

Human Prolactin ELISA Kit ab108679

1 图像

| 概述   |   |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
|--|---|-------|-------|-------|-------|---------|---------|------|------|-------|---------|---------|-------|-------|-------|---------|---------|-------|-------|-------|-------|
| 产品名称   | 人Prolactin ELISA试剂盒   |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 检测方法   | Colorimetric  |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 精确度  | 批次内   |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
|  | <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Level 1</td><td>20</td><td>6.32</td><td>0.4</td><td>6.3%</td></tr><tr><td>Level 2</td><td>20</td><td>19.22</td><td>1.19</td><td>6.18%</td></tr><tr><td>Level 3</td><td>20</td><td>54.52</td><td>2.64</td><td>4.84%</td></tr></table> | 样品    | n     | Mean  | SD    | CV%     | Level 1 | 20   | 6.32 | 0.4   | 6.3%    | Level 2 | 20    | 19.22 | 1.19  | 6.18%   | Level 3 | 20    | 54.52 | 2.64  | 4.84% |
|  | 样品  | n     | Mean  | SD    | CV%   |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
|  | Level 1   | 20    | 6.32  | 0.4   | 6.3%  |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
|  | Level 2   | 20    | 19.22 | 1.19  | 6.18% |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
|  | Level 3   | 20    | 54.52 | 2.64  | 4.84% |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
|  | 批次间   |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Level 1</td><td>20</td><td>6.58</td><td>0.45</td><td>6.77%</td></tr><tr><td>Level 2</td><td>20</td><td>21.34</td><td>1.3</td><td>6.11%</td></tr><tr><td>Level 3</td><td>20</td><td>58.39</td><td>2.93</td><td>5.01%</td></tr></table> | 样品  | n     | Mean  | SD    | CV%   | Level 1 | 20      | 6.58 | 0.45 | 6.77% | Level 2 | 20      | 21.34 | 1.3   | 6.11% | Level 3 | 20      | 58.39 | 2.93  | 5.01% |       |
| 样品   | n   | Mean  | SD    | CV%   |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| Level 1  | 20  | 6.58  | 0.45  | 6.77% |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| Level 2  | 20  | 21.34 | 1.3   | 6.11% |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| Level 3  | 20  | 58.39 | 2.93  | 5.01% |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 样品类型   | Serum   |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 检测类型   | Sandwich (quantitative)   |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 灵敏度  | 0.25 ng/ml  |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 范围   | 5 ng/ml - 100 ng/ml   |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 实验步骤   | Multiple steps standard assay   |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 种属反应性  | 与反应: Human  |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |
| 产品概述   | Abcam's Prolactin <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Prolactin in Human serum.  |       |       |       |       |         |         |      |      |       |         |         |       |       |       |         |         |       |       |       |       |

Microtiterstrip wells are precoated with inactivated specific *Corynebacterium diphtheriae* toxin (toxoid) antigens to bind corresponding antibodies of the specimen. After washing the wells to remove all unbound sample material horseradish peroxidase (HRP) labelled anti-human IgG

conjugate is added. This conjugate binds to the captured C. diphtheriae toxin-specific antibodies. The immune complex formed by the bound conjugate is visualized by adding Tetramethylbenzidine (TMB) substrate which gives a blue reaction product. The intensity of this product is proportional to the amount of C. diphtheriae toxin-specific IgG antibodies in the specimen. Sulphuric acid is added to stop the reaction. This produces a yellow endpoint colour. Absorbance at 450 nm is read using an ELISA microwell plate reader.

**平台** Microplate

## 性能

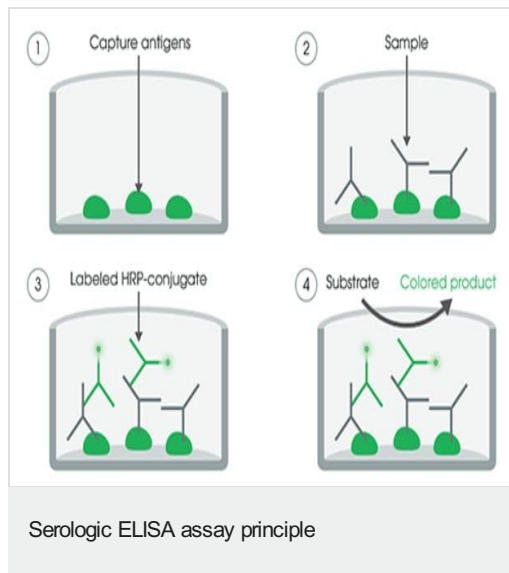
**存放说明** Store at +4°C. Please refer to protocols.

| 组件  | 标识符                  | 1 x 96 tests |
|---|----------------------|--------------|
| 10X Wash Solution                             |                      | 1 x 50ml     |
| Anti-Prolactin-HRP conjugate                  |                      | 1 x 12ml     |
| Cover foil                                    |                      | 1 unit       |
| Prolactin control                             |                      | 1 x 1ml      |
| Prolactin Standard 0: 0 ng/ml                 |                      | 1 x 1ml      |
| Prolactin Standard 1: 5 ng/ml                 |                      | 1 x 1ml      |
| Prolactin Standard 2: 10 ng/ml                |                      | 1 x 1ml      |
| Prolactin Standard 3: 25 ng/ml                |                      | 1 x 1ml      |
| Prolactin Standard 4: 50 ng/ml                |                      | 1 x 1ml      |
| Prolactin Standard 5: 100 ng/ml               |                      | 1 x 1ml      |
| Stop Solution                                 |                      | 1 x 15ml     |
| Streptavidin Coated Microplate (12 x 8 wells) | 12 Strips of 8 wells | 1 unit       |
| Strip holder                                  |                      | 1 unit       |
| TMB Substrate Solution                        |                      | 1 x 15ml     |

**相关性** Various hormones are secreted from the anterior pituitary during development and growth. Prolactin is a growth factor secreted by the anterior pituitary that is necessary for the proliferation and differentiation of the mammary glands. Prolactin acts primarily on the mammary gland by promoting lactation.

**细胞定位** Secreted

## 图片



Specific antigens are coated on the 96-well plate, controls or test samples are added to the well and incubated. The wells are washed to remove any unbound Human anti-antigen antibodies (Ig). A horseradish peroxidase (HRP) labelled anti-Human Ig conjugate is added to the wells. TMB is then catalyzed by the HRP to produce a blue color product that changes to yellow after adding an acidic stop solution. The intensity of yellow coloration is directly proportional to the amount of Human anti-antigen Ig captured on the plate.

**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