



Note that this data sheet is not lot-specific. Please consult the vial label and the certificate of analysis for information on specific lots.

Human RBM3 ELISA Kit

Catalogue Number: BTRBM-001

Package Size: 1 kit (96 wells)

1. Brief Description

The Human RBM3 ELISA provides a fast, highly sensitive and specific quantitative of RNA-binding protein 3 (RBM3) in human serum or plasma samples. Do not use for any other biological sample. One kit contains reagents for 96 determinations, thus allowing the measurement of one standard curve and 40 samples in duplicate.

The calibration curve covers the range from 31.25 pg/ml to 2000 pg/ml. The sensitivity of the assay is 10 pg/ml. The dilution of the serum or plasma samples is recommended 1:10.

2. Kit components

1 Microplate (12 strips with 8 wells) coated with Anti-RBM3-Antibody, 1 Sample dilution buffer, 1 RBM3 Standard concentrate, 1 RBM3 dilution buffer (for preparing RBM3 standards: 0, 31.25, 62.5, 125, 250, 500, 1000 and 2000 pg/ml), 1 RBM3 conjugate concentrate, 1 RBM3 conjugate dilution buffer
Washing buffer concentrate (10x), TMB reagent, Stopping solution (1M H₂SO₄)

3. Other materials and equipment required

Micropipettes, microplate shaker, microplate reader (450 nm / 620 nm), distilled water

4. Sample Preparation

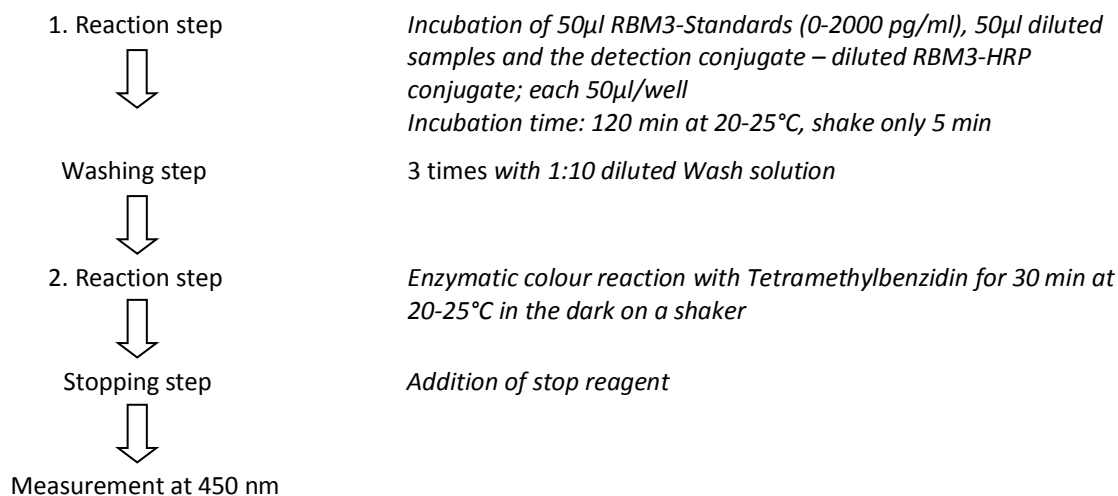
Serum or plasma samples (EDTA or Heparin as anticoagulant) may be stored at -20°C or -80°C. When stored at -20°C or -80°C, it is absolutely necessary to mix the samples thoroughly prior to measuring. Avoid freeze-thaw cycles. Dilute the serum or plasma samples minimum 1:10 with RBM3-SAMPLE-BUF, depending on the possible concentration of the analyte.

5. Principle of the assay

The Human RBM3 ELISA (Enzyme-Linked Immunosorbent Assay) is an immunological quantitative detection method based on a sandwich ELISA.

