



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

IL-2 (Human) Colorimetric ELISA Detection Kit

Catalog #79774-2

Size: 5 x 96 reactions

DESCRIPTION: The cytokine protein Interleukin-2 is an important factor for the maintenance of CD4⁺ regulatory T cells. It plays a critical role in the differentiation of CD4⁺ T cells into a variety of subsets as well as promotes CD8⁺ T-cell and NK cell cytotoxicity. The *IL-2 (Human) Colorimetric ELISA Detection Kit* is designed for detecting and quantifying human interleukin-2 in cell culture medium. This kit comes in a convenient 96-well format, with capture and detection antibodies for IL-2, streptavidin-labeled HRP, blocking buffer, IL-2 standard, and colorimetric HRP substrate for a 96-well plate. Only a few simple steps on a microtiter plate are required for the assay. First, the capture antibody is coated on a 96-well plate. Next, samples containing IL-2 are incubated on the coated plate followed by detecting the captured IL-2 with the detection antibody. Finally, the plate is treated with streptavidin-HRP followed by addition of a colorimetric HRP substrate to produce color, which can then be measured using a UV/Vis spectrophotometer microplate reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
	Interleukin-2 capture antibody	5 x 20 µl	-80°C	(Avoid freeze/thaw cycles!)
	Interleukin-2 detection antibody, biotinylated	5 x 5 µl	-80°C	
	Human IL-2 standard (1 µg/ml)	40 µl	-80°C	
80611	Streptavidin-HRP	5 x 5 µl	+4°C	
79743	Blocking Buffer 3	5 x 50 ml	+4°C	
79651	Colorimetric HRP substrate	5 x 10 ml	+4°C	
	Transparent 96-well microplate	5 x 1	+4°C	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

PBS (Phosphate-buffered saline)

PBST (PBS with 0.05% Tween-20)

1N HCl (aqueous)

Rotating or rocker platform

UV/Vis spectrophotometer microplate reader capable of reading absorbance at 450 nm*

*Alternatively, a spectrophotometer reading at 650 nm may be used without adding 1N HCl, but sensitivity of the assay will be reduced.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

APPLICATIONS: This kit is useful for cytokine detection in cell culture medium.

STABILITY: One year from date of receipt when stored as directed.

REFERENCES:

1. Jiang, T., *et al.* *Role of IL-2 in Cancer Immunotherapy.* *Oncolmmunology*, 2016. **5(6):** e1163462

ASSAY PROTOCOL:

All samples and standards should be tested in duplicate.

Coating the plate with capture Ab:

- 1) Thaw **capture Ab** on ice. Upon first thaw, *briefly* spin tube containing **capture Ab** to recover the full contents of the tube.
- 2) Dilute **capture Ab** to 2 ng/μl in PBS.
- 3) Add 50 μl of diluted **capture Ab** solution to each well and incubate overnight at 4°C.
(Remaining **capture Ab** can be stored at 4°C)

**After overnight coating, it is highly recommended that all remaining steps are completed the following day to obtain the optimal sensitivity.*

- 4) After the overnight incubation, decant to remove the solution. Wash the plate 2 times with 200 μl/well of PBS with 0.05% Tween-20 (PBST). Tap plate onto clean paper towels to remove liquid.
- 5) Block wells by adding 200 μl of **Blocking Buffer** to each well. Incubate for 1 hour at room temperature. Decant to remove the blocking buffer and wash the plate 2 times with 200 μl/well of PBST. Tap plate onto clean paper towels to remove liquid.

Step 1:

- 1) Prepare the sample by diluting in the **Blocking Buffer**. Detection range of the *Colorimetric Human IL-2 Detection Kit* is 5 pg/ml – 500 pg/ml (**Figure 1**). Roughly estimate the amount of human IL-2 in the sample and dilute it accordingly. For quantification, **human IL-2 standard** can be serially diluted (1,000 pg/ml to 5 pg/ml) in the **Blocking Buffer** and run in the same plate. (Aliquot remaining **human IL-2 standard** into single use aliquots. Immediately store remaining undiluted protein in aliquots at -80°C.)

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 2) Add 50 μ l of the diluted samples (or **human IL-2 standard**) to each well and incubate the plate for 2 hours at room temperature.
- 3) After 2 hours incubation, decant to remove the solution and wash the plate 2 times with 200 μ l/well of PBST. Tap plate onto clean paper towels to remove liquid.
- 4) Dilute **biotinylated-detection Ab** 1:1,000 in the Blocking Buffer, and add 50 μ l to the wells. Incubate the plate for 1 hour at room temperature.
- 5) After 1-hour incubation, decant to remove the solution and wash the plate 3 times with 200 μ l/well of PBST. Tap plate onto clean paper towels to remove liquid.
- 6) Dilute **Streptavidin-HRP** 1:1,000 in the Blocking Buffer, and add 50 μ l to the wells. Incubate the plate for 30 minutes at room temperature.
- 7) After 30 minutes incubation, decant to remove the solution and wash the plate 5 times with 200 μ l/well of PBST. Tap plate onto clean paper towels to remove liquid
- 8) Add 100 μ l of the **Colorimetric HRP substrate** to each well and incubate the plate at room temperature until blue color is developed in the positive control well. This usually takes a few to several minutes to fully develop. The optimal incubation time may vary, and should be determined empirically by the user.
- 9) After the blue color is developed, add 100 μ l of 1 M HCl to each well. Read the absorbance at 450 nm using UV/Vis spectrophotometer microplate reader. The blank wells should exhibit an absorbance of \sim 0.05 at 450 nm. *Alternatively, the plate may be read at 650 nm without adding 1N HCl, but the Signal-to-Background ratio will be decreased.*

