

Human CCL28 ELISA Kit ab99988

2 References 画像数 2

製品の概要

製品名	Human CCL28 ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 25 pg/ml
検出範囲	27.4 pg/ml - 20000 pg/ml
添加回収試験	> 65 %

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 supernatant	67.7	59% - 81%
Serum	70.6	60% - 98%
Plasma	68.6	59% - 82%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's CCL28 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human CCL28 in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human CCL28 coated on a 96-well plate. Standards and samples are pipetted into the wells and CCL28 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated antiHuman CCL28 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 CCL28 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
特記事項	Optimization may be required with urine samples.
試験プラットフォーム	Microplate

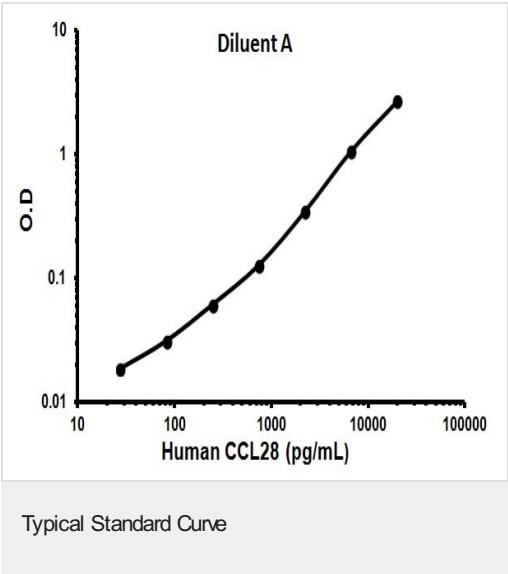
製品の特性

保存方法Store at -20°C. Please refer to protocols.

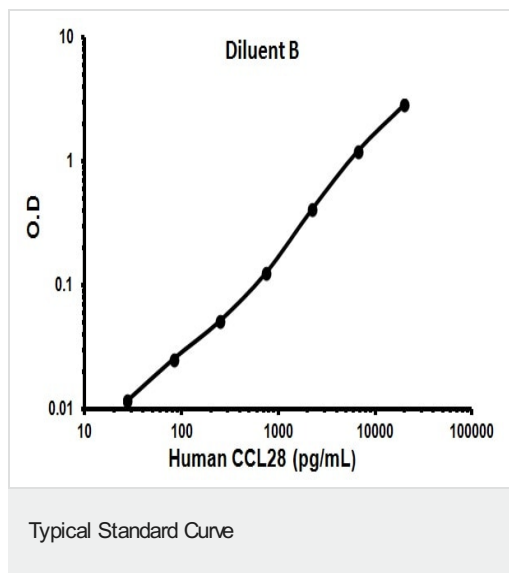
内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human CCL28	2 vials
CCL28 Microplate (12 strips x 8 wells)	1 x 200µl
Recombinant Human CCL28 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

機能	Chemotactic activity for resting CD4, CD8 T-cells and eosinophils. Binds to CCR3 and CCR10 and induces calcium mobilization in a dose-dependent manner.
組織特異性	Preferentially expressed by epithelial cells of diverse tissues including normal and pathological colon, salivary gland, mammary gland, trachea and rectum. Also found in prostate, spleen, thyroid, psoriasis skin and in lower levels in peripheral blood leukocytes, small intestine, Peyer patches, stomach and normal skin.
配列類似性	Belongs to the intercrine beta (chemokine CC) family.
細胞内局在	Secreted.

画像



Representative Standard Curve using ab99988



**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