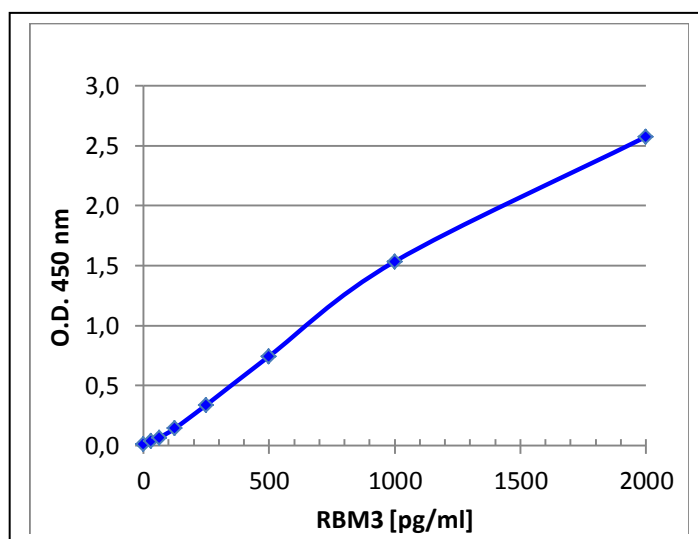
Human RBM3 ELISA is performed in a 96-well microplate. The wells of the microplate are pre-coated with a specific capture antibody against human RBM3 that binds free RBM3 from the patient sample. In one step simultaneous incubation of the RBM3 standards or samples and of the detection conjugate (peroxidase conjugate of an anti-RBM3 antibody) is carried out. Incubation takes place 2 hours at 20-25°C. After a washing step, the colour substrate TMB (tetramethylbenzidine) is added to the wells. The reaction is stopped by adding the Stop Solution and the yellow colour is read in a microtiter plate reader at 450 nm. The concentration of RBM3 in a sample is determined by interpolation from the standard curve. The RBM3 concentration is directly proportional to the colour intensity.

6. Reaction scheme



7. Example of standard value and standard curve

RBM3 Standard (pg/ml)	RBM3 Standard curve Absorbance (450 nm)		
	Value 1	Value 2	Mean Value
0.00	0.0130	0.0110	0.0120
31.25	0.0350	0.0370	0.0360
62.50	0.0650	0.0710	0.0680
125.00	0.1420	0.1480	0.1450
250.00	0.3410	0.3470	0.3440
500.00	0.7250	0.7670	0.7460
1000.00	1.5640	1.5240	1.5440
2000.00	2.5610	2.5630	2.5620



Note: These values are only an example.
 The standard value of RBM3-Standards
 has to be measured in each test.

8. Test evaluation

The evaluation can be performed using commercial ELISA programs. A standard curve is generated by plotting the mean absorbance (x-axis, fig. 2) against pg/ml standard (y-axis, fig.2).
 The pg/ml values of the samples can be read directly from the graph or calculated by the regression coefficients. The recalculated RBM3 values must be further converted by the appropriate sample dilution factor.

9. Assay Parameter

- Target: Human RNA Binding Protein 3 (RBM3)
- Test format: microplate (12 strips with 8 wells)
- RBM3 standard range: 0, 31.25, 62.5, 125, 250, 500, 1000, 2000 pg/ml RBM3
- Sample: serum and plasma
- Sample Dilution: 1:10 is recommended
For dilutions of 1:20 to 1:30, the linearity of the assay results to their dilution is assured.
- Incubation time: 1 x 120 min (sample, standard, conjugate) +1 x 30 min TMB
- Detection limit: <10 pg/ml RBM3
- Recovery rate: 82 – 110 %
- Cross reactivity: No significant cross reactivity to Cold inducible Binding protein (CIRBP), human
- Intra assay variation: <5%
- Inter assay variation: <5%
- Storage: 2-8°C
- Shelf Life: 12 month under storage conditions
- Note: For research use only.

Catalogue Number: BTRBM-001

Gefördert durch:



Bundesministerium
für Wirtschaft
und Technologie

aufgrund eines Beschlusses
des Deutschen Bundestages

distributed in the US/Canada by:
EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063
Phone: 617-419-2019 FAX: 617-419-1110
www.EagleBio.com • info@eaglebio.com



EAGLE
BIOSCIENCES