

# Human Tick-Borne Encephalitis Virus plus IgG ELISA Kit (TBE/FSME) ab108763

[1 Image](#)

### Overview

**Product name** Human Tick-Borne Encephalitis Virus plus IgG ELISA Kit (TBE/FSME)

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Standard B	7	0.32U/ml		6.8%
Standard C	8	0.98U/ml		5.7%
Standard D	8	1.3U/ml		4.7%
Standard E	8	2.15U/ml		4.2%

Inter-assay

Sample	n	Mean	SD	CV%
Standard B	14	0.31U/ml		7.35%
Standard C	16	0.92U/ml		8.3%
Standard D	15	1.37U/ml		7.8%
Standard E	16	2.28U/ml		6.7%

**Sample type** Serum, Hep Plasma, Cit plasma

**Assay type** Indirect

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's anti-Tick-borne encephalitis (TBE) virus IgG plus Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of IgG 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 IgG 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 IgG sample captured in plate.

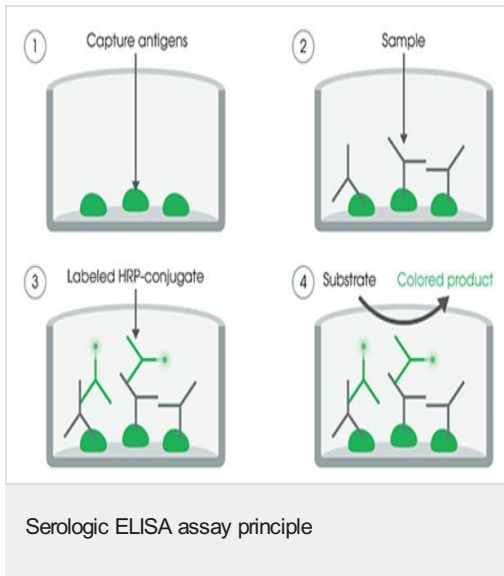
**Platform** Microplate

## Properties

**Storage instructions** Store at +4°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
20X Washing Solution	White cap	1 x 50ml
Cover foil		1 unit
IgG Sample Diluent	Yellow, White cap	1 x 100ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TBE virus (IgG) Coated Microplate (12 x 8 wells)	12 strips of 8 wells	1 unit
TBE virus anti-IgG HRP Conjugate	Blue, Black cap	1 x 20ml
TBE virus IgG Additional Control	Yellow, Orange cap	1 x 2ml
TBE virus IgG Standard A – 0 U/mL	Yellow, Blue cap	1 x 2ml
TBE virus IgG Standard B – 50 U/mL	Yellow, Green cap	1 x 2ml
TBE virus IgG Standard C – 130 U/mL	Yellow, Yellow cap	1 x 2ml
TBE virus IgG Standard D – 200 U/mL	Yellow, Red cap	1 x 2ml
TBE virus IgG Standard E – 300 U/mL	Yellow, White cap	1 x 2ml
TMB Substrate Solution	Yellow cap	1 x 15ml

## Images



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.com/abpromise> or contact our technical team.

### Terms and conditions

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