

Product datasheet

Human IL-4 ELISA Kit ab215089

Recombinant SimpleStep ELISA

★★★★★ 1 Abreviews 12 References 8 Images

Overview

Product name	Human IL-4 ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall	5			6%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall	3			7%
Sample type	Cell culture supernatant, Serum, Plasma, Hep Plasma, EDTA Plasma, Cit plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	1.08 pg/ml				
Range	6.25 pg/ml - 400 pg/ml				
Recovery	Sample specific recovery				
	Sample type		Average %		Range
	Serum		95		91% - 102%
	Cell culture media		117		112% - 120%
	Hep Plasma		95		88% - 98%
	EDTA Plasma		109		103% - 114%
	Cit plasma		100		94% - 104%
Assay time	1h 30m				
Assay duration	One step assay				

Species reactivity**Reacts with:** Human**Product overview**

Human IL-4 ELISA Kit (ab215089) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of IL-4 protein in cit plasma, edta plasma, hep plasma, plasma, serum, and cell culture supernatant. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human IL-4 with 1.08 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-4 is a 129 amino acids long secreted protein. IL-4 is a cytokine produced mostly by activated T lymphocytes, mast cells and basophils. It has variety of immune response modulating functions by acting on number of cell types. IL-4 participates in several B-cell activation processes. It is a co-stimulator of DNA-synthesis. It induces the expression of class II MHC molecules on resting B-cells. It enhances both secretion and cell surface expression of IgE and IgG1. It also regulates the expression of the low affinity Fc receptor for IgE (CD23) on both lymphocytes and monocytes. IL-4 positively regulates IL31RA expression in macrophages.

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 5BI	1 x 6ml
10X Human IL-4 Capture Antibody	1 x 600µl
10X Human IL-4 Detector Antibody	1 x 600µl
Human IL-4 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent 75BS	1 x 20ml

Components	1 x 96 tests
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

Function

Participates in at least several B-cell activation processes as well as of other cell types. It is a costimulator of DNA-synthesis. It induces the expression of class II MHC molecules on resting B-cells. It enhances both secretion and cell surface expression of IgE and IgG1. It also regulates the expression of the low affinity Fc receptor for IgE (CD23) on both lymphocytes and monocytes.

Involvement in disease

Genetic variations in IL4 may be a cause of susceptibility to ischemic stroke (ISCHSTR) [MIM:601367]; also known as cerebrovascular accident or cerebral infarction. A stroke is an acute neurologic event leading to death of neural tissue of the brain and resulting in loss of motor, sensory and/or cognitive function. Ischemic strokes, resulting from vascular occlusion, is considered to be a highly complex disease consisting of a group of heterogeneous disorders with multiple genetic and environmental risk factors.

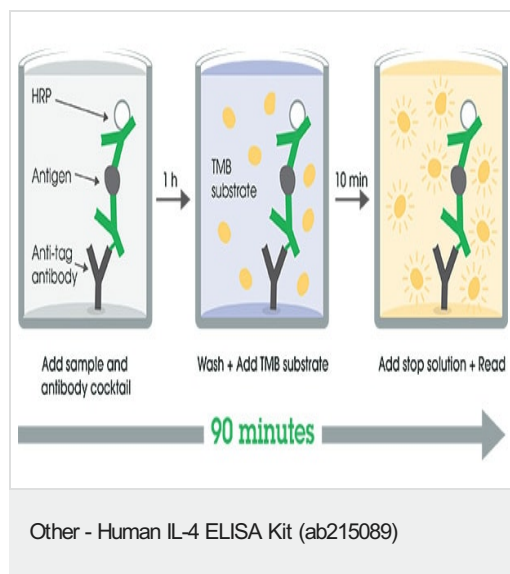
Sequence similarities

Belongs to the IL-4/IL-13 family.

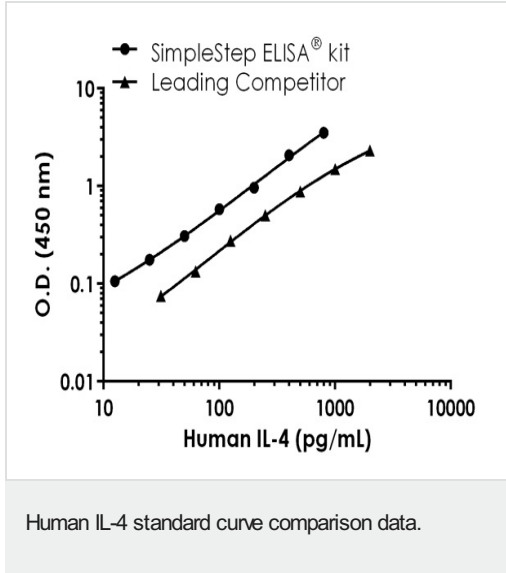
Cellular localization

Secreted.

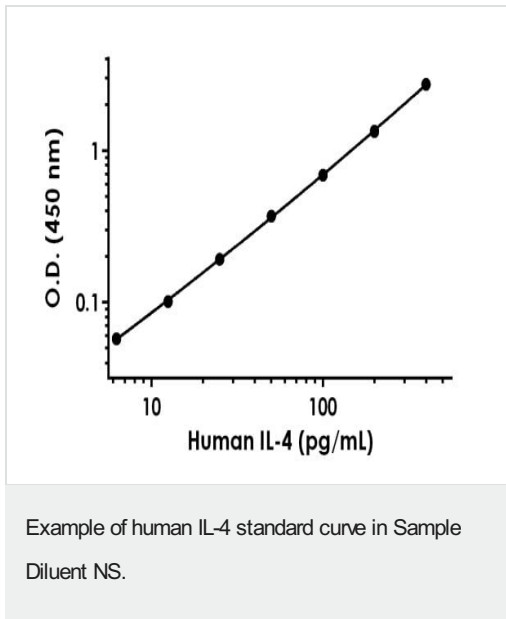
Images



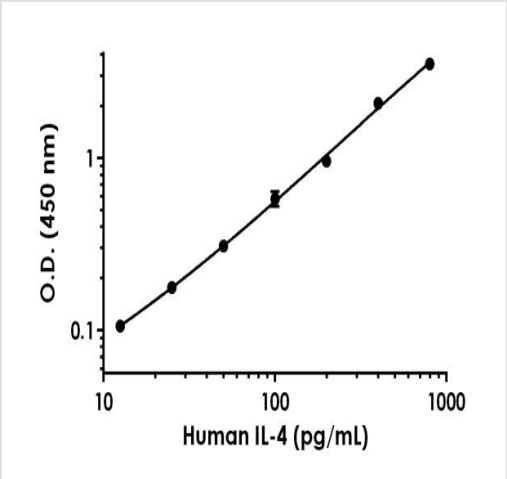
SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



Standard curve comparison between human IL-4 SimpleStep ELISA[®] kit and traditional ELISA kit from leading competitor. SimpleStep ELISA kit shows a 8-fold increase in sensitivity.

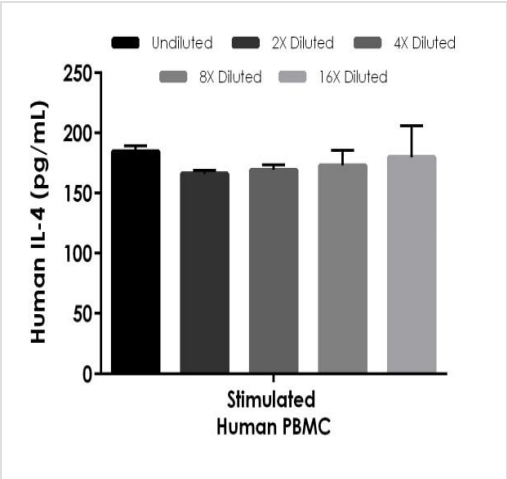


Background-subtracted data values (mean +/- SD) are graphed.



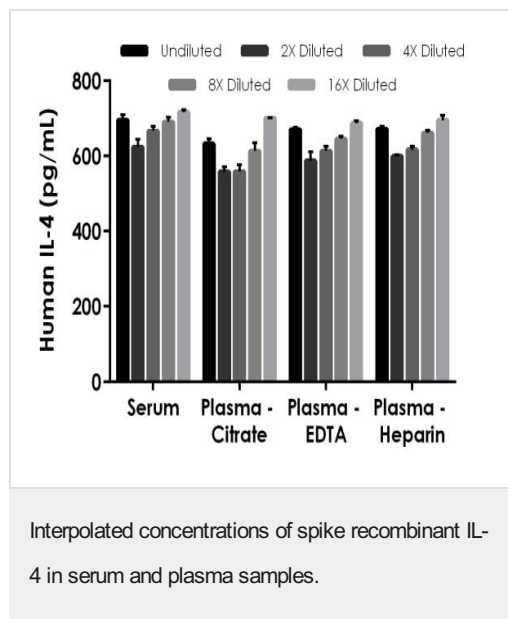
Example of human IL-4 standard curve in Sample Diluent 75BS.

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

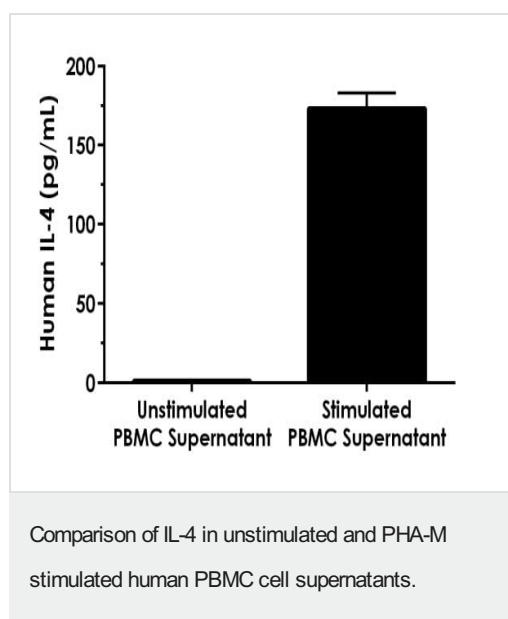


Interpolated concentrations of native IL-4 in PHA-M stimulated human PBMC cell culture supernatant samples.

The concentrations of IL-4 were measured in duplicates, interpolated from the IL-4 standard curves and corrected for sample dilution. Undiluted samples are as follows: stimulated human PBMC cell culture supernatant 100%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean IL-4 concentration was determined to be 174.6 pg/mL in stimulated human PBMC cell culture supernatant.



The concentrations of IL-4 were measured in duplicates, interpolated from the IL-4 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 100%, plasma citrate 100%, plasma EDTA 100%, plasma heparin 100%. The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$).



Human PBMC cells were cultured in the absence or presence of 1.5% PHA-M for 46 hours. The concentrations of IL-4 were measured in neat supernatant samples in duplicates and interpolated from the IL-4 standard curve. The interpolated values are plotted (mean \pm SD, $n=2$). The mean IL-4 concentration was determined to be 174.6 pg/mL in PHA-M stimulated PBMC cell supernatant, 1.4 pg/mL in unstimulated supernatants and undetectable in media (not shown).

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Human IL-4 ELISA Kit (ab215089)

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