

# Human Anti-Varicella-Zoster virus IgG ELISA Kit (VZV) ab108782

### 1 References

#### 概述

**产品名称** 人Anti-Varicella-Zoster virus IgG ELISA试剂盒(VZV)

**检测方法** Colorimetric

**精确度** 批次内

样品	n	Mean	SD	CV%
Pos Serum	21			2.4%

批次间

样品	n	Mean	SD	CV%
Pos. Serum	24			4.9%

**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) IgG Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG 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 IgG 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 IgG sample captured in plate.

**平台** Microplate

**性能**

**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
IgG 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 (IgG) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Varicella-Zoster virus anti-IgG HRP Conjugate	Colored blue; Black cap	1 x 20ml
Varicella-Zoster virus IgG Cut-off Control	Colored yellow; Green cap	1 x 3ml
Varicella-Zoster virus IgG Negative Control	Colored yellow; Blue cap	1 x 2ml
Varicella-Zoster virus IgG Positive Control	Colored yellow; Red cap	1 x 2ml

**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