

Human IL-10 ELISA Kit (Interleukin-10) ab185986

KO 評価済 リコンビナント SimpleStep ELISA

[19 References](#) [画像数 11](#)

製品の概要

製品名 Human IL-10 ELISA Kit (Interleukin-10)

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
supernatant	5			5.2%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
supernatant	3			3.2%

サンプルの種類 Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 1.4 pg/ml

検出範囲 9.4 pg/ml - 3000 pg/ml

添加回収試験 特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	108	105% - 111%
Serum	109	108% - 112%
Cell culture media	106	102% - 111%
Hep Plasma	116	113% - 117%
EDTA Plasma	109	103% - 118%
Cit plasma	111	105% - 119%

全工程の試験時間	1h 30m
ステップ	One step assay
種交差性	交差種: Human 非交差種: Cow
製品の概要	<p>Human IL-10 ELISA Kit (Interleukin-10) (ab185986) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of IL-10 (Interleukin-10) protein in cell culture supernatant, cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human IL-10 (Interleukin-10) with 1.4 pg/ml sensitivity.</p> <p>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:</p> <ul style="list-style-type: none"> - 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 <p>A 384-well SimpleStep ELISA® microplate (ab203359) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.</p>
特記事項	Interleukin 10 (IL-10) is a cytokine involved in immunoregulation through the promotion of Th2 response and inhibition of Th1 responses. It is produced by cells of hematopoietic origin, keratinocytes, placental cytotrophoblasts, and hepatic stellate cells.
試験プラットフォーム	Pre-coated microplate (12 x 8 well strips)

製品の特性

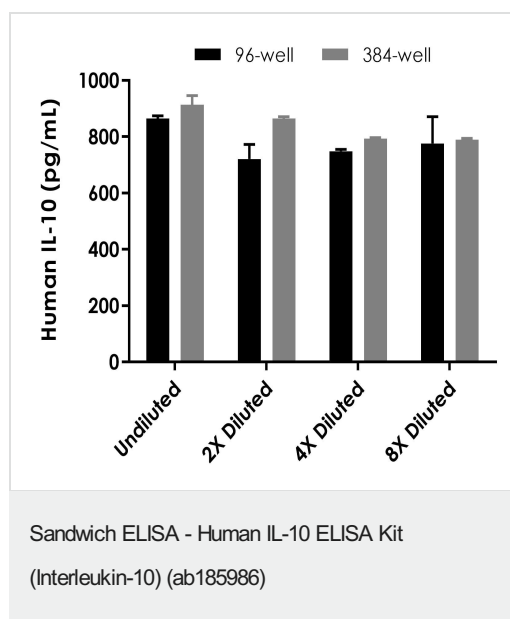
保存方法 Store at +4°C. Please refer to protocols.

内容	1 x 96 tests	1 x 384 tests	1 x 96 tests
10X Human IL-10 Capture Antibody	1 x 600µl	1 x 600µl	1 x 600µl
10X Human IL-10 Detector Antibody	1 x 600µl	1 x 600µl	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 20ml	1 x 20ml
384 well CaptSure™ microplates	0 x 0 unit	1 unit	0 x 0 unit
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml	6 x 1ml	1 x 1ml
Antibody Diluent CPI2	1 x 6ml	1 x 6ml	1 x 6ml
Human IL-10 Lyophilized Recombinant Protein	2 vials	2 vials	2 vials

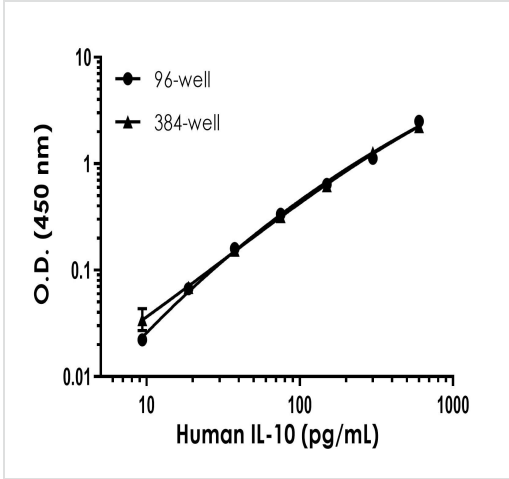
内容	1 x 96 tests	1 x 384 tests	1 x 96 tests
Plate Seals	1 unit	1 unit	1 unit
Sample Diluent NBP	1 x 20ml	1 x 20ml	1 x 20ml
Sample Diluent NS (ab193972)	1 x 50ml	1 x 250ml	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	0 x 0 unit	1 unit
Stop Solution	1 x 12ml	2 x 12ml	1 x 12ml
TMB Development Solution	1 x 12ml	2 x 12ml	1 x 12ml

機能	Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.
組織特異性	Produced by a variety of cell lines, including T-cells, macrophages, mast cells and other cell types.
配列類似性	Belongs to the IL-10 family.
細胞内局在	Secreted.

