

Product datasheet

Human Anti-Tick-borne encephalitis virus IgM ELISA Kit (TBE) ab108774

製品の概要

製品名 Human Anti-Tick-borne encephalitis virus IgM ELISA Kit (TBE)

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%
TBE virus	12			7.5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
TBE virus	24			8.9%

サンプルの種類 Serum, 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-Tick-borne encephalitis (TBE) virus IgM Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Tick-borne encephalitis virus in Human serum and plasma.

A 96-well plate has been precoated with Tick-borne encephalitis 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 Tick-borne encephalitis-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 Tick-borne encephalitis IgM sample captured in plate.

アプリケーション 適用あり: Indirect ELISA

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
Stop Solution		1 x 15ml
Strip holder		1 unit
Tick-borne encephalitis (IgM) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
Tick-borne encephalitis virus anti-IgM HRP Conjugate	Red, Black cap	1 x 20ml
Tick-borne encephalitis virus IgM Cut-off Control	Yellow, Green cap	1 x 3ml
Tick-borne encephalitis virus IgM Negative Control	Yellow, Blue cap	1 x 2ml
Tick-borne encephalitis virus IgM Positive Control	Yellow, Red cap	1 x 2ml
TMB Substrate Solution	Yellow cap	1 x 15ml

アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab108774** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

アプリケーション	Abreviews	特記事項
Indirect ELISA		Use at an assay dependent dilution.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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