

# Human Anti-Herpes simplex virus Type 2 IgG ELISA Kit (HSV2) ab108739

2 References 1 图像

概述

|       |   |    |      |      |
|-------|---|----|------|------|
| 产品名称  | 人Anti-Herpes simplex virus Type 2 IgG ELISA试剂盒(HSV2)  |    |      |      |
| 检测方法  | Colorimetric  |    |      |      |
| 精确度   | 批次内   |    |      |      |
|       | 样品  | n  | Mean | SD   |
|       | Pos.Serum   | 20 |      | 3.4% |
| 精确度   | 批次内   |    |      |      |
|       | 样品  | n  | Mean | SD   |
|       | Pos.Serum   | 24 |      | 3.7% |
| 精确度   | 批次间   |    |      |      |
|       | 样品  | n  | Mean | SD   |
|       | Pos.Serum   | 12 |      | 3.6% |
| 精确度   | 批次间   |    |      |      |
|       | 样品  | n  | Mean | SD   |
|       | Pos.Serum   | 12 |      | 5.1% |
| 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-Herpes simplex virus Type 2 (HSV2) 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 2 in Human serum and plasma.   |    |      |      |
|       | A 96-well plate has been precoated with Herpes simplex virus Type 2 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 2-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 2 IgG sample captured in plate. |    |      |      |
|       |   |    |      |      |
| 平台    | Microplate  |    |      |      |

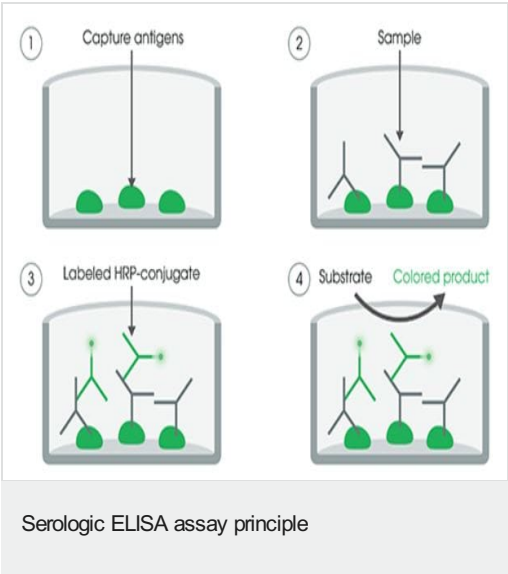
性能

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

图片



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.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