

Mouse IL-15 ELISA Kit ab100701

1 References 1 Image

Overview

Product name	Mouse IL-15 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 0.5 ng/ml
Range	0.33 ng/ml - 80 ng/ml
Recovery	> 100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	137.2	115% - 150%
Plasma	128.5	105% - 148%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mouse
Product overview	

Abcam's IL-15 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse IL-15 in plasma and cell culture supernatants. (Mouse IL-15 concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for mouse 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-mouse 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.

We have not been able to detect endogenous Mouse IL-15 in normal serum with ab100701, only in serum spiked with Mouse IL-15.

**Platform**

Microplate

**Properties****Storage instructions**

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated anti-mouse IL-15	2 vials
IL-15 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IL-15 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

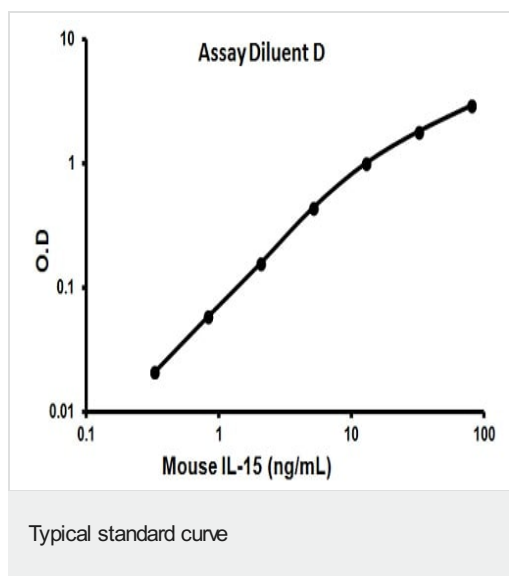
**Relevance**

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.

**Cellular localization**

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.

**Images**



Representative standard curve using ab100701

**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.com/abpromise> or contact our technical team.

#### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors