

Human Anti-Schistosoma mansoni IgG ELISA Kit ab108769

[1 References](#) [1 Image](#)

Overview

Product name	Human Anti-Schistosoma mansoni IgG ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Pos. Serum	24		5.1%
	Inter-assay			
	Sample	n	Mean	SD
	Pos. Serum	4		5.7%
Sample type	Serum, Hep Plasma, Cit plasma			
Assay type	Indirect			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's anti-Schistosoma mansoni IgG Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Schistosoma mansoni in Human serum and plasma.			

A 96-well plate has been precoated with Schistosoma mansoni antigens to bind cognate antibodies. Controls or test samples are added to the wells and incubated. Following washing, a horseradish peroxidise (HRP) labelled Protein A conjugate is added to the wells, which binds to the immobilized Schistosoma mansoni-specific antibodies. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The density of yellow coloration is directly proportional to the amount of Schistosoma mansoni IgG sample captured in plate.

Platform

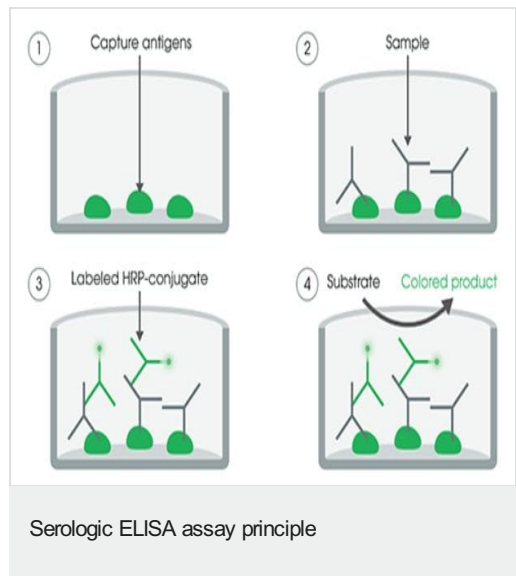
Microplate

Storage instructions

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

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Cover foil		1 unit
IgG Sample Diluent	Yellow, White cap	1 x 100ml
Protein A HRP Conjugate	Blue, Black cap	1 x 20ml
Schistosoma mansoni (IgG) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Schistosoma mansoni IgG Cut-off Control	Yellow, Green cap	1 x 3ml
Schistosoma mansoni IgG Negative Control	Yellow, Blue cap	1 x 2ml
Schistosoma mansoni IgG Positive Control	Yellow, Red cap	1 x 2ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

Images



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