

## Product datasheet

# Anti-Rheumatoid Factor IgM ELISA Kit ab178653

★★★★★ 1 Abreviews 2 References 1 Image

### Overview

<b>Product name</b>	Anti-Rheumatoid Factor IgM ELISA Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Serum, Hep Plasma, Cit plasma
<b>Assay type</b>	Indirect
<b>Sensitivity</b>	< 0.5 U/ml
<b>Assay duration</b>	Multiple steps standard assay
<b>Product overview</b>	<p>Abcam's anti-Rheumatoid Factor IgM ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative determination of autoantibodies to Rheumatoid Factor IgM in Human serum or plasma.</p> <p>The quantitative immunoenzymatic determination of IgM rheumatoid factors is based on the ELISA (Enzyme-linked Immunosorbent Assay) technique. Microtiter strips are precoated with Fc fragments of Human immunoglobulin G to bind corresponding antibodies of the specimen. After washing the wells to remove all unbound sample material horseradish peroxidase (HRP) labelled anti-Human IgM conjugate is added. This conjugate binds to the captured rheumatoid factors. The immune complex formed by the bound conjugate is visualized by adding Tetramethylbenzidine (TMB) substrate which gives a blue reaction product. The intensity of this product is proportional to the amount of IgM rheumatoid factors in the specimen. Sulphuric acid is added to stop the reaction. This produces a yellow endpoint colour. Absorbance at 450 nm is read using an ELISA microwell plate reader.</p>
<b>Platform</b>	Microplate

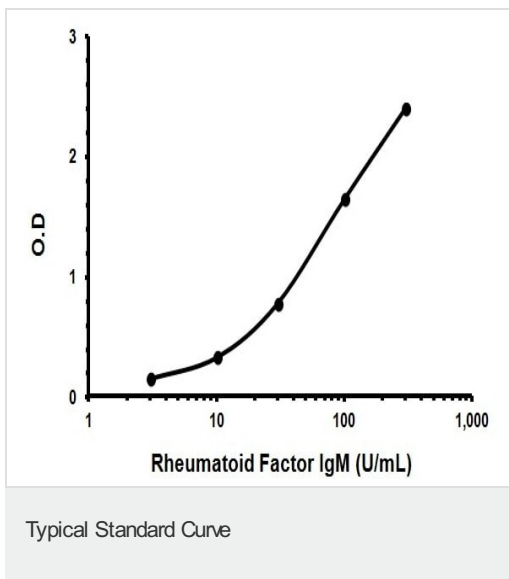
### Properties

**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 Fc Fragment Coated Microplate (12 x 8 wells)		1 unit

Components	Identifier	1 x 96 tests
IgG Sample Diluent		1 x 100ml
Rheumatoid Factor Anti-IgM HRP Conjugate		1 x 20ml
Rheumatoid Factor IgM Negative Control		1 x 2ml
Rheumatoid Factor IgM Positive Control		1 x 2ml
Rheumatoid Factor IgM Standard A - 0 U/mL		1 x 2ml
Rheumatoid Factor IgM Standard B - 3 U/mL		1 x 2ml
Rheumatoid Factor IgM Standard C - 10 U/mL		1 x 2ml
Rheumatoid Factor IgM Standard D - 30 U/mL		1 x 2ml
Rheumatoid Factor IgM Standard E - 100 U/mL		1 x 2ml
Rheumatoid Factor IgM Standard F - 300 U/mL		1 x 2ml
Stop Solution	red cap	1 x 15ml
Strip holder		1 unit
TMB Substrate Solution	Yellow cap	1 x 15ml

**Images**



Representative Standard Curve using ab178653.

**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