

Human Anti-Candida albicans IgA ELISA Kit ab108714

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

Overview

Product name	Human Anti-Candida albicans IgA ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Pos.Serum	10			6.6%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Pos.Serum	33			9%
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 IgA Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgA class antibodies against Candida albicans in Human serum and 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 IgA 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 IgA sample captured in plate.				
Platform	Microplate				

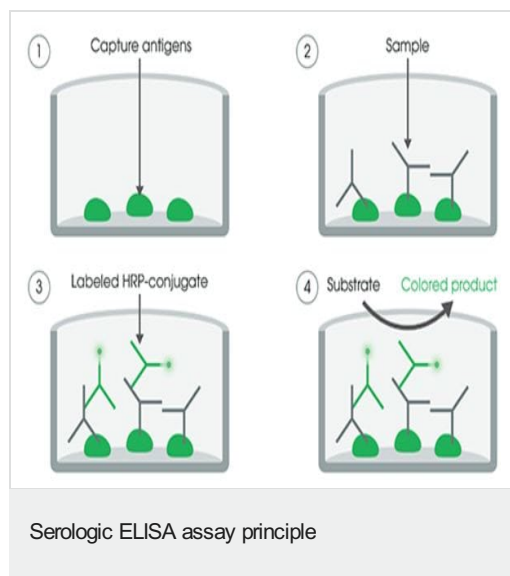
Properties

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 (IgA) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Candida albicans anti-IgA HRP Conjugate	Colored violet; Black cap	1 x 20ml
Candida albicans IgA Cut-off Control	Colored yellow; Green cap	1 x 3ml
Candida albicans IgA Negative Control	Colored yellow; Blue cap	1 x 2ml
Candida albicans IgA Positive Control	Colored yellow; Red cap	1 x 2ml
Cover foil		1 unit
IgA Sample Diluent	Colored yellow; 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