

Mouse IL-23 ELISA Kit ab119545

[3 References](#) [1 图像](#)

概述

产品名称 小鼠IL-23 ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

| 样品 | n | Mean | SD | CV% |
|---------|---|------|----|------|
| Overall | | | | 8.3% |

批次间

| 样品 | n | Mean | SD | CV% |
|---------|---|------|----|------|
| Overall | | | | 2.5% |

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 16 pg/ml

范围 31.3 pg/ml - 2000 pg/ml

回收率 特定样本回收率

| 样品类型 | 平均% | 范围 |
|-------|-----|-----------|
| Serum | 49 | 41% - 53% |

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 Abcam's IL-23 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse IL-23 in cell culture supernatant, serum and plasma (citrate).

IL-23 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated IL-23 detection antibody then incubated at room temperature. Following washing, a Streptavidin-HRP conjugate is added to each well, incubated at room temperature and washed. TMB is added and then catalyzed by HRP to produce a blue color product that changes into yellow after addition of acidic stop solution. The

density of yellow coloration is directly proportional to the amount of IL-23 captured on the plate.

平台 Microplate

性能

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

| 组件 | 1 x 96 tests |
|--|--------------|
| 20X Assay Buffer Concentrate | 1 x 5ml |
| 20X Wash Buffer Concentrate | 1 x 50ml |
| Adhesive Films | 4 units |
| Biotin-Conjugate anti-Mouse IL-23 monoclonal antibody | 1 x 70µl |
| IL-23 Standard lyophilized (4,000 pg/mL upon reconstitution) | 2 vials |
| Microplate coated with monoclonal antibody to Mouse IL-23 (12 x 8 wells) | 1 unit |
| Sample Diluent | 1 x 12ml |
| Stop Solution (1M Phosphoric acid) | 1 x 15ml |
| Streptavidin-HRP | 1 x 150µl |
| TMB Substrate Solution | 1 x 15ml |

功能 Associates with IL12B to form the IL-23 interleukin, an heterodimeric cytokine which functions in innate and adaptive immunity. IL-23 may constitute with IL-17 an acute response to infection in peripheral tissues. IL-23 binds to an heterodimeric receptor complex composed of IL12RB1 and IL23R, activates the Jak-Stat signaling cascade, stimulates memory rather than naive T-cells and promotes production of proinflammatory cytokines. IL-23 induces autoimmune inflammation and thus may be responsible for autoimmune inflammatory diseases and may be important for tumorigenesis.

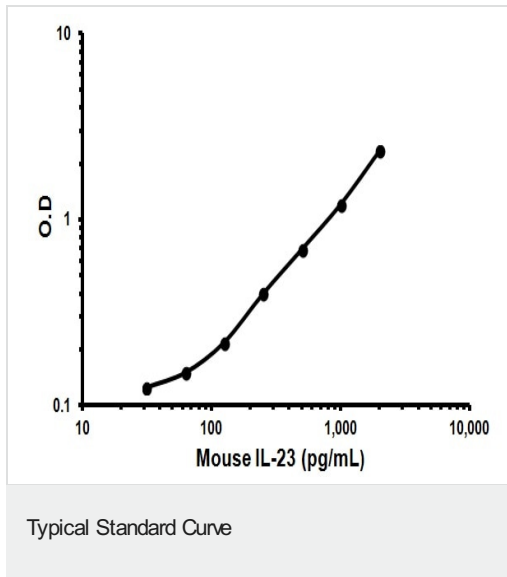
组织特异性 Secreted by activated dendritic and phagocytic cells and keratinocytes. Also expressed by dermal Langerhans cells (at protein level).

序列相似性 Belongs to the IL-6 superfamily.

发展阶段 Expressed by newborns dendritic cells.

细胞定位 Secreted. Secreted upon association with IL12B.

图片



Representative Standard Curve using ab119545.

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