invitrogen USER GUIDE

# Human IL-10 High Sensitivity ELISA Kit

Enzyme-linked Immunosorbent Assay for quantitative detection of human IL-10

Catalog Number BMS215HS

Pub. No. MAN0016564 Rev. D.0 (33)



**WARNING!** Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

# Product description

The Human IL-10 High Sensitivity ELISA Kit is an enzyme-linked immunosorbent assay for the quantitative detection of human IL-10.

### Summary

Interleukin-10 is a pleiotropic cytokine playing an important role as a regulator of lymphoid and myeloid cell function. Due to the ability of IL-10 to block cytokine synthesis and several accessory cell functions of macrophages this cytokine is a potent suppressor of the effector functions of macrophages, T-cells and NK cells. In addition, IL-10 participates in regulating proliferation and differentiation of B-cells, mast cells and thymocytes.

The primary structure of human IL-10 has been determined by cloning the cDNA encoding the cytokine. The corresponding protein exerts 160 amino acids with a predicted molecular mass of 18.5 kDa. Based on its primary structure, IL-10 is a member of the four-helix bundle family of cytokines. In solution human IL-10 is a homodimer with an apparent molecular mass of 39 kDa. Although it contains an N-linked glycosylation site, it lacks detectable carbohydrates. Recombinant protein expressed in E. coli thus retains all known biological activities.

The human IL-10 gene is located on chromosome 1 and is present as a single copy in the genome.

The human IL-10 exhibits strong DNA and amino acid sequence homology to the murine IL-10 and an open reading frame in the Epstein-Barr virus genome, BCRF1, which shares many of the cellular cytokine's biological activities and may therefore play a role in the host-virus interaction.

The immunosuppressive properties of IL-10 suggest a possible clinical use of IL-10 in suppressing rejections of grafts after organ transplantations. IL-10 can furthermore exert strong anti-inflammatory activities.

IL-10 in disease: IL-10 expression was shown to be elevated in parasite infections like in Schistosoma mansoni, Leishmania, Toxoplasma gondii, and Trypanosoma infection. Furthermore, high IL-10 expression was detected in mycobacterial infections as shown for Mycobacterium leprae, Mycobacterium tuberculosis, and Mycobacterium avium infections. High expression levels of IL-10 are also found in retroviral infections inducing immunodeficiency.

For literature update refer to our website.

#### Principles of the test

An anti-human IL-10 coating antibody is adsorbed onto microwells.

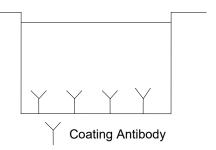


Fig. 1 Coated microwell

Human IL-10 present in the sample or standard binds to antibodies adsorbed to the microwells. A biotin-conjugated anti-human IL-10 antibody is added and binds to human IL-10 captured by the first antibody.

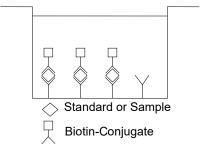


Fig. 2 First incubation

Following incubation unbound biotin-conjugated anti-human IL-10 antibody is removed during a wash step. Streptavidin-HRP is added and binds to the biotin-conjugated anti- human IL-10 antibody.

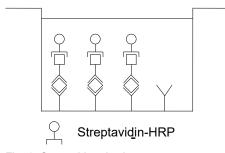


Fig. 3 Second incubation

Following incubation unbound Streptavidin-HRP is removed during a wash step, and amplification reagent I (BiotinyI-Tyramide) is added to the wells.

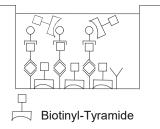
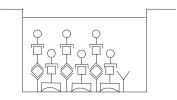


Fig. 4 Third incubation

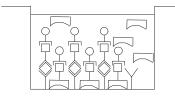
Following incubation unbound amplification reagent I is removed during a wash step and amplification reagent II (Streptavidin-HRP) is added.



Streptavidin-HRP

Fig. 5 Fourth incubation

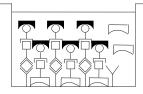
Following incubation unbound amplification reagent II is removed during a wash step and substrate solution reactive with HRP is added.



Substrate

Fig. 6 Fifth incubation

A colored product is formed in proportion to the amount of human IL-10 present in the sample or standard. The reaction is terminated by addition of acid and absorbance is measured at 450 nm. A standard curve is prepared from 7 human IL-10 standard dilutions and human IL-10 sample concentration determined.



Reacted Substrate

Fig. 7 Stop reaction

### Principle of amplification reaction

The amplification reaction is based upon PerkinElmer Life Sciences  $TSA^{\text{TM}}$  technology.

Amplification reagent I contains biotinyl-tyramide. HRP converts multiple biotinyl-tyramide molecules into highly reactive derivates (free radicals). These free radicals bind covalently to any protein in the well.

Thus, the amount of reacted biotinyl-tyramide is proportional to the amount of HRP in the well.

Following incubation unbound biotinyl-tyramide is removed during a wash step. Amplification reagent II contains Streptavidin-HRP, which binds to the biotin sites created during the biotinyl-tyramide reaction, thus multiplying the HRP molecules available at the surface for the substrate reaction.

# Reagents provided

- 1 aluminum pouch with a Microwell Plate (12 strips of 8 wells each) coated with monoclonal antibody to human IL-10
- 1 vial (70 µL) Biotin-Conjugate anti-human IL-10 monoclonal antibody
- 1 vial (150 µL) Streptavidin-HRP
- 2 vials human IL-10 Standard lyophilized, 400 pg/mL upon reconstitution
- 1 vial Control, lyophilized
- 1 vial (5 mL) Assay Buffer Concentrate 20x (PBS with 1% Tween<sup>™</sup> 20 and 10% BSA)
- 2 bottles (12 mL) Sample Diluent
- 1 vial (7 mL) Amplification Diluent Concentrate (2x)
- 2 vials (75 µL) Amplification Reagent I

Note: reagent contains ethyl alcohol

- 1 vial (15 µL) Amplification Reagent II
- 2 bottles (50 mL) Wash Buffer Concentrate 20x (PBS with 1% Tween<sup>™</sup> 20)
- 1 vial (15 mL) Substrate Solution (tetramethyl-benzidine)
- 1 vial (15 mL) Stop Solution (1M Phosphoric acid)
- 8 Adhesive Films

### Storage instructions – ELISA kit

Store kit reagents between  $2^{\circ}$  and  $8^{\circ}$ C except controls. Store lyophilized controls at -20°C.

Immediately after use remaining reagents should be returned to cold storage (2° to 8°C), controls to -20°C, respectively. Expiry of the kit and reagents is stated on labels.

Expiry of the kit components can only be guaranteed if the components are stored properly, and if, in case of repeated use of one component, this reagent is not contaminated by the first handling.

# Sample collection and storage instructions

Cell culture supernatant, serum, and plasma (citrate and heparin) were tested with this assay. Other biological samples might be suitable for use in the assay. Remove serum or plasma from the clot or cells as soon as possible after clotting and separation.

Samples containing a visible precipitate must be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic samples.

Samples should be aliquoted and must be stored frozen at  $-20^{\circ}$ C to avoid loss of bioactive human IL-10. If samples are to be run within 24 hours, they may be stored at 2–8°C (for sample stability refer to "Sample stability" on page 7).

Avoid repeated freeze-thaw cycles. Prior to assay, the frozen sample should be brought to room temperature slowly and mixed gently.

### Materials required but not provided

- 5 mL and 10 mL graduated pipettes
- 5 µL to 1000 µL adjustable single channel micropipettes with disposable tips
- 50 μL to 300 μL adjustable multichannel micropipette with disposable tips
- · Multichannel micropipette reservoir
- Beakers, flasks, cylinders necessary for preparation of reagents
- Device for delivery of wash solution (multichannel wash bottle or automatic wash system)
- · Microplate shaker
- Microwell strip reader capable of reading at 450 nm (620 nm as optional reference wave length)
- Glass-distilled or deionized water
- Statistical calculator with program to perform regression analysis

### Precautions for use

- All chemicals should be considered as potentially hazardous. We
  therefore recommend that this product is handled only by those
  persons who have been trained in laboratory techniques and that
  it is used in accordance with the principles of good laboratory
  practice. Wear suitable protective clothing such as laboratory
  overalls, safety glasses, and gloves. Care should be taken to
  avoid contact with skin or eyes. In the case of contact with skin
  or eyes wash immediately with water. See material safety data
  sheet(s) and/or safety statement(s) for specific advice.
- Reagents are intended for research use only and are not for use in diagnostic or therapeutic procedures.
- Do not mix or substitute reagents with those from other lots or other sources.
- Do not use kit reagents beyond expiration date on label.
- Do not expose kit reagents to strong light during storage or incubation.
- Do not pipet by mouth.
- Do not eat or smoke in areas where kit reagents or samples are handled.
- Avoid contact of skin or mucous membranes with kit reagents or samples.
- Rubber or disposable latex gloves should be worn while handling kit reagents or samples.
- Avoid contact of substrate solution with oxidizing agents and metal.
- Avoid splashing or generation of aerosols.
- To avoid microbial contamination or cross-contamination of reagents or samples that may invalidate the test, use disposable pipette tips and/or pipettes.
- Use clean, dedicated reagent trays for dispensing the conjugate and substrate reagent.
- Exposure to acid inactivates the conjugate.
- Glass-distilled water or deionized water must be used for reagent preparation.
- Substrate solution must be at room temperature prior to use.
- Decontaminate and dispose samples and all potentially contaminated materials as if they could contain infectious agents.
   The preferred method of decontamination is autoclaving for a minimum of 1 hour at 121.5°C.
- Liquid wastes not containing acid and neutralized waste may be mixed with sodium hypochlorite in volumes such that the final mixture contains 1.0% sodium hypochlorite. Allow 30 minutes for effective decontamination. Liquid waste containing acid must be neutralized prior to the addition of sodium hypochlorite.

