

Human Retinol binding protein 4 ELISA Kit (RBP4) ab108897

5 References 画像数 1

製品の概要

製品名	Human Retinol binding protein 4 ELISA Kit (RBP4)			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				3.8%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				8.5%
サンプルの種類	Cell culture supernatant, Saliva, Milk, Urine, Serum, Plasma, Tissue, Cell Lysate, Cerebral Spinal Fluid			
アッセイタイプ	Sandwich (quantitative)			
検出感度	= 8.3 ng/ml			
検出範囲	25 ng/ml - 200 ng/ml			
添加回収試験	97 %			
全工程の試験時間	4h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	Abcam's Retinol binding protein 4 (RBP4) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Retinol binding protein 4 concentrations in plasma, serum, saliva, urine, milk, cerebrospinal fluid and cell culture supernatants.			

A Retinol binding protein 4 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a Retinol binding protein 4 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 Retinol binding protein 4 captured in plate.

Get higher sensitivity in only 90 minutes with Human Retinol Binding Protein 4 ELISA Kit ([ab196264](#)) from our SimpleStep ELISA® range.

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

#### 試験プラットフォーム

Microplate

#### 製品の特性

#### 保存方法

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

内容	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
40X Biotinylated Human Retinol binding Protein 4 Antibody	1 x 150µl
Chromogen Substrate	1 x 7ml
Retinol binding Protein 4 Microplate (12 x 8 well strips)	1 unit
Retinol binding Protein 4 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

#### 機能

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.

#### 関連疾患

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.

#### 配列類似性

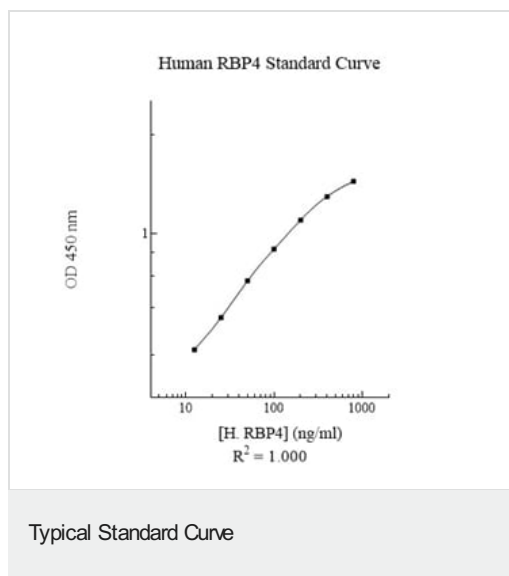
Belongs to the calycin superfamily. Lipocalin family.

#### 細胞内局在

Secreted.

#### 画像





Representative Standard Curve Using ab108897.

**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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors