

### Human IL-2 ELISA Kit ab174444

SimpleStep ELISA

[9 References](#) [7 Images](#)

#### Overview

**Product name** Human IL-2 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall	6			6%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	24			6%

**Sample type**

Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

**Assay type**

Sandwich (quantitative)

**Sensitivity**

9 pg/ml

**Range**

15.63 pg/ml - 2000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	95	93% - 97%
Cell culture media	94	83% - 109%
Hep Plasma	91	82% - 95%
EDTA Plasma	100	95% - 105%
Cit plasma	83	77% - 98%

**Assay time**

1h 30m

**Assay duration**

One step assay

## Species reactivity

## Product overview

**Reacts with:** Human

Human IL-2 ELISA kit has been re-developed. We have identified new recombinant monoclonal antibodies to provide improved performance and consistency. This version will be discontinued when inventory is depleted. The new version is available as [ab270883](#).

Human IL-2 ELISA kit (ab174444) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of IL-2 protein in human cell culture supernatant, plasma and serum samples. It uses our proprietary SimpleStep ELISA® technology. Quantitate human IL-2 with 9 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.

## ASSAY SPECIFICITY

This kit recognizes both native and recombinant Human IL-2 protein in serum, plasma, and cell culture supernatant samples only.

Extract samples have not been tested with this kit.

## SPECIES REACTIVITY

This kit recognizes Human IL-2 protein.

No cross reaction observed with recombinant mouse IL-2 tested at 10 ng/mL.

Other species reactivity not determined.

## Notes

IL-2, also known as T cell growth factor (TCGF), is a glycosylated alpha-helical polypeptide, synthesized as a 153 amino acid (aa) precursor with a 20 aa signal peptide and a 133 aa mature chain. It is secreted by activated CD4+ and CD8+ T cells, neurons, microglia and hematopoietic stem cells in response to antigenic or mitogenic stimulation. IL-2 is required for T-cell proliferation, Natural Killer cells (NK) cytolytic activity, differentiation of regulatory T cells, modulation of T helper (Th) cell differentiation and activation-induced cell death. In particular, IL-2 modulates the expression of receptors for other cytokines and transcription factors, therefore regulating cytokine cascades that correlate with each of the Th differentiation states.

Complete deficiency of IL-2 has been implicated in severe combined immunodeficiency, whereas reduction of the IL-2 correlates with reduced function of CD4+CD25+ regulatory T cells and destabilization of immune homeostasis leading to autoimmune disease. Increased expression of IL-2 has also been implicated in inflammatory conditions such as inflammatory bowel disease and chronic liver diseases. IL-2 therefore is both a immune stimulator and immune suppressor cytokine which efficiently controls the immune system to deal with autoimmunity and adaptive immune response.

**Platform** Microplate

**Properties**

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

Components	
Plate Seals	1 x unit
10X Wash Buffer PT ( <a href="#">ab206977</a> )	1 x 20ml
TMB&nbsp;Development Solution	1 x 12ml
Stop Solution	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate ( <a href="#">ab206978</a> )	1 x unit
10X Human IL-2 Detector Antibody	1 x 600µl
Human IL-2 Lyophilized Recombinant Protein	2 x 1vial
10X Human IL-2 Capture Antibody	1 x 600µl
Sample Diluent NS ( <a href="#">ab193972</a> )	1 x 50ml
Sample Diluent 25BS	1 x 20ml
Antibody Diluent 4BI	1 x 6ml

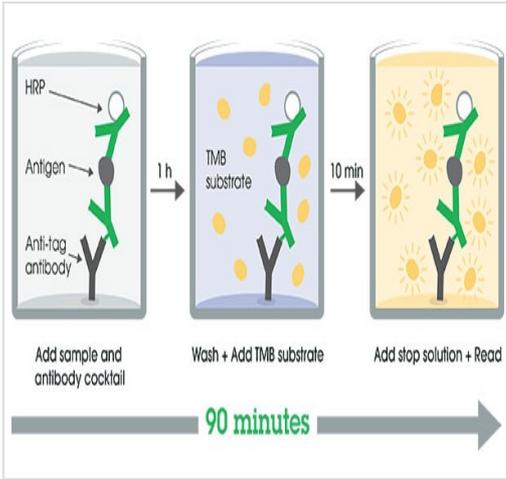
**Function** Produced by T-cells in response to antigenic or mitogenic stimulation, this protein is required for T-cell proliferation and other activities crucial to regulation of the immune response. Can stimulate B-cells, monocytes, lymphokine-activated killer cells, natural killer cells, and glioma cells.

**Involvement in disease** Note=A chromosomal aberration involving IL2 is found in a form of T-cell acute lymphoblastic leukemia (T-ALL). Translocation t(4;16)(q26;p13) with involves TNFRSF17.

