

Product datasheet

# Mouse Retinol Binding Protein 4 Matched Antibody Pair Kit ab214479

1 Image

Overview

---

<b>Product name</b>	Mouse Retinol Binding Protein 4 Matched Antibody Pair Kit
<b>Detection method</b>	Colorimetric
<b>Assay type</b>	ELISA set
<b>Range</b>	15.62 pg/ml - 1000 pg/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Mouse
<b>Product overview</b>	<p>Matched Antibody Pair kit is a titrated unlabeled capture antibody, a titrated biotin-labeled detector and a calibrated protein standard. The Matched Antibody Pair Kit can be used to quantify native and recombinant mouse Retinol Binding Protein 4. Both capture and detector antibodies are rabbit monoclonal antibodies.</p> <p>Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.</p> <p>Additional protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our <a href="#">website</a>.</p> <p>For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA® (<a href="#">ab202404</a>) which uses the same antibody pair.</p> <p><b>To receive an electronic copy of the Certificate of Analysis, please send an <a href="#">email</a> with "CoA for matched antibody pair kit" in the subject line and the desired product number and lot number in the body of the email.</b></p> <p>Buffer information:            The capture antibody is glycerol free.            The detector antibody contains glycerol.</p>
<b>Tested applications</b>	<b>Suitable for:</b> ELISA
<b>Platform</b>	Reagents

## Properties

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

Components	10 x 96 tests	2 x 96 tests
Mouse Retinol Binding Protein 4 Capture Antibody	1 x 100µg	1 x 20µg
Mouse Retinol Binding Protein 4 Detector Antibody	1 x 25µg	1 x 5µg
Mouse Retinol Binding Protein 4 Lyophilized Protein	1 vial	1 vial

**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.

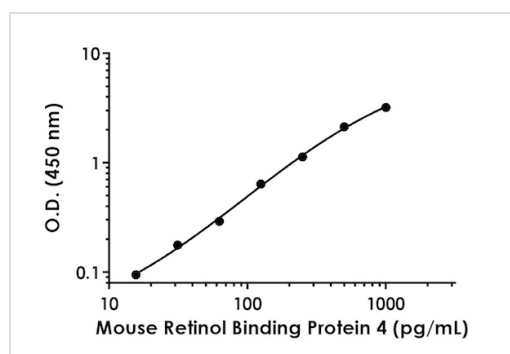
## Applications

Our [Abpromise guarantee](#) covers the use of **ab214479** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
ELISA		Use at an assay dependent concentration.

## Images



Mouse Retinol Binding Protein 4 standard curve.

Standard calibration curve. Background subtracted values are graphed.

**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