

# Human Anti-Herpes simplex virus Type 1 IgG ELISA Kit (HSV1) ab108737

3 References    画像数 1

製品の概要

|            |   |    |     |      |
|------------|---|----|-----|------|
| 製品名        | Human Anti-Herpes simplex virus Type 1 IgG ELISA Kit (HSV1)   |    |     |      |
| 検出方法       | Colorimetric  |    |     |      |
| 再現性        | Intra-Assay (同時再現性)   |    |     |      |
|            | サンプル  | N  | 平均値 | SD   |
|            | Pos.Serum   | 20 |     | CV%  |
|            | Pos.Serum   | 24 |     | 3.4% |
|            |   |    |     | 3.7% |
|            | Inter-Assay (日差再現性)   |    |     |      |
|            | サンプル  | N  | 平均値 | SD   |
|            | Pos.Serum   | 12 |     | CV%  |
|            | Pos.Serum   | 12 |     | 2.7% |
|            |   |    |     | 4.9% |
|            |   |    |     |      |
| 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      | Abcam's anti-Herpes simplex virus Type 1 (HSV1) IgG Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Herpes simplex virus Type 1 in Human serum and plasma.   |    |     |      |
|            | A 96-well plate has been precoated with Herpes simplex virus Type 1 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 Herpes simplex virus Type 1-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 Herpes simplex virus Type 1 IgG sample captured in plate. |    |     |      |
| 試験プラットフォーム | Microplate  |    |     |      |

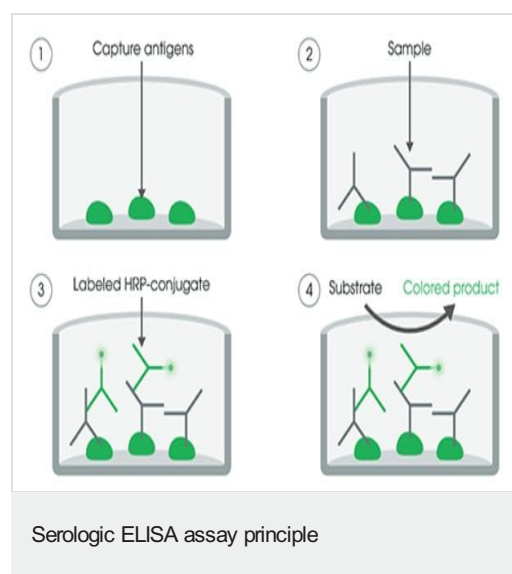
## 製品の特性

### 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       |
| Herpes simplex virus Type 1 (IgG) Coated Microplate (12 x 8 wells) | 12 strips of 8 wells | 1 unit       |
| Herpes simplex virus Type 1 anti-IgG HRP Conjugate                 | Blue, Black cap      | 1 x 20ml     |
| Herpes simplex virus Type 1 IgG Cut-off Control                    | Yellow, Green cap    | 1 x 3ml      |
| Herpes simplex virus Type 1 IgG Negative Control                   | Yellow, Blue cap     | 1 x 2ml      |
| Herpes simplex virus Type 1 IgG Positive Control                   | Yellow, Red cap      | 1 x 2ml      |
| IgG Sample Diluent   | 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     |

## 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