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

Example of Detection Results:

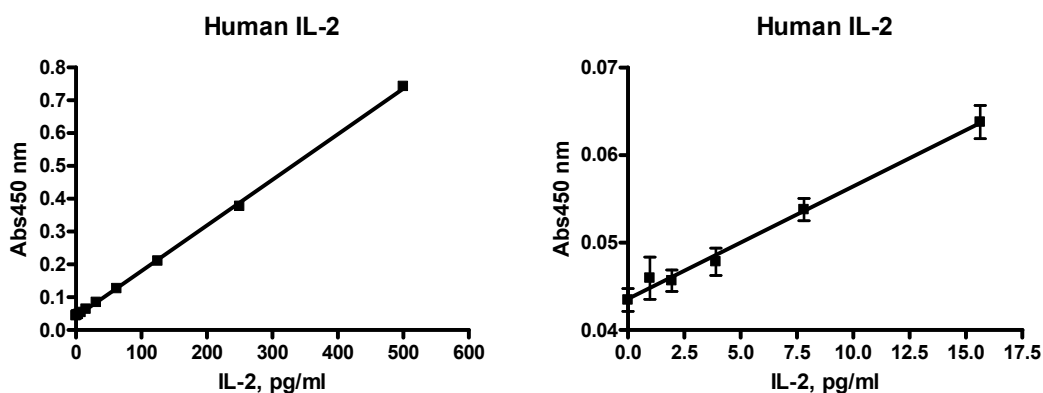


Figure 1. Detection of the human IL-2 (left: 1 pg/ml – 500 pg/ml, right: 1 pg/ml – 15 pg/ml) standard using the *Colorimetric Human IL-2 Detection Kit*. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com

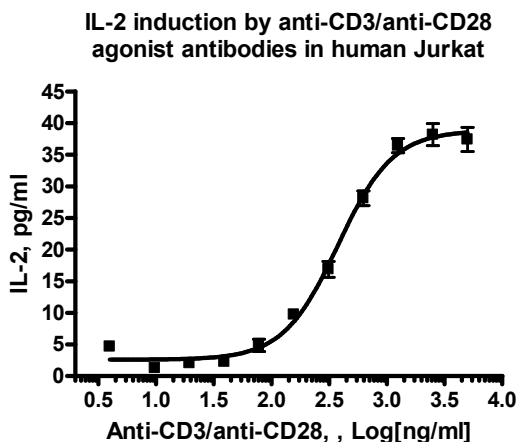


Figure 2. Induction of the human IL-2 in Jurkat cells by anti-CD3/anti-CD28 agonist antibodies. Jurkat cells were treated with the 1:1 mixture of the anti-CD3 (**BPS cat# 71274**) /anti-CD28 (**BPS cat# 100186**) agonist antibodies for 24 hours, and the amount of IL-2 secreted was measured by the *Colorimetric Human IL-2 Detection Kit*

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS:

Product Name

Catalog #

Size

Human Interleukin-1 beta
Human Interleukin-1 beta
Mouse Interleukin-1 beta
Mouse Interleukin-1 beta
Streptavidin-HRP
Blocking Buffer 3

90168-A
90168-B
90172-A
90172-B
80611
79743

2 µg
10 µg
2 µg
10 µg
100 µl
50 ml

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: [**info@bpsbioscience.com**](mailto:info@bpsbioscience.com)

Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)



6042 Cornerstone Court W, Ste B
 San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Colorimetric signal of positive control reaction is weak	Antibodies have lost activity	Antibodies and IL-1 β standard may lose activity upon repeated freeze/thaw cycles. Use fresh protein. Store proteins in single-use aliquots. Increase time of incubation. Increase protein or antibody concentration.
	Incorrect settings on instruments	Refer to instrument instructions for settings to increase sensitivity.
	Colorimetric HRP substrate was not incubated long enough	Increase the amount of time that the colorimetric HRP substrate is incubated in the wells. Avoid azides.
Colorimetric signal is erratic or varies widely among wells	Inaccurate pipetting/technique	Run duplicates of all reactions. Use a multichannel pipettor. Use master mixes to minimize errors.
	Bubbles in wells	Pipette slowly to avoid bubble formation. Tap plate lightly to disperse bubbles; be careful not to splash between wells.
	Signal is out of range of detection (too high)	Decrease the amount of time that the colorimetric HRP substrate is incubated in the wells
Background (signal to noise ratio) is high	Insufficient washes or blocking	Be sure to include blocking steps after wash steps. Increase number of washes. Increase wash volume. Increase Tween-20 concentration to 0.1% in TBST. Be sure to dilute Streptavidin-HRP in blocking buffer, not assay buffer.
	Sample solvent is inhibiting the protein	Run negative control assay including solvent. Maintain DMSO level at <1%. Increase time of protein incubation.
	Results are outside the linear range of the assay	Use different concentrations of IL-1 β standard to create a standard curve.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com