# Preparation of reagents

- 1. Buffer concentrates should be brought to room temperature and should be diluted before starting the test procedure.
- 2. If crystals have formed in the Buffer Concentrates, warm them gently until they have completely dissolved.

#### Wash buffer (1x)

- Pour entire contents (50 mL) of the Wash Buffer Concentrate (20x) into a clean 1000 mL graduated cylinder. Bring to final volume of 1000 mL with glass-distilled or deionized water. Mix gently to avoid foaming.
- 2. Transfer to a clean wash bottle and store at 2° to 25°C. Please note that Wash Buffer (1x) is stable for 30 days.
- Wash Buffer (1x) may also be prepared as needed according to the following table:

Number of Strips	Wash Buffer Concentrate (20x) (mL)	Distilled Water (mL)
1 - 6	25	475
1 - 12	50	950

### Assay buffer (1x)

- Pour the entire contents (5 mL) of the Assay Buffer Concentrate (20x) into a clean 100 mL graduated cylinder. Bring to final volume of 100 mL with distilled water. Mix gently to avoid foaming.
- 2. Store at 2–8°C. The Assay Buffer (1x) is stable for 30 days.
- Assay Buffer (1x) may also be prepared as needed according to the following table:

Number of Strips	Assay Buffer Concentrate (20x) (mL)	Distilled Water (mL)
1–6	2.5	47.5
1–12	5.0	95.0

### Biotin-Conjugate

Note: The Biotin-Conjugate should be used within 30 minutes after dilution

Make a 1:100 dilution of the concentrated Biotin-Conjugate solution with Sample Diluent in a clean plastic tube as needed according to the following table:

Number of Strips	Biotin-Conjugate (mL)	Sample Diluent (mL)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

#### Streptavidin-HRP

Note: The Streptavidin-HRP should be used within 30 minutes after dilution.

Make a 1:200 dilution of the concentrated Streptavidin-HRP solution with Assay Buffer (1x) in a clean plastic tube as needed according to the following table:

Number of Strips	Streptavidin-HRP (mL)	Assay Buffer (1x) (mL)
1 - 6	0.03	5.97
1 - 12	0.06	11.94

### Human IL-10 standard

- Reconstitute human IL-10 standard by addition of distilled water. Reconstitution volume is stated on the label of the standard vial. Swirl or mix gently to insure complete and homogeneous solubilization (concentration of reconstituted standard = 400 pg/mL).
- Allow the standard to reconstitute for 10-30 minutes. Mix well prior to making dilutions.
- After usage remaining standard cannot be stored and has to be discarded.
- 4. The concentrated human IL-10 standard must be diluted 1:8 with Sample Diluent just prior to use in a clean plastic test tube according to the following dilution scheme:
  - 100  $\mu$ L concentrated human IL-10 standard + 700  $\mu$ L Sample Diluent. Shake gently to mix (concentration of standard = 50 pg/mL).
- Standard dilutions can be prepared directly on the microwell plate (see "Test protocol" on page 4) or alternatively in tubes (see "External standard dilution" on page 3).

#### External standard dilution

- Label 7 tubes, one for each standard point: S1, S2, S3, S4, S5, S6, S7
- 2. Prepare 1:2 serial dilutions for the standard curve as follows: Pipette 225  $\mu L$  of Sample Diluent into each tube.
- Pipette 225 μL of diluted standard (concentration of standard = 50 pg/mL) into the first tube, labeled S1, and mix (concentration of standard 1 = 25 pg/mL).
- Pipette 225 μL of this dilution into the second tube, labeled S2, and mix thoroughly before the next transfer.

Repeat serial dilutions 5 more times thus creating the points of the standard curve (see Figure 8).

Sample Diluent serves as blank.

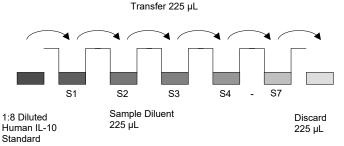


Fig. 8 Dilute standards - tubes

#### Control

- Reconstitute lyophilized controls by adding distilled water (10-30 minutes). Reconstitution volume is stated on the label of the control vial. Swirl or mix gently to ensure complete and homogeneous solubilization. Predilute solubilized control 1:9 in Sample Diluent: 100 μL control + 800 μL Sample Diluent.
- Further treat the controls like your samples in the assay. For control range please refer to certificate of analysis or vial label. Store reconstituted control aliquoted at -20°C. Avoid repeated freeze and thaw cycles.

