

Human Anti-Measles virus IgM ELISA Kit ab108751

画像数 1

製品の概要

製品名 Human Anti-Measles virus IgM 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%
Pos. Serum	20			7.7%
Pos. Serum	24			6.9%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Pos. Serum	14			4.2%
Pos. Serum	14			5.6%

サンプルの種類 Serum, Hep Plasma, Cit plasma

アッセイタイプ Indirect

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's anti-Measles virus IgM Human *in vitro* 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.

試験プラットフォーム Microplate

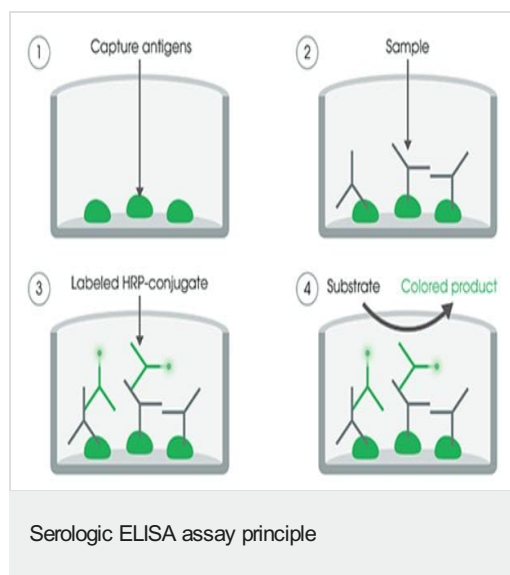
製品の特性

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
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

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