

Strongyloides IgG/IgM ELISA kit ab247203

1 Image

Overview

Product name Strongyloides IgG/IgM ELISA kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
sample	24	0.499nM		7.27%
sample	24	1.21nM		4.42%
sample	24	1.106nM		3.89%

Inter-assay

Sample	n	Mean	SD	CV%
sample	12	31.77nM		10.72%
sample	12	24.62nM		9.84%
sample	12	8.68nM		9.69%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Sandwich (qualitative)

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Strongyloides IgG/IgM ELISA Kit (ab247203) is designed for the qualitative determination of IgG/IgM class antibodies against Strongyloides in human serum or plasma (citrate, heparin).

The qualitative immunoenzymatic determination of specific antibodies is based on the ELISA (Enzyme-linked Immunosorbent

Assay) technique. Microplates are coated with specific antigens to bind corresponding antibodies of the sample. After washing the wells to remove all unbound sample material a horseradish peroxidase (HRP) labelled conjugate is added. This conjugate binds to the captured antibodies.

In a second washing step unbound conjugate is removed. The immune complex formed by the bound conjugate is visualized by adding Tetramethylbenzidine (TMB) substrate which gives a blue reaction product. The intensity of this product is proportional to the amount of specific antibodies in the sample. Sulphuric acid is added to stop the reaction. This produces a yellow endpoint colour. Absorbance at 450/620 nm is read using an ELISA microwell plate reader.

Notes

Strongyloides is a genus containing some 50 species of obligate gastrointestinal parasites of vertebrates. *Strongyloides stercoralis* is the scientific name of a human parasitic roundworm causing the disease of strongyloidiasis. The *Strongyloides stercoralis* nematode can parasitize humans. The adult parasitic stage lives in tunnels in the mucosa of the small intestine. *S. stercoralis* can be found in areas with tropical and subtropical climates but cases also occur in temperate area, more frequently in rural areas. *S. stercoralis* has a very low prevalence in societies where fecal contamination of soil or water is rare. Many people infected are usually asymptomatic at first. Symptoms include dermatitis: swelling, itching, larva currens, and mild hemorrhage at the site where the skin has been penetrated. If the parasite reaches the lungs, the chest may feel as if it is burning, and wheezing and coughing may result, along with pneumonia-like symptoms (Löfler's syndrome). The intestines could eventually be invaded, leading to burning pain, tissue damage, sepsis, and ulcers. In severe cases, edema may result in obstruction of the intestinal tract, as well as loss of peristaltic contractions. *Strongyloides* infection in immunocompromised individuals (particularly following the administration of steroids, for example following transplant surgery) can result in disseminated strongyloidiasis, in which worms move beyond the confines of the gut into other organs.

Platform

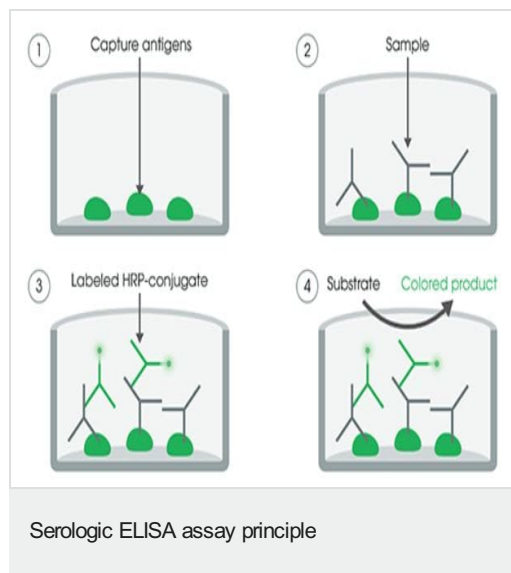
Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions

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

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
anti-human IgG/IgM HRP conjugate		1 x 20ml
Cover Foil		1 unit
Sample Diluent		1 x 100ml
Stop Solution	red cap	1 x 15ml
Strongyloides Coated Microplate		1 unit
Strongyloides Cut-off Control:		1 x 3ml
Strongyloides Negative Control		1 x 2ml
Strongyloides Positive Control		1 x 2ml
TMB Substrate Solution	Yellow cap	1 x 15ml



Specific antigens are coated on the 96-well plate, controls or test samples are added to the well and incubated. The wells are washed to remove any unbound Human anti-antigen antibodies (Ig). A horseradish peroxidase (HRP) labelled anti-Human Ig conjugate is added to the wells. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The intensity of yellow coloration is directly proportional to the amount of Human anti-antigen Ig captured on the plate.

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