



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-1 β (Human) Colorimetric ELISA Detection Kit

Catalog #79771-2

Size: 5 x 96 reactions

DESCRIPTION: The cytokine protein Interleukin-1 β is a mediator in the inflammatory response and is a therapeutic target for autoinflammatory diseases such as Alzheimer's disease and Schnitzler syndrome. The *IL-1 β (Human) Colorimetric ELISA Detection Kit* is designed for detecting and quantifying human interleukin-1 β in cell culture medium. This kit comes in a convenient 96-well format, with capture and detection antibodies for IL-1 β , streptavidin-labeled HRP, blocking buffer, IL-1 β 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-1 β are incubated on the coated plate followed by detecting the captured IL-1 β 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-1 β capture antibody	5 x 20 μ l	-80 $^{\circ}$ C	(Avoid freeze/thaw cycles!)
	Interleukin-1 β detection antibody, biotinylated	5 x 5 μ l	-80 $^{\circ}$ C	
	Human IL-1 β standard (1 μ g/ml)	40 μ l	-80 $^{\circ}$ C	
80611	Streptavidin-HRP	5 x 5 μ l	+4 $^{\circ}$ C	
79743	Blocking Buffer 3	5 x 50 ml	+4 $^{\circ}$ C	
79651	Colorimetric HRP substrate	5 x 10 ml	+4 $^{\circ}$ C	
79964	Transparent 96-well microplate	5 x 1	+4 $^{\circ}$ 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. Sciacca, FL., *et al.* *Interleukin-1B polymorphism is associated with age at onset of Alzheimer's disease.* Neurobiology of Aging 2003. **24:** 927-931.
2. Roerink, M., *et al.* *Interleukin-1 as a mediator of fatigue in disease.* Journal of Neuroinflammation. 2017, **14:**16

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/ μ l in PBS.
- 3) Add 50 μ l 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 μ l/well of PBS with 0.05% Tween-20 (PBST). Tap plate onto clean paper towels to remove liquid.
- 5) Block wells by adding 200 μ l 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 μ l/well of PBST. Tap plate onto clean paper towels to remove liquid.

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

Step 1:

- 1) Prepare the sample by diluting in the **Blocking Buffer**. Detection range of the *Colorimetric Human IL-1 β Detection Kit* is 2 pg/ml – 500 pg/ml (**Figure 1**). Roughly estimate the amount of human IL-1 β in the sample and dilute it accordingly. For quantification, **human IL-1 β standard** can be serially diluted (500 pg/ml to 2 pg/ml) in the **Blocking Buffer** and run in the same plate. (Aliquot remaining **human IL-1 β** into single use aliquots. Immediately store remaining undiluted protein in aliquots at -80°C.)
- 2) Add 50 μ l of the diluted samples (or **human IL-1 β 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. 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 ~ 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:

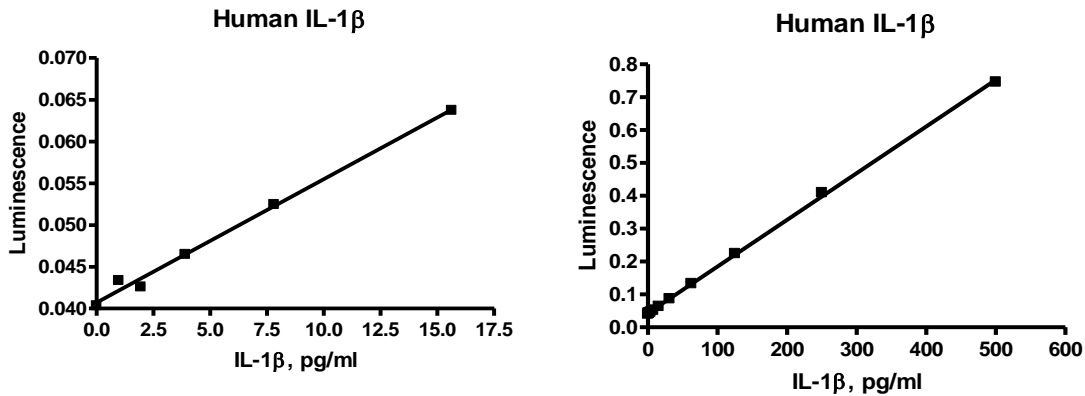


Figure 1. Detection of the human IL-1β (1 pg/ml – 15 pg/ml (Left), 1 pg/ml – 500 pg/ml (right) standard using the *Colorimetric™ Human IL-1β Detection Kit*. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com

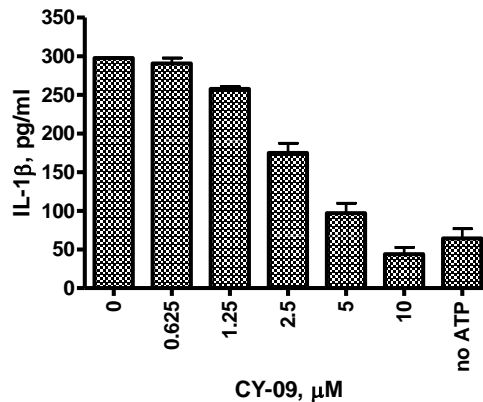


Figure 2. Detection of the human IL-1β from THP-1 cells after inflammasome activation. THP-1 cells were differentiated with PMA followed by priming and activation of the inflammasome with LPS and ATP respectively. Cell growing medium was removed and diluted 1:40 in Blocking Buffer, and the amount of IL-1β secreted by THP-1 cells was measured by the *Colorimetric Human IL-1β 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:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
Human Interleukin-1 beta	90168-A	2 µg
Human Interleukin-1 beta	90168-B	10 µg
Mouse Interleukin-1 beta	90172-A	2 µg
Mouse Interleukin-1 beta	90172-B	10 µg
Streptavidin-HRP	80611	100 µl
Blocking Buffer 3	79743	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