

Human IL-3 ELISA Kit ab100569

2 图像

概述

| | |
|------|---|
| 产品名称 | 人IL-3 ELISA试剂盒 |
| 检测方法 | Colorimetric |
| 样品类型 | Cell culture supernatant, Serum, Plasma |
| 检测类型 | Sandwich (quantitative) |
| 灵敏度 | < 20 pg/ml |
| 范围 | 16.46 pg/ml - 12000 pg/ml |
| 回收率 | 100 % |

特定样本回收率

| 样品类型 | 平均 % | 范围 |
|--------------------------|-------|-------------|
| Cell culture supernatant | 118 | 110% - 127% |
| Serum | 114.4 | 103% - 122% |
| Plasma | 98.82 | 86% - 118% |

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's IL-3 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-3 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human IL-3 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-3 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of IL-3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

说明 Optimization may be required with urine samples.

平台 Microplate

性能

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

| 组件 | 1 x 96 tests |
|---|--------------|
| 20X Wash Buffer | 1 x 25ml |
| 5X Assay Diluent B | 1 x 15ml |
| 700X HRP-Streptavidin Concentrate | 1 x 200µl |
| Assay Diluent A | 1 x 30ml |
| Biotinylated anti-Human IL-3 | 2 vials |
| IL-3 Microplate (12 x 8 wells) | 1 unit |
| Recombinant Human IL-3 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

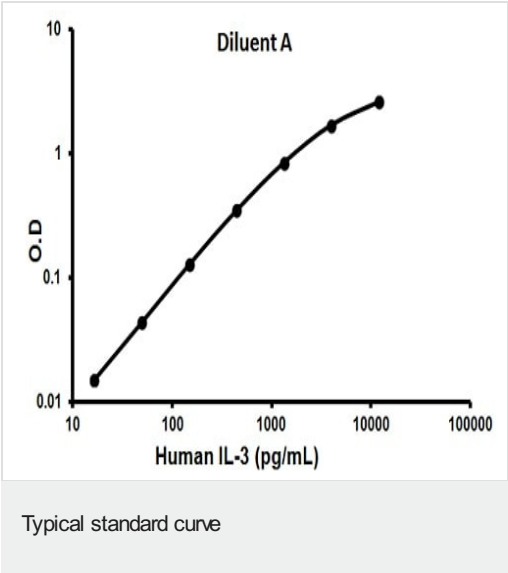
功能 Granulocyte/macrophage colony-stimulating factors are cytokines that act in hematopoiesis by controlling the production, differentiation, and function of 2 related white cell populations of the blood, the granulocytes and the monocytes-macrophages.
This CSF induces granulocytes, macrophages, mast cells, stem cells, erythroid cells, eosinophils and megakaryocytes.

组织特异性 Activated T-cells, mast cells, natural killer cells.

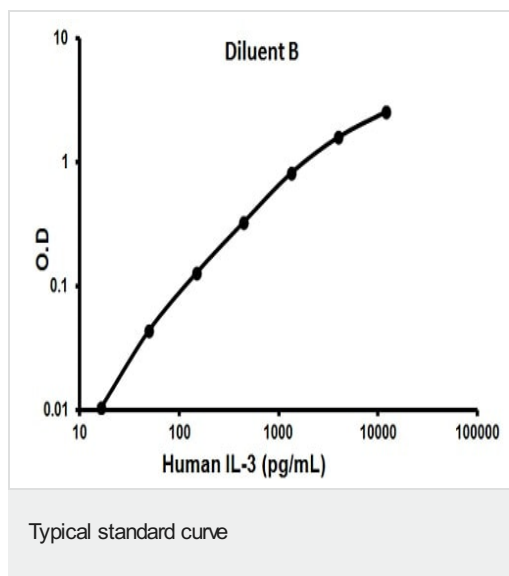
序列相似性 Belongs to the IL-3 family.

细胞定位 Secreted.

图片



Representative standard curve using ab100569



Representative standard curve using ab100569

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