### Amplification diluent (1x)

Preparation of Amplification Diluent (1x) has to be done immediately prior to use.

Make a 1:2 dilution of the concentrated Amplification Diluent (2x) as needed according to the following table:

Number of Strips	Amplification Diluent (2x) (mL)	Distilled Water (mL)
1–6	3	3
1–12	6	6

### Amplification Solution I

- 1. Prepare Amplification Solution I immediately prior to application on the plate.
- 2. Dilute Amplification Reagent I in Amplification Diluent (1x) as indicated in the Certificate of Analysis.
- Discard immediately any prediluted Amplification Solution I after usage.

#### Amplification solution II

- Prepare Amplification Solution II immediately prior to application on the plate.
- Centrifuge vial for a few seconds in a microcentrifuge before opening to collect liquid trapped in the lid.
- 3. Dilute Amplification Reagent II in Assay Buffer (1x) as as indicated in the Certificate of Analysis.
- 4. Discard immediately any prediluted Amplification Solution II after usage.

### Test protocol

### IMPORTANT!

- Because this ELISA is a high sensitivity system, it is extremely important to stick exactly to the manual (washing procedure, chronology of and preparation of solutions, incubation time) to obtain optimal test performance.
- Amplification Solutions must be prepared immediately prior to application on the plate! It is extremely important to wash the wells properly to obtain a good test performance.
- Shaking is absolutely necessary for an optimal test performance.
   Protect microwell plate from light during incubation steps

- Determine the number of microwell strips required to test
  the desired number of samples plus appropriate number of
  wells needed for running blanks and standards. Each sample,
  standard, blank and optional control sample should be assayed
  in duplicate. Remove extra microwell strips from holder and store
  in foil bag with the desiccant provided at 2°-8°C sealed tightly.
- Wash the microwell strips twice with exactly 400 μL Wash Buffer per well with thorough aspiration of microwell contents between washes. Allow the Wash Buffer to sit in the wells for about 10 15 seconds before aspiration. Soaking is highly recommended between the washes to obtain a good test performance! Take care not to scratch the surface of the microwells.

After the last wash step, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess Wash Buffer. Use the microwell strips immediately after washing. Do not allow wells to dry.

3. Standard dilution on the microwell plate (Alternatively the standard dilution can be prepared in tubes - see "External standard dilution" on page 3.):

Add 100  $\mu$ L of Sample Diluent in duplicate to all standard wells. Pipette 100  $\mu$ L of prepared standard (see Preparation of Standard "Human IL-10 standard" on page 3, concentration = 50.00 pg/mL) in duplicate into well A1 and A2 (see Table 1). Mix the contents of wells A1 and A2 by repeated aspiration and ejection (concentration of standard 1, S1 = 25.00 pg/mL), and transfer 100  $\mu$ L to wells B1 and B2, respectively. (see Figure 9). Take care not to scratch the inner surface of the microwells. Continue this procedure 5 times, creating two rows of human IL-10 standard dilutions ranging from 25.00 to 0.39 pg/mL. Discard 100  $\mu$ L of the contents from the last microwells (G1, G2) used.

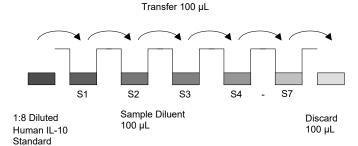


Fig. 9 Dilute standards - microwell plate

Table 1 Example of the arrangement of blanks, standards and samples in the microwell strips.

	1	2	3	4
А	Standard 1 25.00 pg/mL	Standard 1 25.00 pg/mL	Sample 1	Sample 1
В	Standard 2 12.50 pg/mL	Standard 2 12.50 pg/mL	Sample 2	Sample 2
С	Standard 3 6.25 pg/mL	Standard 3 6.25 pg/mL	Sample 3	Sample 3
D	Standard 4 3.13 pg/mL	Standard 4 3.13 pg/mL	Sample 4	Sample 4
Е	Standard 5 1.56 pg/mL	Standard 5 1.56 pg/mL	Sample 5	Sample 5
F	Standard 6 0.78 pg/mL	Standard 6 0.78 pg/mL	Sample 6	Sample 6
G	Standard 7 0.39 pg/mL	Standard 7 0.39 pg/mL	Sample 7	Sample 7
Н	Blank	Blank	Sample 8	Sample 8

In case of an external standard dilution (see "External standard dilution" on page 3.), pipette  $100~\mu L$  of these standard dilutions (S1 - S7) in the standard wells according to Table 1.

- 4. Add 100 µL of Sample Diluent in duplicate to the blank wells.
- 5. Add 50 µL of Sample Diluent to the sample wells.
- 6. Add 50  $\mu$ L of each sample in duplicate to the sample wells.

- Prepare Biotin-Conjugate (see Preparation of Biotin-Conjugate "Biotin-Conjugate" on page 3).
- 8. Add 50  $\mu$ L of Biotin-Conjugate to all wells.
- Cover with an adhesive film and incubate at room temperature (18° to 25°C) for 2 hours, if available on a microplate shaker (Shaking is absolutely necessary for an optimal test performance).
- **10.** Prepare Streptavidin-HRP (refer to Preparation of Streptavidin-HRP "Streptavidin-HRP" on page 3).
- 11. Remove adhesive film and empty wells. Wash microwell strips 6 times according to point 2. of the test protocol. Proceed immediately to the next step.
- 12. Add 100  $\mu L$  of diluted Streptavidin-HRP to all wells, including the blank wells.
- 13. Cover with an adhesive film and incubate at room temperature (18° to 25°C) for 1 hour on a microplate shaker (Shaking is absolutely necessary for an optimal test performance).
- Prepare Amplification Solution I diluted in Amplification Diluent (1x) (see "Amplification Solution I" on page 4) immediately prior to use.
- 15. Remove adhesive film and empty wells. Wash microwell strips 6 times according to point 2. of the test protocol. Proceed immediately to the next step.
- 16. Add 100  $\mu L$  of Amplification Solution I to all wells, including the blank wells.
- 17. Cover with an adhesive film and incubate at room temperature (18° to 25°C) for exactly 15 minutes on a microplate shaker. (Shaking is absolutely necessary for an optimal test performance.)
- **18.** Prepare Amplification Solution II diluted in Assay Buffer (1x) (see "Amplification solution II" on page 4) immediately prior to use.
- 19. Remove adhesive film and empty wells. Wash microwell strips 6 times according to point 2. of the test protocol. Proceed immediately to the next step.
- 20. Add 100  $\mu L$  of Amplification Solution II to all wells, including the blank wells.
- Cover with an adhesive film and incubate at room temperature (18° to 25°C) for exactly 30 minutes, if available on a microplate shaker.
- 22. Remove adhesive film and empty wells. Wash microwell strips 6 times according to point 2. of the test protocol. Proceed immediately to the next step.
- 23. Pipette 100  $\mu$ L of TMB Substrate Solution to all wells.
- 24. Incubate the microwell strips at room temperature (18° to 25°C) for about 10-20 minutes. Avoid direct exposure to intense light.

The color development on the plate should be monitored and the substrate reaction stopped (see next point of this protocol) before positive wells are no longer properly recordable. Determination of the ideal time period for color development has to be done individually for each assay.

It is recommended to add the stop solution when the highest standard has developed a dark blue color. Alternatively the color development can be monitored by the ELISA reader at 620 nm. The substrate reaction should be stopped as soon as Standard 1 has reached an OD of 0.9 – 0.95.

- 25. Stop the enzyme reaction by quickly pipetting 100 μL of Stop Solution into each well. It is important that the Stop Solution is spread quickly and uniformly throughout the microwells to completely inactivate the enzyme. Results must be read immediately after the Stop Solution is added or within one hour if the microwell strips are stored at 2 8°C in the dark.
- 26. Read absorbance of each microwell on a spectro-photometer using 450 nm as the primary wave length (optionally 620 nm as the reference wave length; 610 nm to 650 nm is acceptable). Blank the plate reader according to the manufacturer's instructions by using the blank wells. Determine the absorbance of both the samples and the standards.

Note: In case of incubation without shaking the obtained O.D. values may be lower than indicated below. Nevertheless the results are still valid

### Calculation of results

- Calculate the average absorbance values for each set of duplicate standards and samples. Duplicates should be within 20 percent of the mean value.
- Create a standard curve by plotting the mean absorbance for each standard concentration on the ordinate against the human IL-10 concentration on the abscissa. Draw a best fit curve through the points of the graph (a 5-parameter curve fit is recommended).
- To determine the concentration of circulating human IL-10 for each sample, first find the mean absorbance value on the ordinate and extend a horizontal line to the standard curve. At the point of intersection, extend a vertical line to the abscissa and read the corresponding human IL-10 concentration.
- If instructions in this protocol have been followed samples have been diluted 1:2 (50 μL sample + 50 μL Sample Diluent) and controls 1:18 (50 μL of 1:9 prediluted control + 50 μL Sample Diluent). Thus concentrations read from the standard curve must be multiplied by the dilution factor (x 2 for samples, x 18 for controls).
- Calculation of samples with a concentration exceeding standard 1 may result in incorrect, low human IL-10 levels. Such samples require further external predilution according to expected human IL-10 values with Sample Diluent in order to precisely quantitate the actual human IL-10 level.
- It is suggested that each testing facility establishes a control sample of known human IL-10 concentration and runs this additional control with each assay. If the values obtained are not within the expected range of the control, the assay results may be invalid.
- A representative standard curve is shown in Figure 10.

