

Product datasheet

Anti-Rubella Virus IgM ELISA Kit ab178654

Overview

Product name Anti-Rubella Virus IgM ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Pos Serum	23			2.7%
Pos Serum	24			4%

Inter-assay

Sample	n	Mean	SD	CV%
Pos. Serum	3			10.3%
Pos. Serum	3			11.5%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Sandwich (qualitative)

Assay duration Multiple steps standard assay

Product overview

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

A 96-well plate has been precoated with anti-Human IgM antibodies to bind to corresponding antibodies of the sample. Controls or test samples are added to the wells and incubated. Following washing, a horseradish peroxidase (HRP) labelled Rubella Virus antigen conjugate is added to the wells, which binds to the captured Rubella Virus-specific IgM 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 Rubella 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
Cover Foil		1 unit
IgM Sample Diluent		1 x 100ml
Rubella Virus HRP Conjugate		1 x 12ml
Rubella Virus IgM Cut-off Control		1 x 3ml
Rubella Virus IgM Negative Control		1 x 2ml
Rubella Virus IgM Positive Control		1 x 2ml
Stop Solution	red cap	1 x 15ml
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