

Product datasheet

Novel Coronavirus IgM Antibody Detection Kit (SARS-CoV-2) ab272244

1 Image

Overview

Product name	Novel Coronavirus IgM Antibody Detection Kit (SARS-CoV-2)
Sample type	Serum, Plasma, Whole Blood
Species reactivity	Reacts with: Human
Product overview	Novel Coronavirus IgM Antibody Detection Kit (SARS-CoV-2) (ab272244) is suitable for the qualitative detection of Novel Coronavirus (SARS-CoV-2) IgM antibodies in human serum, plasma, and whole blood.

Note: Novel Coronavirus IgG Antibody Detection Kit (SARS-CoV-2) - [ab272243](#) and Novel Coronavirus IgM Antibody Detection Kit (SARS-CoV-2) - [ab272244](#) are intended for Research Use Only, not for use in diagnostic procedures.

The detection kit uses the principle of immunochromatography: the separation of components in a mixture through a medium using capillary force and the specific and rapid binding of an antibody to its antigen. Each cassette is a dry medium that has been coated separately with novel coronavirus N protein (“T” test line) and goat antichickn IgY antibody (“C” control line) (Figure 1). Two free colloidal gold-labeled antibodies, mouse anti-human IgM (mIgM) and chicken IgY, are in the release pad section (S). Once diluted serum, plasma, or whole blood is applied to the release pad section, the mIgM antibody will bind to coronavirus IgM antibodies if they are present, forming an IgM-IgM complex. The sample and antibodies will then move across the cassette’s medium via capillary action. If coronavirus IgM antibody is present in the sample, the test line (T) will be bound by the IgM-IgM complex and develop color. If there is no coronavirus IgM antibody in the sample, free mIgM will not bind to the test line (T) and no color will develop. The free chicken IgY antibody will bind to the control line (C); this control line should be visible after the detection step as this confirms that the kit is working properly.

The following Detection Kits are available:

Novel Coronavirus IgG Antibody Detection Kit (SARS-CoV-2) - [ab272243](#)

Novel Coronavirus IgM Antibody Detection Kit (SARS-CoV-2) - this kit

Notes [ab272243](#) and [ab272244](#) were tested on confirmed positive Covid-19 samples and known negative samples. The following data set was the outcome of two different screening batches

compiled together. Confirmed positive and negative control samples were tested for the presence of IgG and IgM novel coronavirus antibodies. The kit was used as instructed by the manual, and positive or negative indications were listed as such. The results of that testing indicates a sensitivity to viral antibodies of 87%, with a specificity of 84%.

Sample Type	Sample	IgM-Ab272244	IgG - Ab272243	Overall Positive	Overall Negative	Sensitivity	Spe
Positive	30	24	24	26	4	87%	
Negative	90	12	3	14	76		84

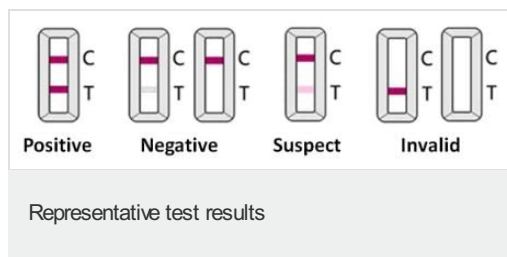
Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	20 tests
Detection Cassette	20 units
Dilution Buffer	20 vials

Relevance Coronavirus specific IgM

Images



Positive for Coronavirus Both the test line (T) and the quality control line (C) are colored dark pink.

Negative for Coronavirus: The test line (T) does not develop color, or a faint gray band may be visible, but the quality control line (C) is colored.

Suspect: A light pink band is an inconclusive result. The sample requires an alternate testing method (such as RT-qPCR) to determine positivity.

Invalid: There is no colored control line (C) band. The results are invalid regardless of whether a red band appears on the test line (T); additional testing is required.

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