

Human Anti-Candida albicans IgM ELISA Kit ab108716

1 图像

概述

产品名称 人Anti-Candida albicans IgM ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

| 样品 | n | Mean | SD | CV% |
|-----------|---|------|----|------|
| Pos.Serum | 6 | | | 4.5% |

批次间

| 样品 | n | Mean | SD | CV% |
|-----------|----|------|----|------|
| Pos.Serum | 38 | | | 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-Candida albicans IgM Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Candida albicans in Human serum or plasma.

A 96-well plate has been precoated with Candida albicans 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 Candida albicans-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 Candida albicans IgM sample captured in plate.

平台 Microplate

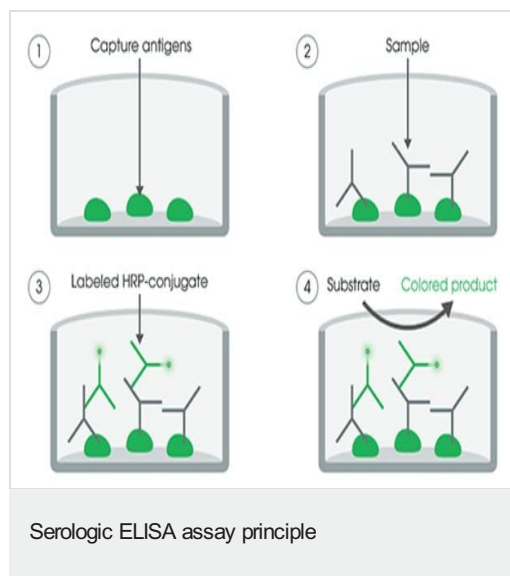
性能

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 |
| Candida albicans (IgM) Coated Microplate (12 x 8 wells) | | 1 unit |
| Candida albicans anti-IgM HRP Conjugate | Red Bottle, Black Cap | 1 x 20ml |
| Candida albicans IgM Cut-off Control | Yellow Bottle, Green Cap | 1 x 3ml |
| Candida albicans IgM Negative Control | Yellow Bottle, Blue Cap | 1 x 2ml |
| Candida albicans IgM Positive Control | Yellow Bottle, Red Cap | 1 x 2ml |
| Foil Cover | | 1 unit |
| IgM Sample Diluent | Green Bottle, 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