



Product datasheet

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

1 References

Overview

| | | | | |
|--------------------|--|----|------|------|
| Product name | Human Anti-Varicella-Zoster virus IgG ELISA Kit (VZV) | | | |
| Detection method | Colorimetric | | | |
| Precision | Intra-assay | | | |
| | Sample | n | Mean | SD |
| | Pos Serum | 21 | | 2.4% |
| | Inter-assay | | | |
| | Sample | n | Mean | SD |
| | Pos. Serum | 24 | | 4.9% |
| 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-Varicella-Zoster virus (VZV) IgG Human <i>in vitro</i> 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.

Platform

Microplate

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 |
| 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.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