
Positive Control	1 vial
SARS-CoV-2 N protein coated Microplate	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

## Images



Example Positive Control calibration curve.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish

- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors