

Human IL-15 ELISA Kit ab100554

★★★★☆ [1 Abreviews](#) [4 References](#) [画像数 2](#)

製品の概要

製品名	Human IL-15 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	< 3 pg/ml
検出範囲	3.3 pg/ml - 800 pg/ml
添加回収試験	94 %

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	96.37	83% - 104%
Serum	92.26	80% - 102%
Plasma	94.53	82% - 103%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's IL-15 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-15 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human IL-15 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-15 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IL-15 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-15 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 Optimization may be required with urine samples

製品の特性

保存方法

Store at -20°C. Please refer to protocols.

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

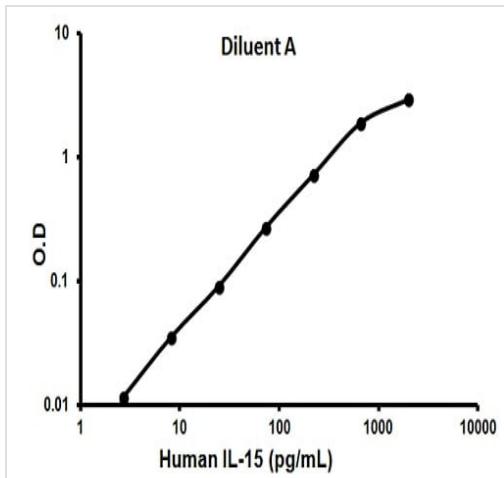
関連性

IL-15 (114 amino acids) is a cytokine that regulates T and natural killer cell activation and proliferation. It has a predicted molecular mass of approximately 12.5 kDa. Human IL-15 shares approximately 97% and 73% amino acid sequence identity with simian and mouse IL-15, respectively. Both human and simian IL-15 are active on mouse cells. IL-15 was initially isolated from the simian kidney epithelial cell line CV1/EBNA. It has also been isolated from mouse and human cell sources. The cytokines IL-15 and IL2 share many biological properties and stimulatory activities (T, B, and NK cells). Both IL-15 and IL-2 stimulate mouse CTLL2 cells. In activated peripheral blood T lymphocytes, IL-2 is highly expressed but the expression of IL-15 is not detectable. There is no sequence homology between IL-15 and IL-2, though computer modeling indicates both possess a four alpha helical bundle structure. IL-15 competes for binding sites with IL-2, as both IL2 and IL-15 stimulate the growth of cells through the IL-2 receptor. IL-15 mRNA is expressed in many cell types and tissues including adherent peripheral blood mononuclear cells, fibroblasts, and epithelial cells, monocytes, placenta, and skeletal muscle.

細胞内局在

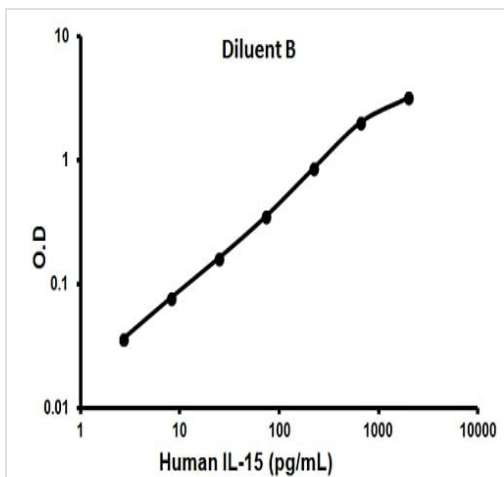
Isoform IL-15-S48AA: Secreted. Isoform IL-15-S21AA: Cytoplasm. Nucleus. Note: IL-15-S21AA is not secreted, but rather is stored intracellularly, appearing in the nucleus and cytoplasmic components.

画像



Representative standard curve using ab100554

Typical standard curve



Representative standard curve using ab100554

Typical standard curve

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