

Product datasheet

COVID-19 S-Protein (S1RBD) Human IgA ELISA Kit  
ab276185

[1 Image](#)

Overview

<b>Product name</b>	COVID-19 S-Protein (S1RBD) Human IgA 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>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	<p>COVID-19 S-Protein (S1RBD) Human IgA ELISA Kit (ab276185) is an <i>in vitro</i> indirect ELISA for the quantitative measurement of human IgA antibody against SARS-CoV-2 S1 RBD protein in human serum.</p> <p>Standard 96-well plates (12 strips with 8 wells/strip) are coated with the SARS-CoV-2 S1 RBD 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 IgA 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 S1 RBD protein human IgA antibody bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p> <p>The Positive Controls are from an inactivated serum sample which contains SARS-COV-2 S1 RBD protein human IgA antibody. We do not know the exact amount of SARS-COV-2 S1 RBD protein human IgA antibody in the Positive Control sample. The Positive Control can be used as a calibration curve for interpretation purposes in different assays.</p>
<b>Notes</b>	<p>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.</p> <p>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.</p>
<b>Tested applications</b>	<b>Suitable for:</b> ELISA
<b>Platform</b>	Pre-coated microplate (12 x 8 well strips)

## Properties

---

### Storage instructions

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
SARS-CoV-2 S1 RBD protein coated Microplate	1 unit
Albumin protein coated 96 well-Microplate	1 unit
20X Wash Buffer	1 x 40ml
Positive Control	2 vials
Biotinylated Anti-Human IgA antibody	2 vials
HRP-Streptavidin Concentrate	1 vial
TMB One-Step Substrate Reagent	1 x 24ml
Stop Solution	1 x 16ml
5X Assay Diluent B	1 x 15ml
5X Sample Diluent	1 x 25ml

## Applications

---

### The Abpromise guarantee

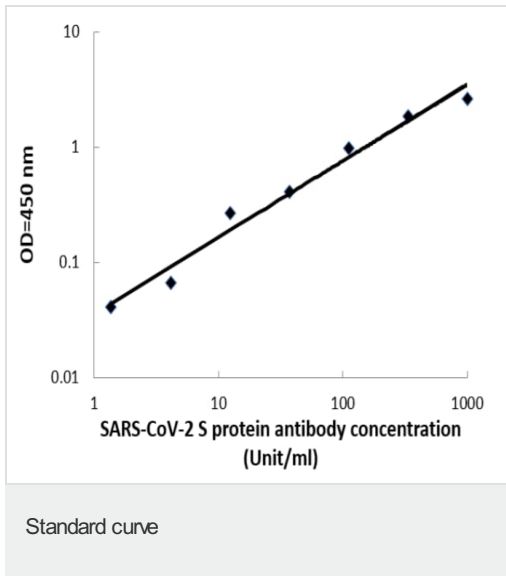
Our [Abpromise guarantee](#) covers the use of ab276185 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

## Images

---



COVID-19 S-Protein (S1RBD) Human IgA ELISA Kit (ab276185)  
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