

### Rat IL-10 ELISA Kit ab100765

★★★★★ [1 Abreviews](#) [25 References](#) [画像数 1](#)

#### 製品の概要

|         |                                        |   |     |       |
|---------|----------------------------------------|---|-----|-------|
| 製品名     | Rat IL-10 ELISA Kit                    |   |     |       |
| 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%   |
|         |                                        |   |     | < 10% |
|         | Inter-Assay (日差再現性)                    |   |     |       |
|         | サンプル                                   | N | 平均値 | SD    |
|         | Overall                                |   |     | CV%   |
|         |                                        |   |     | < 12% |
| サンプルの種類 | Cell culture extracts, Tissue Extracts |   |     |       |
| アッセイタイプ | Sandwich (quantitative)                |   |     |       |
| 検出感度    | < 10 pg/ml                             |   |     |       |
| 検出範囲    | 8.23 pg/ml - 6000 pg/ml                |   |     |       |
| 添加回収試験  | 140 %                                  |   |     |       |

|                       |              |             |
|-----------------------|--------------|-------------|
|                       | 特定サンプルでの回収試験 |             |
| サンプルの種類               | 平均 %         | 測定範囲        |
| Cell culture extracts | 142.1        | 133% - 149% |
| Tissue Extracts       | 138.7        | 127% - 146% |

|       |                                                                                                                                                                                                           |
|-------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| ステップ  | Multiple steps standard assay                                                                                                                                                                             |
| 種交差性  | 交差種: Rat                                                                                                                                                                                                  |
| 製品の概要 | Abcam's IL-10 Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of rat IL-10 in cell lysates and tissue lysates. |

This assay employs an antibody specific for rat IL-10 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-10 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-rat IL-10 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-10 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

#### 特記事項

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

#### 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 |
|----------------------------------------------|--------------|
| 140X HRP-Streptavidin Concentrate            | 1 x 200µl    |
| 20X Wash Buffer                              | 1 x 25ml     |
| 2X Cell Lysis Buffer                         | 1 x 5ml      |
| 5X Assay Diluent                             | 1 x 15ml     |
| 5X Sample Diluent Buffer                     | 1 x 10ml     |
| Biotinylated anti-Rat IL-10                  | 2 vials      |
| IL-10 Microplate (12 strips x 8 wells)       | 1 unit       |
| Recombinant rat IL-10 Standard (lyophilized) | 2 vials      |
| Stop Solution                                | 1 x 8ml      |
| TMB One-Step Substrate Reagent               | 1 x 12ml     |

#### 機能

Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

#### 組織特異性

Produced by a variety of cell lines, including T-cells, macrophages, mast cells and other cell types.

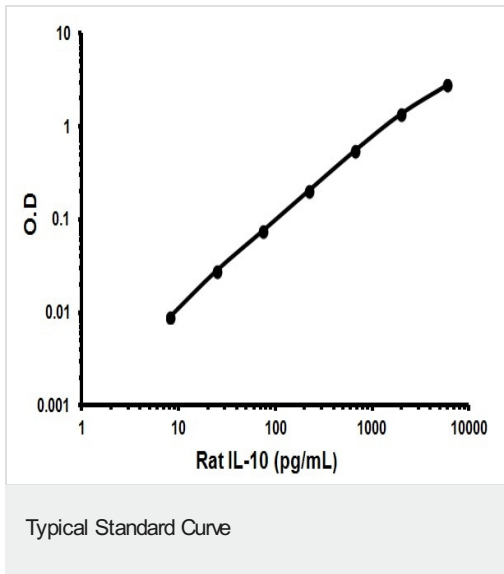
#### 配列類似性

Belongs to the IL-10 family.

#### 細胞内局在

Secreted.

#### 画像



Representative standard curve using ab100765.

**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