

Human Anti-Varicella-Zoster virus IgA ELISA Kit (VZV) ab108781

1 References 1 Image

Overview

Product name	Human Anti-Varicella-Zoster virus IgA ELISA Kit (VZV)			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Pos. Serum	3		3.8%
	Inter-assay			
	Sample	n	Mean	SD
	Pos. Serum	7		5.3%
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-Varicella-Zoster virus (VZV) IgA Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgA class antibodies against Varicella-Zoster virus (VZV) in Human serum and plasma.			
	A 96-well plate has been precoated with Varicella-Zoster virus (VZV) 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 Varicella-Zoster virus (VZV)-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 Varicella-Zoster virus (VZV) IgA sample captured in plate.			
Platform	Microplate			

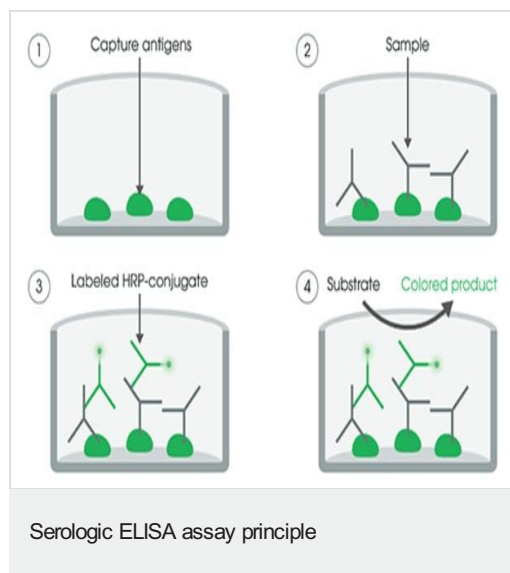
Properties

Storage instructions

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

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
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
Varicella-Zoster virus (IgA) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Varicella-Zoster virus anti-IgA HRP Conjugate	Colored violet; Black cap	1 x 20ml
Varicella-Zoster virus IgA Cut-off Control	Colored yellow; Green cap	1 x 3ml
Varicella-Zoster virus IgA Negative Control	Colored yellow; Blue cap	1 x 2ml
Varicella-Zoster virus IgA Positive Control	Colored yellow; Red cap	1 x 2ml

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