



Product datasheet

# Human Anti-Chikungunya Virus IgM ELISA Kit ab177848

11 References

Overview

Product name	Human Anti-Chikungunya Virus IgM ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Low + serum	23			4.8%
	High + serum	24			3.4%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Low + serum	3			7.3%
	High + serum	3			2.9%
Sample type	Serum, Hep Plasma, Cit plasma				
Assay type	Sandwich (qualitative)				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	<p>Abcam's anti-Chikungunya Virus IgM Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of IgM class antibodies against Chikungunya Virus in Human serum and plasma.</p> <p>A 96-well plate has been precoated with anti-Human antibodies to bind to corresponding antibodies of the sample. Controls or test samples are added to the wells and incubated. Following washing, the Chikungunya Virus antigen is added to the wells and incubated. The wells are then washed as before and the biotinylated Chikungunya Virus antibodies are added to each well followed by incubation. After a last wash step, the streptavidin peroxidase (SP) conjugate which binds to the biotinylated Chikungunya Virus-specific antibodies is added to the wells. TMB is then catalyzed by the SP 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</p>				

Chikungunya Virus 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
Anti-Human (IgM) Coated Microplate (12 x 8 wells)		1 unit
Chikungunya IgM Negative Control		1 x 1.5ml
Chikungunya Virus antigen Solution 1 (Lyophilized)		6 vials
Chikungunya Virus biotinylated antibody Solution 2		1 x 6ml
Chikungunya Virus IgM Cut-off Control		1 x 2ml
Chikungunya Virus IgM Positive Control		1 x 1.5ml
Cover Foil		1 unit
Sample Diluent		1 x 100ml
Stop Solution	red cap	1 x 15ml
Streptavidin Conjugate		1 x 6ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

**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