

Human Anti-Candida albicans IgM ELISA Kit ab108716

1 Image

Overview

Product name	Human Anti-Candida albicans IgM ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Pos.Serum	6		CV%
				4.5%
	Inter-assay			
	Sample	n	Mean	SD
	Pos.Serum	38		CV%
				5.1%
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-Candida albicans IgM Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Candida albicans in Human serum or plasma.			

A 96-well plate has been precoated with Candida albicans antigens to bind cognate antibodies. Controls or test samples are added to the wells and incubated. Following washing, a horseradish peroxidase (HRP) labelled anti-Human IgM conjugate is added to the wells, which binds to the immobilized Candida albicans-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 Candida albicans IgM sample captured in plate.

Platform

Microplate

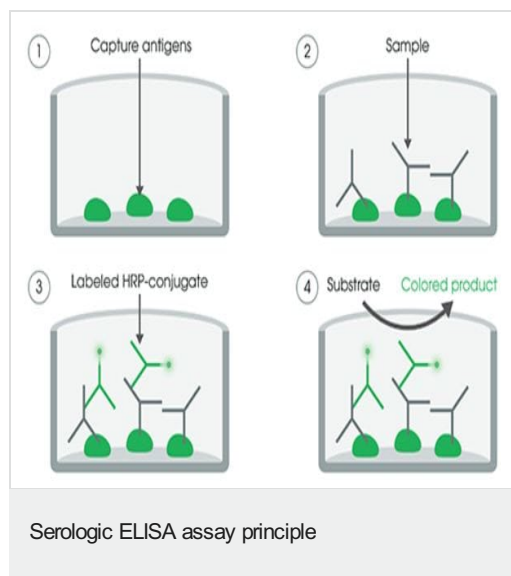
Properties

## Storage instructions

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

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Candida albicans (IgM) Coated Microplate (12 x 8 wells)		1 unit
Candida albicans anti-IgM HRP Conjugate	Red Bottle, Black Cap	1 x 20ml
Candida albicans IgM Cut-off Control	Yellow Bottle, Green Cap	1 x 3ml
Candida albicans IgM Negative Control	Yellow Bottle, Blue Cap	1 x 2ml
Candida albicans IgM Positive Control	Yellow Bottle, Red Cap	1 x 2ml
Foil Cover		1 unit
IgM Sample Diluent	Green Bottle, White Cap	1 x 100ml
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