**Sequence similarities** Belongs to the IL-2 family.

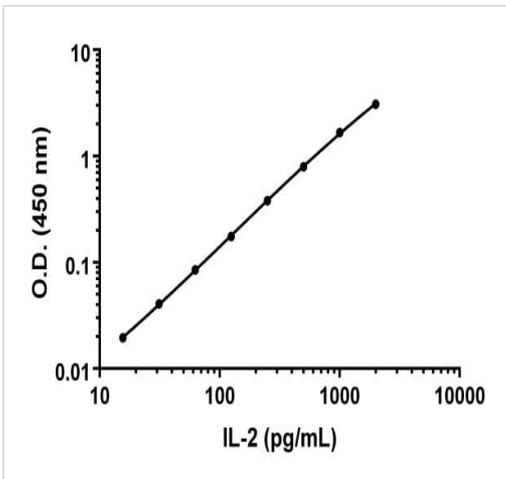
**Cellular localization** Secreted.

**Images**



Other - Human IL-2 ELISA Kit (ab174444)

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.



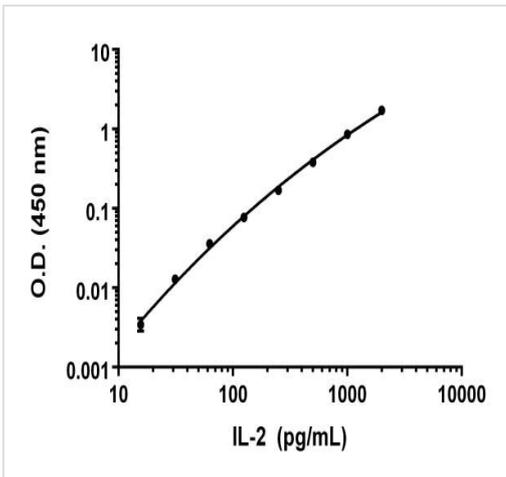
Example IL-2 standard curve in sample diluents NS.

Example IL-2 standard curve in sample diluents NS.

Standard Curve Measurements			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.061	0.059	0.0602
15.63	0.088	0.087	0.088
31.3	0.112	0.106	0.109
62.5	0.164	0.162	0.163
125	0.285	0.246	0.265
250	0.556	0.482	0.519
500	1.081	0.908	0.995
1,000	1.938	2.014	1.976
2,000	3.559	3.548	3.553

The IL-2 standard curve was prepared as described. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example IL-2 standard curve in Sample Diluents NS.



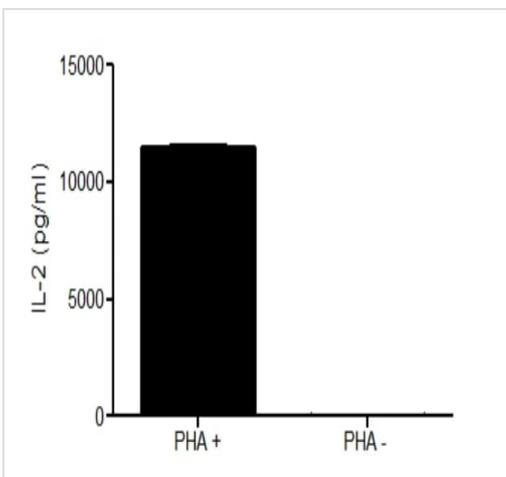
Example IL-2 standard curve in standard diluent 25BS.

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

Standard Curve Measurements			
Conc. (pg/mL)	O.D. 450 nm		Mean O.D.
	1	2	
0	0.058	0.057	0.058
15.63	0.061	0.061	0.061
31.3	0.069	0.072	0.070
62.5	0.093	0.094	0.093
125	0.136	0.133	0.134
250	0.232	0.222	0.227
500	0.439	0.433	0.436
1,000	0.912	0.914	0.913
2,000	1.720	1.811	1.765

Example IL-2 standard curve in Sample Diluent 25BS.

The curve was prepared by loading standard curve as described. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Specificity of IL-2 signal on stimulated and non stimulated media supernatants.

Human PBMCs were cultured in RPMI supplemented with 10% fetal calf serum, 2mM L-glutamine, 100U/mL penicillin, and 100 µg/mL streptomycin. Cells were cultured for 2 days at 37°C in the presence or absence of PHA. The concentrations of IL-2 were interpolated from the calibration curve and corrected for sample dilution. The mean IL-2 concentration was determined to be 16 pg/mL in unstimulated PBMC supernatants and 11,460 pg/mL in stimulated PBMC supernatants.

	Interpolated value	NHS 50%	NHP Citrate 50%	NHP EDTA 50%	NHP Heparin 50%	Media 10%
1:1	pg/mL	2,065	1,893	1,959	1,802	2,214
	% Expected value	100	100	100	100	100
1:2	pg/mL	1,010	964	964	920	1,061
	% Expected value	102	102	98	102	105
1:4	pg/mL	543	532	543	510	498
	% Expected value	112	113	111	113	98
1:8	pg/mL	293	293	305	281	243
	% Expected value	124	125	125	124	94

Linearity of dilution.

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Recombinant Human IL-2 was spiked into the following biological samples and diluted in a 2-fold dilution series in Sample Diluent NS.

**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