

Anti-Malaria ELISA Kit ab178649

画像数 1

製品の概要

製品名

Anti-Malaria ELISA Kit

検出方法

Colorimetric

再現性

Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Positive	22			2.8%
Positive	22			3.9%
Negative	22			5.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Positive	24			3.2%
Positive	24			4.8%
Negative	24			10.3%

サンプルの種類

Serum, Hep Plasma, Cit plasma

アッセイタイプ

Indirect

検出感度

= 95.9 %

ステップ

Multiple steps standard assay

種交差性

交差種: Human

製品の概要

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.

試験プラットフォーム

Microplate

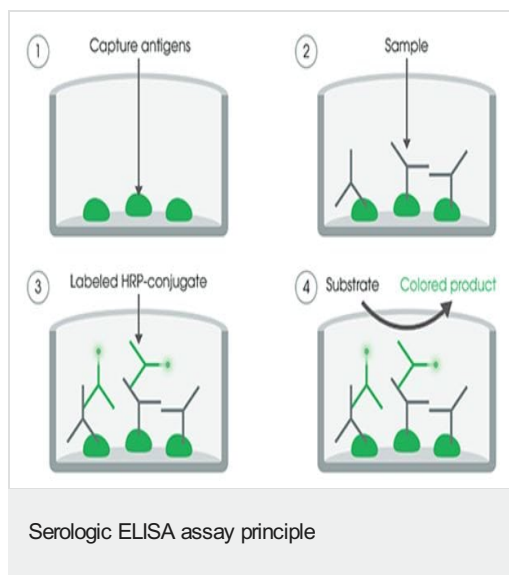
製品の特性

保存方法

Store at +4°C. Please refer to protocols.

内容	ラベル	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

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors