

Anti-Malaria ELISA Kit ab178649

1 图像

概述

产品名称

检测方法

精确度

Anti-Malaria ELISA 试剂盒

Colorimetric

批次内

样品	n	Mean	SD	CV%
Positive	22			2.8%
Positive	22			3.9%
Negative	22			5.5%

批次间

样品	n	Mean	SD	CV%
Positive	24			3.2%
Positive	24			4.8%
Negative	24			10.3%

样品类型

检测类型

灵敏度

实验步骤

种属反应性

产品概述

Serum, Hep Plasma, Cit plasma

Indirect

= 95.9 %

Multiple steps standard assay

与反应: Human

Abcam's anti-Malaria Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgG and IgM class antibodies against Malaria in Human serum and plasma.

A 96-well plate has been precoated with Malaria 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 and anti-Human IgM conjugate is added to the wells, which binds to the immobilized Malaria antigens. 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 Malaria IgG or Malaria IgM sample captured in plate.

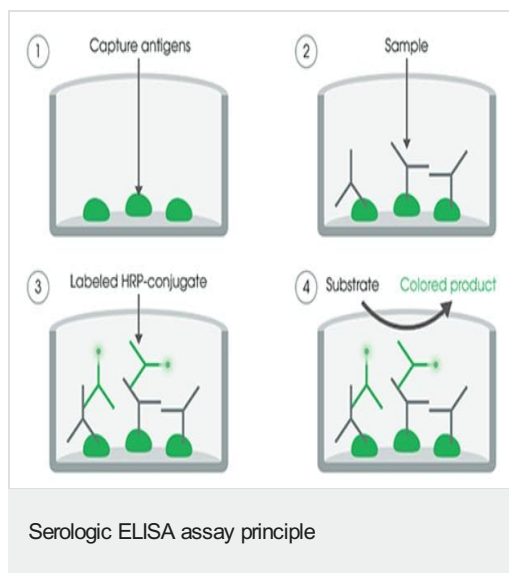
平台 Microplate

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
IgG Sample Diluent		1 x 100ml
Malaria Anti-IgG and Anti-IgM HRP Conjugate		1 x 20ml
Malaria Coated Microplate (12 x 8 wells)		1 unit
Malaria Cut-off Control		1 x 3ml
Malaria Negative Control		1 x 2ml
Malaria Positive Control		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