画像

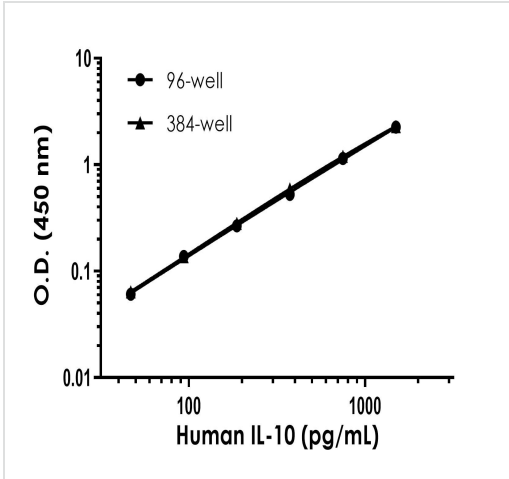


Interpolated concentration of native IL-10 was measured in duplicate at different sample concentrations in 96-well vs. 384-well plates. Undiluted samples are 50% stimulated PBMC supernatant (1.5% PHA). The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). Sample dilutions are made in Sample Diluent NS + 1X Enhancer.



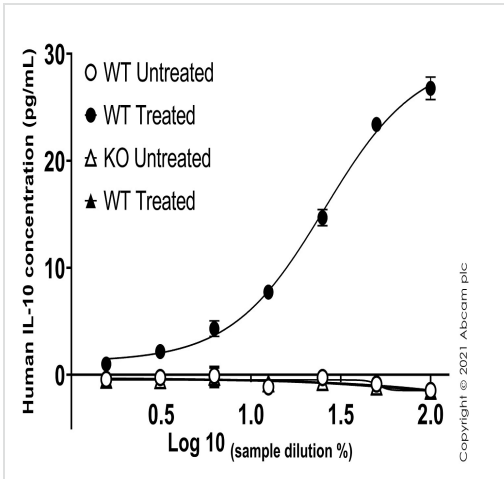
Example of human IL-10 standard curve in 96-well vs. 384-well plate. Background-subtracted data values (mean +/- SD) are graphed.

Sandwich ELISA - Human IL-10 ELISA Kit
(Interleukin-10) (ab185986)



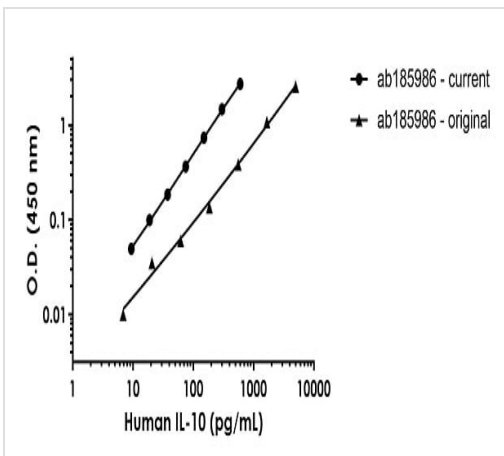
Example of human IL-10 standard curve in 96-well vs. 384-well plate. Background-subtracted data values (mean +/- SD) are graphed.

Sandwich ELISA - Human IL-10 ELISA Kit
(Interleukin-10) (ab185986)



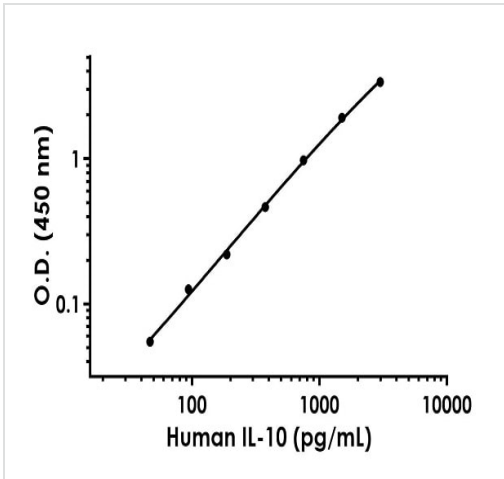
KO data

Human IL-10 concentration was interpolated from the IL-10 standard curve. Supernatants from cell culture samples were serially diluted and assessed by the Human IL-10 ELISA Kit (Interleukin-10) (ab185986). Wild-type and IL-10 knockout THP-1 ([ab273763](#)) vehicle control and PMA (100 ng/mL, 56 h) + LPS (1 µg/mL, 24 h) treated cells were assessed in duplicate (n=2). Data are represented as the mean and error bars represent standard deviation.”



