

Mouse IL-11 ELISA Kit ab100698

1 References 1 Image

Overview

Product name	Mouse IL-11 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 1 ng/ml
Range	1.85 ng/ml - 450 ng/ml
Recovery	99 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	77.35	70% - 87%
Plasma	117.3	95% - 113%

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

Abcam's IL-11 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse IL-11 in plasma and cell culture supernatants. (Mouse IL-11 concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for mouse IL-11 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-11 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse IL-11 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-11 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 the endogenous Mouse IL-11 in normal serum with ab100698, only in serum spiked with Mouse IL-11.

Platform	Microplate
----------	------------

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

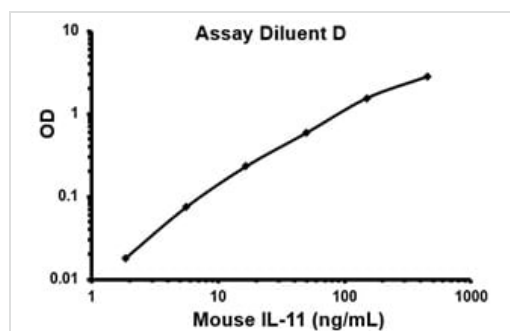
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
5X Assay Diluent D	1 x 15ml
Biotinylated anti-mouse IL-11	2 vials
IL-11 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IL-11 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Function** Directly stimulates the proliferation of hematopoietic stem cells and megakaryocyte progenitor cells and induces megakaryocyte maturation resulting in increased platelet production.

**Sequence similarities** Belongs to the IL-6 superfamily.

**Cellular localization** Secreted.

## Images



Typical standard curve

Representative standard curve using ab100698.

**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