

Human Retinol binding protein ELISA Kit (RBP) ab108899

★★★★★ 1 Abreviews 2 图像

概述

|       |   |   |      |      |
|-------|---|---|------|------|
| 产品名称  | 人Retinol binding蛋白ELISA试剂盒(RBP)   |   |      |      |
| 检测方法  | Colorimetric  |   |      |      |
| 精确度   | 批次内   |   |      |      |
|       | 样品  | n | Mean | SD   |
|       | Overall   |   |      | 4.6% |
|       | 批次间   |   |      |      |
|       | 样品  | n | Mean | SD   |
|       | Overall   |   |      | 7.2% |
| 样品类型  | Cell culture supernatant, Saliva, Milk, Urine   |   |      |      |
| 检测类型  | Sandwich (quantitative)   |   |      |      |
| 灵敏度   | 0.0078 µg/ml  |   |      |      |
| 范围    | 0.008 µg/ml - 0.5 µg/ml   |   |      |      |
| 回收率   | 98 %  |   |      |      |
| 检测时间  | 4h 00m  |   |      |      |
| 实验步骤  | Multiple steps standard assay   |   |      |      |
| 种属反应性 | 与反应: Human  |   |      |      |
| 产品概述  | Abcam's Retinol binding protein (RBP) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Retinol binding protein concentrations in Human milk, urine, CSF, saliva and cell culture supernatants. |   |      |      |

A Retinol binding protein 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 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 captured in plate.

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

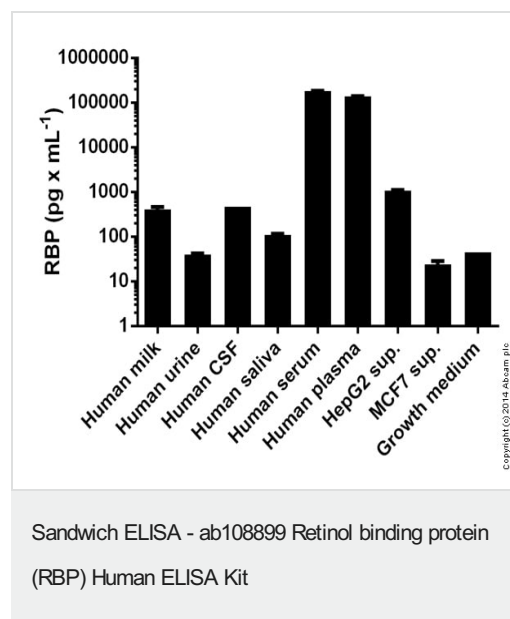
平台 Microplate

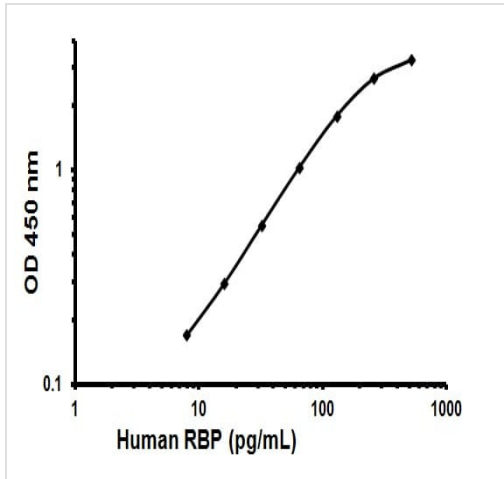
## 性能

存放说明 Store at -20°C. Please refer to protocols.

| 组件  | 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 RBP Antibody                     | 1 x 150µl    |
| Chromogen Substrate                                     | 1 x 8ml      |
| Retinol binding protein Microplate (12 x 8 well strips) | 1 unit       |
| Retinol binding protein Standard                        | 1 vial       |
| Sealing Tapes   | 3 units      |
| Stop Solution   | 1 x 12ml     |

## 图片





Representative Standard Curve using ab108899.

Typical Standard Curve

**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.cn/abpromise> or contact our technical team.

### Terms and conditions

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