

Mouse Retinol Binding Protein ELISA Kit (RBP) ab190535

1 Image

Overview

Product name Mouse Retinol Binding Protein ELISA Kit (RBP)

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Sample type Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.152 ng/ml

Range 0.625 ng/ml - 20 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	> 85	% - %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview

Abcam's Retinol Binding Protein (RBP) Mouse ELISA kit (ab190535) is an *in vitro* enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of Retinol Binding Protein (RBP) in mouse serum and plasma.

In this assay the Retinol Binding Protein (RBP) present in samples reacts with the anti-Retinol Binding Protein (RBP) antibodies which have been adsorbed to the surface of polystyrene microtitre wells. After the removal of unbound proteins by washing, anti-RBP antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound RBP. Following another washing step, the enzyme bound to the immunosorbent is assayed by the addition of a chromogenic substrate, 3,3',5,5'-

tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of RBP in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of RBP in the test sample. The quantity of RBP in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

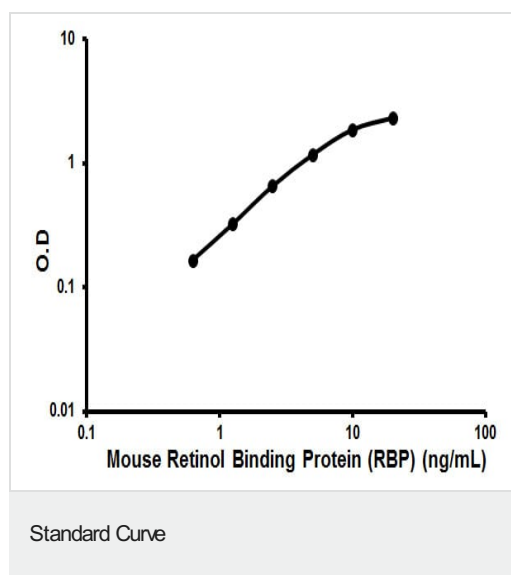
Platform Microplate

Properties

Storage instructions Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 50ml
5X Diluent Concentrate	1 x 50ml
RBP Mouse Antibody Coated Microwells	1 unit
RBP Mouse Calibrator (Lyophilised)	1 vial
RBP Mouse HRP Conjugate	1 vial
Stop Solution	1 x 12ml
Chromogen Substrate Solution	1 x 12ml

Images



Representative standard curve using ab190535 Retinol Binding Protein (RBP) Mouse ELISA Kit.

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