

Mouse Retinol binding protein 4 ELISA Kit (RBP4)
ab195459

1 Image

Overview

Product name	Mouse Retinol binding protein 4 ELISA Kit (RBP4)				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall				3.2%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall				9.2%
Sample type	Cell culture supernatant, Urine, Serum, Plasma, Tissue, Cell Lysate				
Assay type	Sandwich (quantitative)				
Sensitivity	0.44 ng/ml				
Range	1.25 ng/ml - 20 ng/ml				
Recovery	97 %				
Assay time	4h 00m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Mouse				
Product overview	Abcam's Retinol binding protein 4 (RBP4) Mouse <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of RBP4 concentrations in mouse plasma, serum, urine, cell lysate, and cell culture supernatants.				

An RBP4 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an RBP4 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop

solution. The density of yellow coloration is directly proportional to the amount of RBP4 captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

Platform Microplate

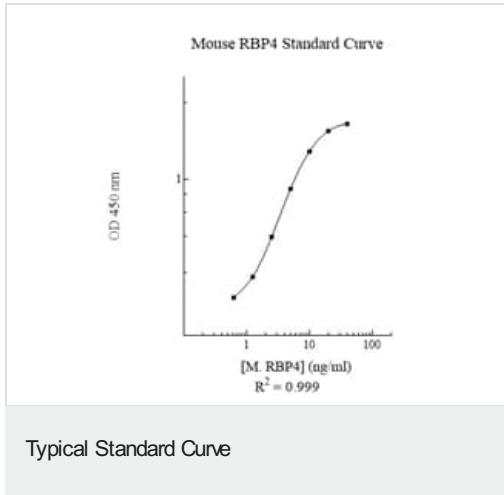
Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
20X Wash Buffer Concentrate	2 x 30ml
25X Biotinylated Mouse RBP4 Antibody	1 x 200µl
Chromogen Substrate	1 x 7ml
Mouse RBP4 Microplate (12 x 8 well strips)	1 unit
Mouse RBP4 Standard (Lyophilised)	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

Function	Delivers retinol from the liver stores to the peripheral tissues. In plasma, the RBP-retinol complex interacts with transthyretin, this prevents its loss by filtration through the kidney glomeruli.
Involvement in disease	Defects in RBP4 are a cause of retinol-binding protein deficiency (RBP deficiency) [MIM:180250]. This condition causes night vision problems. It produces a typical 'fundus xerophthalmicus', featuring a progressed atrophy of the retinal pigment epithelium.
Sequence similarities	Belongs to the calycin superfamily. Lipocalin family.
Cellular localization	Secreted.

Images



Representative standard curve using ab195459.

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