

# Anti-Epstein-Barr Virus IgG Avidity ELISA kit (VCA) ab222940

画像数 1

### 製品の概要

|         |                                                                                                                                                                                                                                                                                                                  |
|---------|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| 製品名     | Anti-Epstein-Barr Virus IgG Avidity ELISA kit (VCA)                                                                                                                                                                                                                                                              |
| 検出方法    | Colorimetric                                                                                                                                                                                                                                                                                                     |
| サンプルの種類 | Serum, Hep Plasma, Cit 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   | <p>The Anti-Epstein-Barr Virus IgG ELISA (Enzyme-Linked Immunosorbent Assay) kit (VCA) (ab222940) is designed for the qualitative determination of Epstein-Barr virus viral capsid (VCA)-specific IgG avidity in human serum or plasma (citrate, heparin) to differentiate between acute and past infection.</p> |

Microplates are coated with specific antigens to bind the corresponding antibodies of the sample (dual pipetting). After washing the wells to remove all unbound sample material, one well is incubated with reagent and the corresponding well with washing buffer. The reagent removes the low-avidity antibodies from the antigens whereas the high-avidity ones are still bound to the specific antigens. After a second washing step to remove the rest of reagent and low-avidity antibodies, a horseradish peroxidase (HRP) labeled conjugate is added. This conjugate binds to the captured antibodies. In a third washing step unbound conjugate is removed. The immune complex formed by the bound conjugate is visualized by adding Tetramethylbenzidine (TMB) substrate which gives a blue reaction product.

The intensity of this product is proportional to the amount of specific antibodies in the specimen. Sulphuric acid is added to stop the reaction. This produces a yellow endpoint color.

Absorbance at 450/620 nm is read using an ELISA microwell plate reader.

### 特記事項

The presence of IgG antibodies to Epstein-Barr Virus indicates the occurrence of the infection but does not distinguish between recent and past infection. Specific IgM antibodies are first detected approximately in ten days and peak at about four weeks post infection. They may persist for several months after acute infections. Based on the evidence that antibody avidity gradually increases after exposure to an immunogen, avidity of IgG antibodies can be used as a marker for distinguishing recent primary from long-term infections. Avidity describes the binding strength of a specific antibody to its antigen. Low-avidity IgG antibodies indicate a primary infection, whereas the presence of IgG antibodies with high avidity points to persistency or reactivation of infection.

**試験プラットフォーム**

Pre-coated microplate (12 x 8 well strips)

**製品の特 性****保存方法**

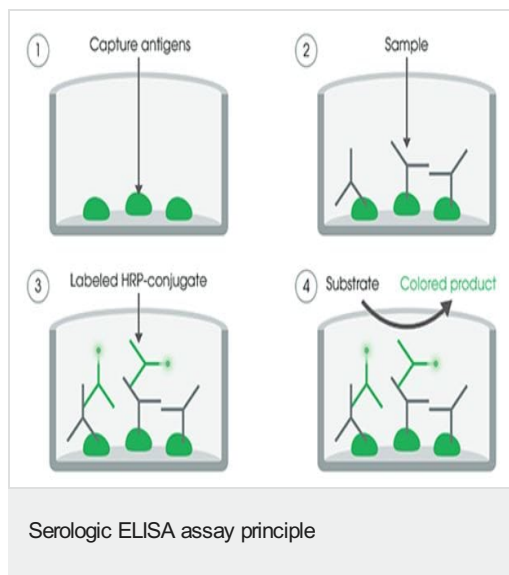
Store at +4°C. Please refer to protocols.

| 内 容                                                       | ラベル                       | 1 x 96 tests |
|-----------------------------------------------------------|---------------------------|--------------|
| 20X Washing Solution                                      | White cap                 | 1 x 50ml     |
| Bottle                                                    |                           | 1 unit       |
| Cover foil                                                |                           | 1 unit       |
| Epstein Barr virus (IgG) Coated Microplate (12 x 8 wells) |                           | 1 unit       |
| Epstein Barr virus anti-IgG HRP Conjugate                 | colored blue; black cap   | 1 x 20ml     |
| Epstein-Barr Virus (VCA) IgG Control High                 |                           | 1 x 2ml      |
| Epstein-Barr Virus (VCA) IgG Control Low                  |                           | 1 x 2ml      |
| Epstein-Barr Virus IgG Cut-off Control                    | colored yellow; green cap | 1 x 3ml      |
| Epstein-Barr Virus IgG Negative Control                   | colored yellow; blue cap  | 1 x 2ml      |
| Epstein-Barr Virus IgG Positive Control                   | colored yellow; red cap   | 1 x 2ml      |
| IgG Sample Diluent                                        | colored yellow; white cap | 1 x 100ml    |
| Reagent                                                   |                           | 1 x 15ml     |
| Stop Solution                                             | red cap                   | 1 x 15ml     |
| Strip holder                                              |                           | 1 unit       |
| TMB Substrate Solution                                    | Yellow cap                | 1 x 15ml     |

**細胞内局在**

Secreted

**画 像**



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.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