

## Human Anti-Epstein Barr virus IgG ELISA Kit (EBV-EBNA) ab108731

[1 References](#) [1 图像](#)

### 概述

**产品名称** 人Anti-Epstein Barr virus IgG ELISA试剂盒(EBV-EBNA)

**检测方法** Colorimetric

**精确度** 批次内

样品	n	Mean	SD	CV%
Pos. Serum	12			2%

批次间

样品	n	Mean	SD	CV%
Pos. Serum	26			2.4%

**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-Epstein Barr virus (EBV-EBNA) IgG Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG class antibodies against Epstein Barr virus nuclear antigen (EBNA) in Human serum and plasma.

A 96-well plate has been precoated with Epstein Barr virus 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 Epstein Barr virus-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 Epstein Barr virus IgG sample captured in plate.

**平台** Microplate

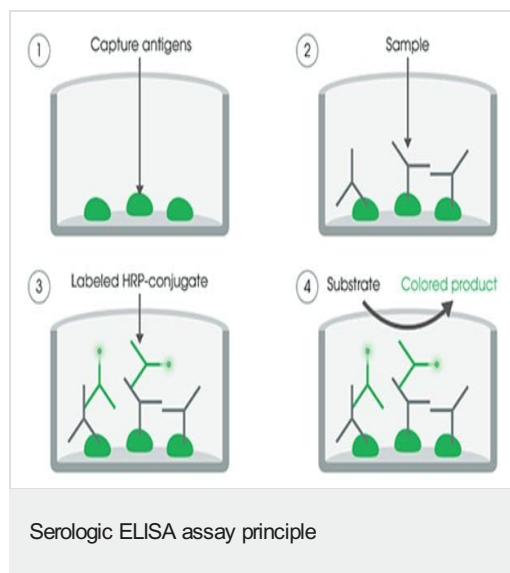
**性能**

## 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
Epstein Barr virus (IgG) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Epstein Barr virus anti-IgG HRP Conjugate	colored blue, black cap	1 x 20ml
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
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