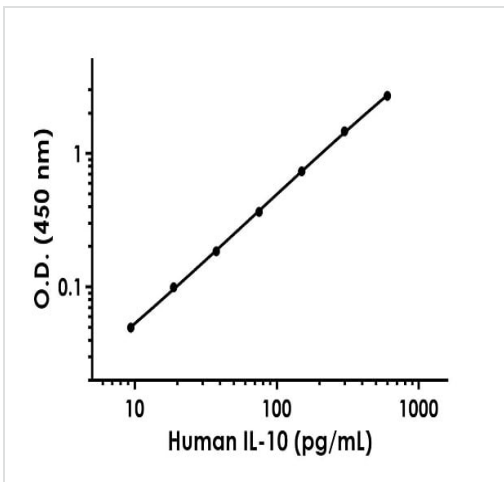
Human IL-10 Standard Curve Comparison

Standard Curve comparison between human IL-10 current (new) SimpleStep ELISA kit and original ELISA kit. The current SimpleStep ELISA kit shows increased sensitivity.



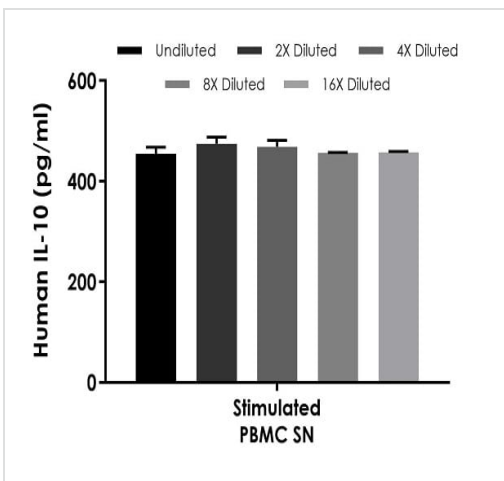
Example of human IL-10 standard curve in Sample Diluent 50BP + Enhancer.

The IL-10 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



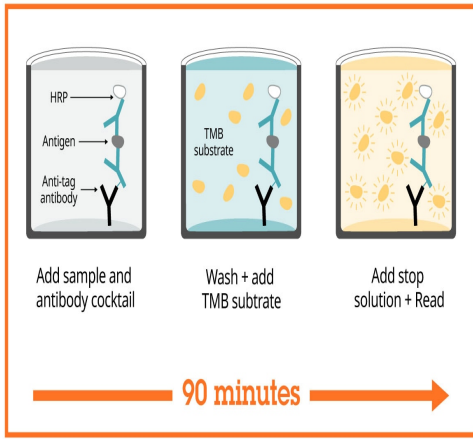
Example of human IL-10 standard curve in Sample Diluent NS + Enhancer.

The IL-10 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Interpolated concentrations of native IL-10 in human PBMC cell culture supernatant samples.

The concentrations of IL-10 were measured in duplicates, interpolated from the IL-10 standard curves and corrected for sample dilution. Undiluted samples are as follows: stimulated PBMC cell culture supernatant 50%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean IL-10 concentration was determined to be 462 pg/ml in stimulated PBMC cell culture supernatant. PBMC cell culture supernatant samples were stimulated for 46 hours with 1.5% PHA. Unstimulated PBMC cell culture supernatant samples were also measured; all values were below the detectable range of the assay.



Sandwich ELISA - Human IL-10 ELISA Kit
(Interleukin-10) (ab185986)

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.

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-10 ELISA Kit
(Interleukin-10) (ab185986)

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

Get more done with
SimpleStep ELISA

- Easy to use**
Single-wash 90-minute protocol
- Flexible**
Matched antibody pairs available
- Precision antibodies**
High sensitivity, specificity and reproducibility
- Scalable**
Now in 10-pack and 384-well formats

Sandwich ELISA - Human IL-10 ELISA Kit
(Interleukin-10) (ab185986)

To learn more about the advantages of SimpleStep ELISA[®] kits 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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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