**Note:** Do not use this standard curve to derive test results. Each laboratory must prepare a standard curve for each group of microwell strips assayed.

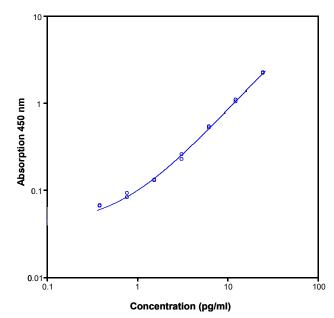


Fig. 10 Representative standard curve for human IL-10 ELISA. Human IL-10 was diluted in serial 2-fold steps in Sample Diluent.

Table 2 Typical data using the human IL-10 ELISA.

Measuring wavelength: 450 nm Reference wavelength: 620 nm

Standard	human IL-10 Concentration (pg/mL)	O.D. at 450 nm	Mean O.D. at 450 nm	C.V. (%)
1	25.00	2.193 2.201	2.197	0.3
2	12.50	1.020 1.070	1.045	3.4
3	6.25	0.531 0.514	0.523	2.3
4	3.13	0.253 0.225	0.239	8.3
5	1.56	0.129 0.127	0.128	1.1
6	0.78	0.091 0.081	0.086	8.2
7	0.39	0.065 0.066	0.066	1.1
Blank	0	0.037 0.036	0.037	1.4

The OD values of the standard curve may vary according to the conditions of assay performance (e.g., operator, pipetting technique, washing technique, or temperature effects). Furthermore, shelf life of the kit may affect enzymatic activity and thus color intensity. Values measured are still valid.

### Limitations

- Since exact conditions may vary from assay to assay, a standard curve must be established for every run.
- Bacterial or fungal contamination of either screen samples or reagents or cross-contamination between reagents may cause erroneous results.
- Disposable pipette tips, flasks or glassware are preferred, reusable glassware must be washed and thoroughly rinsed of all detergents before use.
- Improper or insufficient washing at any stage of the procedure will
  result in either false positive or false negative results. Empty wells
  completely before dispensing fresh wash solution, fill with Wash
  Buffer as indicated for each wash cycle and do not allow wells to
  sit uncovered or dry for extended periods.
- The use of radioimmunotherapy has significantly increased the number of subjects with human anti-mouse IgG antibodies (HAMA). HAMA may interfere with assays utilizing murine monoclonal antibodies leading to both false positive and false negative results. Serum samples containing antibodies to murine immunoglobulins can still be analyzed in such assays when murine immunoglobulins (serum, ascitic fluid, or monoclonal antibodies of irrelevant specificity) are added to the sample.

# Performance characteristics

### Sensitivity

The limit of detection of human IL-10 defined as the analyte concentration resulting in an absorbance significantly higher than that of the dilution medium (mean plus 2 standard deviations) was determined to be 0.05 pg/mL (mean of 6 independent assays).

### Reproducibility

Intra-assay

Reproducibility within the assay was evaluated in 3 independent experiments. Each assay was carried out with 6 replicates of 8 serum samples containing different concentrations of human IL-10. 2 standard curves were run on each plate. Data below show the mean human IL-10 concentration and the coefficient of variation for each

sample (see Table 3). The calculated overall intra-assay coefficient of variation was 6.8%.

Table 3 The mean human IL-10 concentration and the coefficient of variation for each sample

Sample	Experiment	Mean Human IL-10 Concentration (pg/mL)	Coefficient of Variation (%)
	1	7.64	10
1	2	7.99	8
	3	7.97	8
	1	1.27	8
2	2	1.16	8
	3	1.47	8
	1	1.16	7
3	2	1.12	9
	3	1.09	5
	1	5.11	1
4	2	5.23	3
	3	6.39	4
	1	1.55	7
5	2	1.38	8
	3	1.64	8
	1	3.65	4
6	2	3.27	8
	3	4.41	6
	1	2.45	8
7	2	2.68	7
	3	3.04	7
	1	3.09	9
8	2	3.13	7
	3	3.60	6

#### Inter-assay

Assay to assay reproducibility within one laboratory was evaluated in 3 independent experiments. Each assay was carried out with 6 replicates of 8 serum samples containing different concentrations of human IL-10. 2 standard curves were run on each plate. Data below show the mean human IL-10 concentration and the coefficient of variation calculated on 18 determinations of each sample (see Table 4). The calculated overall inter-assay coefficient of variation was 7.5%.

