

Human Anti-Measles virus IgM ELISA Kit ab108751

1 图像

概述

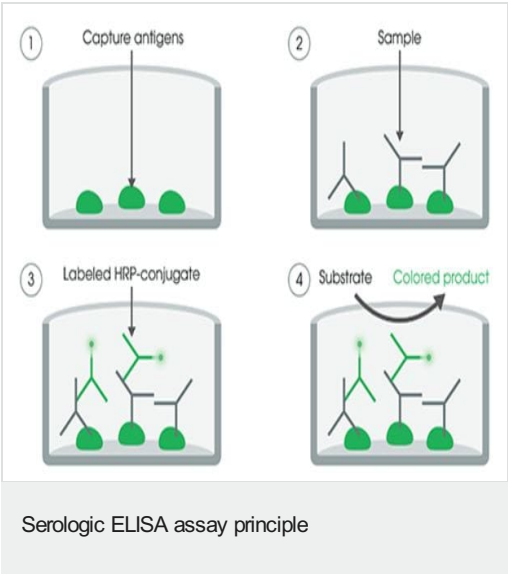
| | | | | |
|-------|---|----|------|------|
| 产品名称 | 人Anti-Measles virus IgM ELISA试剂盒 | | | |
| 检测方法 | Colorimetric | | | |
| 精确度 | 批次内 | | | |
| | 样品 | n | Mean | SD |
| | Pos. Serum | 20 | | 7.7% |
| | 批次间 | | | |
| | 样品 | n | Mean | SD |
| | Pos. Serum | 14 | | 4.2% |
| | 批次间 | | | |
| | 样品 | n | Mean | SD |
| | Pos. Serum | 14 | | 5.6% |
| 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-Measles virus IgM Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Measles virus in Human serum and plasma. | | | |
| | A 96-well plate has been precoated with Measles 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 IgM conjugate is added to the wells, which binds to the immobilized Measles 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 Measles virus IgM 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 |
| IgM Sample Diluent | Green, White cap | 1 x 100ml |
| Measles virus (IgM) Coated Microplate (12 x 8 wells) | 12 strips of 8 wells | 1 unit |
| Measles virus anti-IgM HRP Conjugate | Red, Black cap | 1 x 20ml |
| Measles virus IgM Cut-off Control | Yellow, Green cap | 1 x 3ml |
| Measles virus IgM Negative Control | Yellow, Blue cap | 1 x 2ml |
| Measles virus IgM Positive Control | Yellow, Red cap | 1 x 2ml |
| 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