

Product datasheet

Mouse IL-11 ELISA Kit ab215084

Recombinant SimpleStep ELISA

[5 Images](#)

Overview

Product name Mouse IL-11 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	8			5.4%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	3			10.3%

Sample type Cell culture supernatant, Serum, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity 17.4 pg/ml

Range 78.12 pg/ml - 5000 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	102	98% - 105%
Serum	100	98% - 101%
Cit plasma	91	86% - 97%

Assay time 1h 30m

Assay duration One step assay

Species reactivity

Reacts with: Mouse

Does not react with: Cow

Product overview

Mouse IL-11 ELISA Kit (ab215084) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of IL-11 protein in cell culture supernatant, cit plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse IL-11 with 17.4 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes

IL-11 is a multifunctional cytokine that is a key regulator of hematopoiesis. It stimulates hematopoietic stem cell and megakaryocyte progenitor cell proliferation. It is also involved in the response to liver damage by inducing proliferation of hepatocytes.

Platform

Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPR	1 x 6ml
10X Mouse IL-11 Capture Antibody	1 x 600µl
10X Mouse IL-11 Detector Antibody	1 x 600µl
Mouse IL-11 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
Stop Solution	1 x 12ml
TMB Development Solution	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.

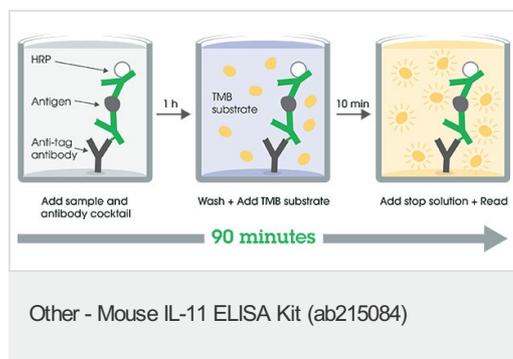
Sequence similarities

Belongs to the IL-6 superfamily.

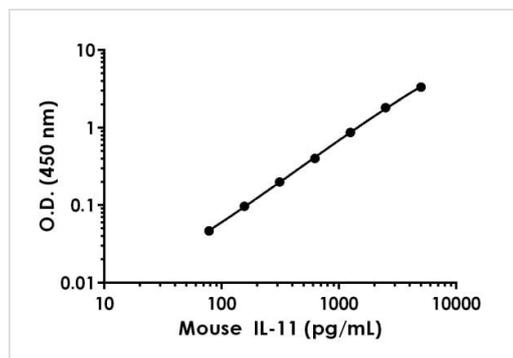
Cellular localization

Secreted.

Images

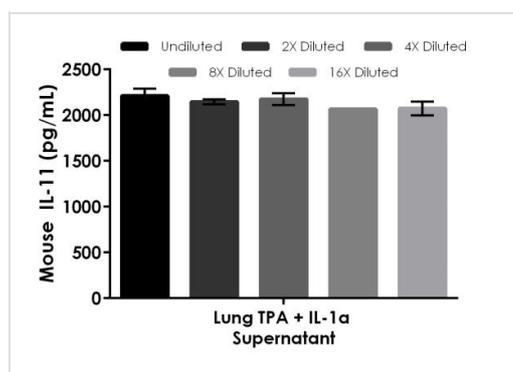


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



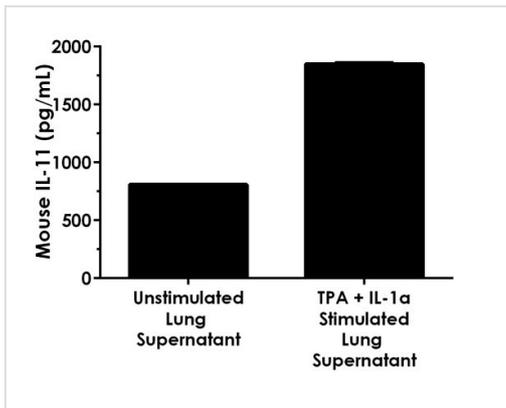
Background-subtracted data values (mean +/- SD) are graphed.

Example of mouse IL-11 standard curve in Sample Diluent NS.



The concentrations of IL-11 were measured in duplicates and interpolated from the IL-11 standard curves. Undiluted samples are as follows: lung supernatant stimulated with TPA and IL-1a 100%. The interpolated values are plotted (mean +/- SD, n=2). The mean IL-11 concentration was determined to be 2143 pg/mL in lung supernatant stimulated with TPA and IL-1a.

Interpolated concentrations of native IL-11 in mouse lung cell culture supernatant samples.



Mouse lung cells were cultured unstimulated or stimulated with 10 ng/mL TPA and 1 ng/mL IL-1 α .

Conditioned media was harvested after 48 hours. IL-11 was measured in 100% unstimulated and stimulated lung supernatant. The concentrations of IL-11 were measured in duplicate, interpolated from the IL-11 standard curves. The interpolated values are plotted (mean \pm SD, n=2). The mean IL-11 concentration was determined to be 1849 pg/mL in TPA and IL-1 α stimulated lung supernatant and 809 pg/mL in unstimulated lung supernatant.

Powered by recombinant antibodies

 <p>Research with confidence Consistent and reproducible results</p>	 <p>Long-term and scalable supply Recombinant technology</p>
 <p>Success from the first experiment Confirmed specificity</p>	 <p>Ethical standards compliant Animal-free production</p>

Sandwich ELISA - Mouse IL-11 ELISA Kit
(ab215084)

To learn more about the advantages of recombinant antibodies see [here](#).

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