Table 4 The mean human IL-10 concentration and the coefficient of variation of each sample

Sample	Mean Human IL-10 Concentration (pg/mL)	Coefficient of Variation (%)
1	7.87	9.0
2	1.30	8.0
3	1.12	7.1
4	5.58	2.5
5	1.53	7.7
6	3.78	6.1
7	2.73	7.1
8	3.27	7.1

#### Spike recovery

The spike recovery was evaluated by spiking 4 levels of human IL-10 into pooled normal human serum samples. Recoveries were determined in 2 independent experiments with 8 replicates each. The amount of endogenous human IL-10 in unspiked serum was subtracted from the two spike values. The recovery ranged from 78–111% with an overall mean recovery of 94%.

#### Dilution parallelism

Four serum samples with different levels of human IL-10 were analyzed at serial 2-fold dilutions with 4 replicates each. The recovery ranged from 98% to 118% with an overall recovery of 111%.

Sample	Dilution	Expected human IL-10 concentratio n (pg/mL)	Observed human IL-10 concentratio n (pg/mL)	Recovery of expected human IL-10 concentratio n (%)
	1	_	40.4	_
1	2	20.2	22.3	111
'	4	11.2	13.2	118
	8	6.6	7.4	112
	1	-	38.2	-
2	2	19.1	18.8	98
2	4	9.4	11.1	118
	8	5.5	6.4	115
	1	_	17.7	_
3	2	8.9	9.3	105
3	4	4.7	5.5	117
	8	2.7	3.0	109
	1	_	40.4	_
4	2	20.2	22.3	111
4	4	11.2	13.2	118
	8	6.6	7.4	112

### Sample stability

### Freeze-Thaw stability

Aliquots of serum samples (spiked or unspiked) were stored at -20°C and thawed 5 times, and the human IL-10 levels determined. There was no significant loss of human IL-10 immunoreactivity detected by freezing and thawing.

#### Storage stability

Aliquots of serum samples (spiked or unspiked) were stored at –20°C, 2–8°C, room temperature, and at 37°C, and the human IL-10 level determined after 24 hours. There was no significant loss of human IL-10 immunoreactivity detected during storage at –20°C, 2–8°C and room temperature.

A significant loss of human IL-10 immunoreactivity (30%) was detected during storage at 37°C after 24 hours.

# Comparison of serum and plasma

From 2 individuals each, serum, citrate and heparin plasma was obtained at the same time and tested for human IL-10. Concentrations were not significantly different and therefore all these blood preparations are suitable for use in the assay. It is nevertheless highly recommended to assure the uniformity of blood preparations.

#### Specificity

The interference of circulating factors of the immune system was evaluated by spiking these proteins at physiologically relevant concentrations into a human IL-10 positive serum. No cross-reactivity was detected.

#### **Expected values**

A panel of 16 serum samples from randomly selected apparently healthy donors (males and females) was tested for human IL-10. The detected human IL-10 sera levels ranged between 0.16 and 12.70 pg/mL with a mean level of 4.70 pg/mL and a standard deviation of 3.40 pg/mL. The normal levels measured may vary with the sample collective used.

#### Calibration

This immunoassay is calibrated with highly purified recombinant human IL-10, which has been evaluated against the International Reference Standard NIBSC 93/722 and has been shown to be equivalent.

NIBSC 93/722 is quantitated in International Units (IU), 1IU corresponding to 200 pg human IL-10.

# Reagent preparation summary

### Wash buffer (1x)

Add Wash Buffer Concentrate 20x (50 mL) to 950 mL distilled water.

Number of Strips	Wash Buffer Concentrate (mL)  Distilled Wat	
1 - 6	25	475
1 - 12	50	950

#### Assay buffer (1x)

Add Assay Buffer Concentrate 20x (5 mL) to 95 mL distilled water.

Number of Strips	Assay Buffer Concentrate (mL)	Distilled Water (mL)
1–6	2.5	47.5
1–12	5.0	95.0

### Biotin-Conjugate

Make a 1:100 dilution of Biotin-Conjugate in Sample Diluent:

Number of Strips	Biotin-Conjugate (mL)	Sample Diluent (mL)
1 - 6	0.03	2.97
1 - 12	0.06	5.94

### Streptavidin-HRP

Make a 1:200 dilution of Streptavidin-HRP in Assay Buffer (1x):

Number of Strips	Streptavidin-HRP (mL)	Assay Buffer (1x) (mL)
1 - 6	0.03	5.97
1 - 12	0.06	11.94

#### Human IL-10 standard

Reconstitute lyophilized human IL-10 standard with distilled water. (Reconstitution volume is stated on the label of the standard vial.) The concentrated human IL-10 standard must be diluted 1:8 with Sample Diluent.

