

Human Anti-Brucella IgM ELISA Kit ab108713

1 Image

Overview

Product name Human Anti-Brucella IgM ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Pos.Serum	24			5.7%
	Pos.Serum	24			3.9%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Pos.Serum	12			3.2%
	Pos.Serum	12			2.2%

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

A 96-well plate has been precoated with Brucella 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 Brucella-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 Brucella IgM sample captured in plate.

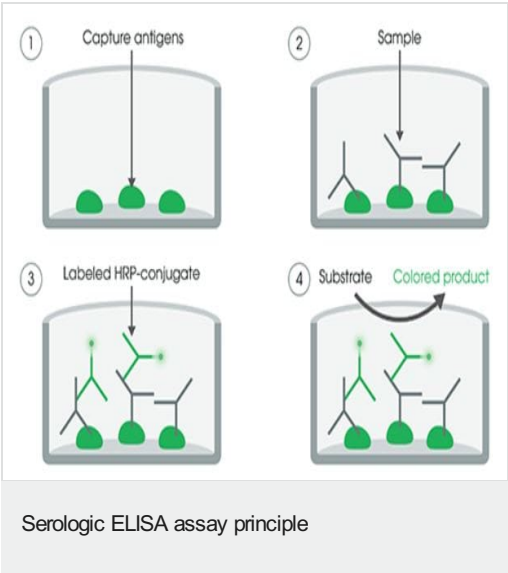
Platform Microplate

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
Brucella (IgM) Coated Microplate (12 x 8 wells)		1 unit
Brucella anti-IgM HRP Conjugate		1 x 20ml
Brucella IgM Cut-off Control		1 x 3ml
Brucella IgM Negative Control		1 x 2ml
Brucella IgM Positive Control		1 x 2ml
Cover foil		1 unit
IgM Sample Diluent		1 x 100ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
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