

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

[3 References](#) [1 图像](#)

概述

产品名称 人Anti-Herpes simplex virus Type 1 IgG ELISA试剂盒(HSV1)

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Pos.Serum	20			3.4%
Pos.Serum	24			3.7%

批次间

样品	n	Mean	SD	CV%
Pos.Serum	12			2.7%
Pos.Serum	12			4.9%

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 1 (HSV1) IgG Human *in vitro* 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.

平台 Microplate

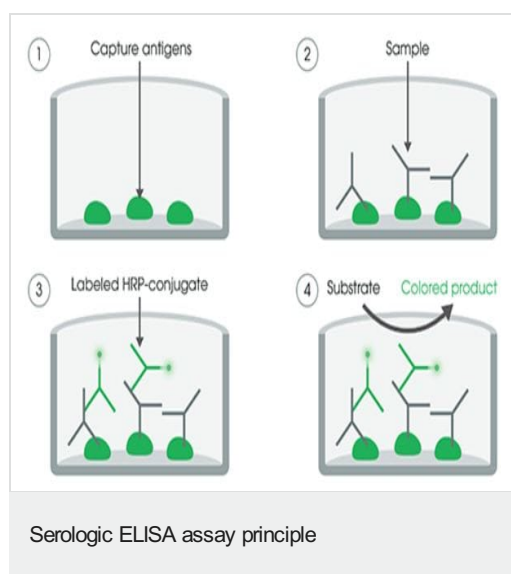
性能

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 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

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