### Amplification diluent (1x)

Prepare Amplification Diluent (1x) immediately prior to use.

Number of Strips	Amplification Diluent (2x) (mL)	Distilled Water (mL)
1–6	3	3
1–12	6	6

# Amplification Solution I

Dilute Amplification Reagent I in Amplification Diluent (1x) immediately prior to application on the plate as indicated in the Certificate of Analysis.

### Amplification solution II

Centrifuge vial for a few seconds in a micro-centrifuge before opening to collect liquid trapped in the lid. Dilute Amplification Solution II in Assay Buffer (1x) immediately prior to application on the plate as indicated in the Certificate of Analysis.

#### Contro

Reconstitute lyophilized controls by adding distilled water (10-30 minutes). Reconstitution volume is stated on the label of the control vial. Predilute reconstituted control 1:9 with Sample Diluent.

### Test protocol summary

**Note:** Prepare Amplification Solutions immediately prior to application on the plate. It is extremely important to wash the wells properly to obtain a good test performance.

- 1. Determine the number of microwell strips required.
- 2. Wash microwell strips twice with Wash Buffer.
- Standard dilution on the microwell plate: Add 100 μL Sample Diluent, in duplicate, to all standard wells. Pipette 100 μL prepared standard into the first wells and create standard dilutions by transferring 100 μL from well to well. Discard 100 μL from the last wells.

Alternatively external standard dilution in tubes (see "External standard dilution" on page 3): Pipette 100  $\mu$ L of these standard dilutions in the microwell strips.

- 4. Add 100 µL Sample Diluent in duplicate, to the blank wells.
- 5. Add 50 µL Sample Diluent to sample wells.
- 6. Add 50 µL sample in duplicate, to designated sample wells.
- 7. Prepare Biotin-Conjugate.
- 8. Add 50 µL Biotin-Conjugate to all wells.
- 9. Cover microwell strips and incubate 2 hours at room temperature (18° to 25°C). (Shaking is absolutely necessary for an optimal test performance.)
- 10. Prepare Streptavidin-HRP.
- 11. Empty and wash microwell strips 6 times with Wash Buffer.
- 12. Add 100 µL diluted Streptavidin-HRP to all wells.
- 13. Cover microwell strips and incubate 1 hour at room temperature (18° to 25°C). (Shaking is absolutely necessary for an optimal test performance.)
- 14. Prepare Amplification Solution I diluted in Amplification Diluent (1x) immediately prior to application on the plate.
- 15. Empty and wash microwell strips 6 times with Wash Buffer.
- 16. Add 100 µL Amplification Solution I to all wells.
- 17. Cover microwell strips and incubate for exactly 15 minutes at room temperature (18° to 25°C). (Shaking is absolutely necessary for an optimal test performance.)
- **18.** Prepare Amplification Solution II diluted in Assay Buffer (1x) immediately prior to application on the plate.
- 19. Empty and wash microwell strips 6 times with Wash Buffer

- 20. Add 100 µL Amplification Solution II to all wells.
- 21. Cover microwell strips and incubate for exactly 30 minutes at room temperature (18° to 25°C).
- 22. Empty and wash microwell strips 6 times with Wash Buffer.
- 23. Add 100 µL of TMB Substrate Solution to all wells.
- 24. Incubate the microwell strips for about 10-20 minutes at room temperature (18°to 25°C).
- 25. Add 100 µL Stop Solution to all wells.
- 26. Blank microwell reader and measure color intensity at 450 nm.

**Note:** If instructions in this protocol have been followed samples have been diluted 1:2 (50  $\mu$ L sample + 50  $\mu$ L Sample Diluent) and controls 1:18 (50  $\mu$ L of 1:9 prediluted control + 50  $\mu$ L Sample Diluent). Thus concentrations read from the standard curve must be multiplied by the dilution factor (x 2 for samples, x 18 for controls).

### Customer and technical support

Visit thermofisher.com/support for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
  - Product FAQs
  - Software, patches, and updates
  - Training for many applications and instruments
- Order and web support
- Product documentation
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at <a href="https://www.thermofisher.com/us/en/home/global/terms-and-conditions.html">www.thermofisher.com/us/en/home/global/terms-and-conditions.html</a>. If you have any questions, please contact Life Technologies at <a href="https://www.thermofisher.com/support">www.thermofisher.com/support</a>.



Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history: Pub. No. MAN0016564

Revision	Date	Description
D.0 (33)	29 April 2022	EDTA was removed from the document (see "Comparison of serum and plasma" on page 7).
		The control preparation instructions were updated (see "Control" on page 4).
		Initial release with revision history table.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2022 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. TSA is a trademark of PerkinElmer Life Sciences.

