

Product datasheet

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

製品の概要

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

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Pos. Serum	3			3.8%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Pos. Serum	7			5.3%

サンプルの種類 Serum, Heparin Plasma, Citrate Plasma

アッセイタイプ Indirect

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's anti-Varicella-Zoster virus (VZV) IgA Human *in vitro* 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.

アプリケーション 適用あり: Indirect ELISA

試験プラットフォーム Microplate

製品の特性

保存方法

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

内容	ラベル	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		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

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab108781** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Indirect ELISA		Use at an assay dependent concentration.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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