

# Human Anti-Varicella-Zoster virus IgM ELISA Kit (VZV) ab108783

1 [References](#) 1 [图像](#)

概述

产品名称	人Anti-Varicella-Zoster virus IgM ELISA试剂盒(VZV)			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Pos. Serum	8		5.2%
	批次间			
	样品	n	Mean	SD
	Pos. Serum	16		4.2%
样品类型	Serum, Hep Plasma, Cit plasma			
检测类型	Indirect			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Human			
产品概述	Abcam's anti-Varicella-Zoster virus (VZV) IgM Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Varicella-Zoster virus in Human serum and plasma.			

A 96-well plate has been precoated with Varicella-Zoster virus 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 Varicella-Zoster virus-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 IgM sample captured in plate.

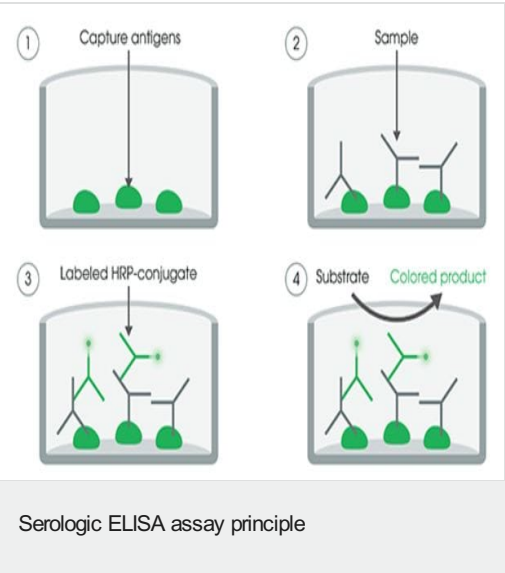
平台 Microplate

存放说明

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

组件	标识符	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Cover foil		1 unit
IgM Sample Diluent	Colored green; 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 (IgM) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Varicella-Zoster virus (VZV) IgM Cut-off Control	Colored yellow; Green cap	1 x 3ml
Varicella-Zoster virus (VZV) IgM Negative Control	Colored yellow; Blue cap	1 x 2ml
Varicella-Zoster virus (VZV) IgM Positive Control	Colored yellow; Red cap	1 x 2ml
Varicella-Zoster virus anti-IgM HRP Conjugate	Colored red; Black cap	1 x 20ml

图片



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.cn/abpromise> or contact our technical team.

#### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors