



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data Sheet

CTLA4[Biotinylated]:B7-2 Inhibitor Screening Assay Kit

Catalog # 72024

Size: 96 reactions

BACKGROUND: B7-2 (CD86) signaling through CTLA4(CD152) has been shown to inhibit T cell activation. This co-inhibitory pathway can be overactive in many tumors, enabling cancers to escape the host's immune system. CTLA4-blocking antibodies, including Ipilimumab (Yervoy) and Tremelimumab, have shown clinical efficacy in treating cancer.

DESCRIPTION: The *CTLA4[Biotinylated]:B7-2 Inhibitor Screening Assay Kit* is designed for screening and profiling inhibitors of CTLA4:B7-2 interaction. This kit comes in a convenient 96-well format, with biotin-labeled CTLA4 (CD152), purified B7-2 (CD86), streptavidin-labeled HRP, and assay buffer for 100 binding reactions. The key to this kit is the high sensitivity of detection of biotin-labeled CTLA4 by streptavidin-HRP. Only a few simple steps on a microtiter plate are required for the assay. First, B7-2 is coated on a 96-well plate. Next, CTLA4-biotin is incubated with B7-2 on the plate. Finally, the plate is treated with streptavidin-HRP followed by addition of an HRP substrate to produce chemiluminescence, which can then be measured using a chemiluminescence reader.

COMPONENTS:

| Catalog # | Component | Amount | Storage | |
|-----------|--|--------|---------|---|
| 71152 | CTLA4 (CD152), Fc fusion, Biotin-labeled | 3 µg | -80°C | (Avoid freeze/ thaw cycles!) |
| 71150 | B7-2 (CD86), Fc fusion | 10 µg | -80°C | |
| | Streptavidin-HRP | 15 µl | +4°C | |
| 79311 | 3x Immuno Buffer 1 | 50 ml | -20°C | |
| | Blocking Buffer | 50 ml | +4°C | |
| | HRP chemiluminescent substrate A (transparent bottle) | 6 ml | +4°C | |
| | HRP chemiluminescent substrate B (brown bottle) | 6 ml | +4°C | |
| | White 96-well microplate | 1 | +4°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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

PBS (Phosphate buffered saline)
Luminometer or microplate reader capable of reading chemiluminescence
Rotating or rocker platform

APPLICATIONS: This kit is useful for screening for inhibitors of CTLA4 binding to B7-2.

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

REFERENCES: 1. Ohtani, H., *et al.*, *Lab Invest.* 1997; **77(3)**: 231-241.
2. Rovert, C., *et al.*, *N. Engl. J. Med.* 2011; **364**: 2517-25262.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Coating the plate with B7-2:

- 1) Thaw **B7-2** on ice. Upon first thaw, briefly spin tube containing **B7-2** to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining B7-2 in aliquots at -80°C. *Note: B7-2 is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.*
- 2) Dilute **B7-2** to 2 µg/ml in PBS.
- 3) Add 50 µl of diluted **B7-2** solution to each well and incubate overnight at 4°C. Leave a couple of wells empty (uncoated), for use with the "Ligand Control" (see below).
- 4) Dilute **3x Immuno Buffer 1** to **1x Immuno Buffer 1** in water.
- 5) Decant to remove supernatant. Wash the plate 3 times with 100 µl **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 6) Block wells by adding 100 µl of **Blocking Buffer** to each well. Incubate for 1 hour at room temperature. Remove supernatant as described in step 4.

Step 1:

- 1) Prepare the master mixture: N wells × (10 µl **3x Immuno Buffer 1** + 15 µl H₂O).
- 2) Add 25 µl of master mixture to each well. Use uncoated wells for the "Ligand Control".

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

- 3) Add 5 µl of inhibitor solution to each well designated "Test Inhibitor". For the "Positive Control", "Ligand Control" and "Blank", add 5 µl of the same solution without inhibitor (inhibitor buffer). Incubate at room temperature for one hour.

| | Blank | Ligand Control | Positive Control | Test Inhibitor |
|---------------------------------|--------------|----------------|------------------|----------------|
| 3x Immuno Buffer 1 | 10 µl | 10 µl | 10 µl | 10 µl |
| H ₂ O | 15 µl | 15 µl | 15 µl | 15 µl |
| Test Inhibitor/Activator | – | – | – | 5 µl |
| Inhibitor buffer (no inhibitor) | 5 µl | 5 µl | 5 µl | – |
| 1x Immuno Buffer 1 | 20 µl | – | – | – |
| CTLA4-biotin (1 µg/ml) | – | 20 µl | 20 µl | 20 µl |
| Total | 50 µl | 50 µl | 50 µl | 50 µl |

- 4) Thaw **CTLA4-biotin** on ice. Upon first thaw, briefly spin tube containing protein to recover full contents of the tube. Aliquot **CTLA4-biotin** into single use aliquots. Immediately store remaining undiluted protein in aliquots at -80°C. *Note: CTLA4-biotin is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 5) Dilute **CTLA4-biotin** in **1x CTLA Assay Buffer** at 1 µg/ml. Keep diluted protein on ice until use. Discard any unused diluted protein after use.
- 6) Add 20 µl of **1x Immuno Buffer 1** to the well designated "Blank".
- 7) Initiate reaction by adding 20 µl of diluted **CTLA4-biotin** (see Step 1-5) to wells labeled "Positive Control", "Ligand Control" and "Test Inhibitor". Incubate at room temperature for two hours.
- 8) Decant to remove supernatant. Wash the plate 3 times with 100 µl/well **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 9) Block wells by adding 100 µl of **Blocking Buffer** to each well. Incubate for 10 minutes at room temperature. Remove supernatant as in Step 1-8.

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: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Step 2:

- 1) Dilute **Streptavidin-HRP** 1000-fold with **Blocking Buffer**.
- 2) Add 100 μ l to each well. Incubate for 1 hour at room temperature with slow shaking.
- 3) Wash plate three times with **1x Immuno Buffer 1**. Tap onto clean paper towels to remove liquid.
- 4) Block wells by adding 100 μ l of **Blocking Buffer** to each well. Incubate for 10 minutes at room temperature. Decant to remove supernatant. Tap plate onto clean paper towels to remove liquid.
- 5) Just before use, mix on ice 50 μ l **HRP Chemiluminescent Substrate A** and 50 μ l **HRP Chemiluminescent Substrate B** per well of the reaction, then add 100 μ l to each well. Discard any unused chemiluminescent reagent after use.
- 6) Immediately read sample in a luminometer or microtiter-plate capable of reading chemiluminescence. "Blank" value is subtracted from all readings.

Reading Chemiluminescence:

Chemiluminescence is the emission of light (luminescence) which results from a chemical reaction. The detection of chemiluminescence requires no wavelength selection because the method used is emission photometry and is not emission spectrophotometry.

To properly read chemiluminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second, delay after plate movement is 100 msec. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader are: use the "hole" position on the filter wheel; Optics position: Top; Read type: endpoint. Sensitivity may be adjusted based on the luminescence of a control assay without binding partner (typically we set this value as 100).

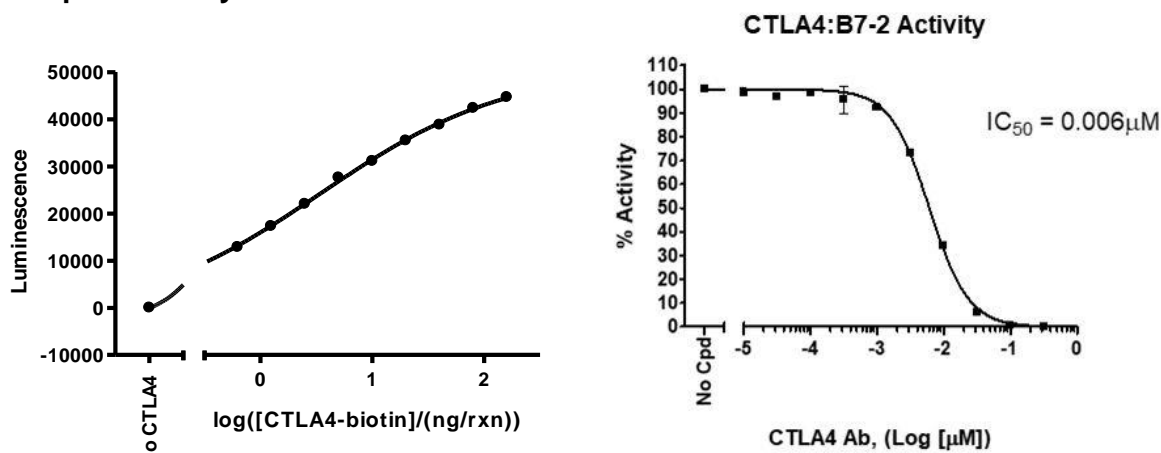
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
 San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Example of Assay Results:



CTLA4[Biotinylated]:B7-2 binding (left) and inhibition (right), measured using the using the *CTLA4[Biotinylated]:B7-2 Inhibitor Screening Assay Kit*, BPS Bioscience, Catalog #72024. Luminescence was measured using a Bio-Tek fluorescent microplate reader. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at info@bpsbioscience.com.*

RELATED PRODUCTS:

| <u>Product Name</u> | <u>Catalog #</u> | <u>Size</u> |
|--|------------------|-------------|
| CTLA4 (CD152), Fc fusion | 71149 | 100 μg |
| CTLA4 (CD152) Neutralizing Antibody | 71212 | 50 μg |
| B7-1, Biotin labeled | 71114 | 50 μg |
| B7-1 | 71125 | 100 μg |
| B7-2 | 71150 | 100 μg |
| CD28 | 71113 | 200 μg |
| CTLA4:B7-1[Biotinylated] Inhibitor Screening Assay Kit | 72009 | 96 rxns |
| CD28:B7-1[Biotinylated] Inhibitor Screening Assay Kit | 72007 | 96 rxns |
| PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit | 72003 | 96 rxns |
| PD-1:PD-L2[Biotinylated] Inhibitor Screening Assay Kit | 72004 | 96 rxns |
| PD-1 | 71106 | 100 μg |
| PD-1, Biotin labeled | 71109 | 50 μg |
| PD-L1 | 71104 | 100 μg |
| PD-L1, Biotin-labeled | 71105 | 50 μg |
| PD-L2 | 71107 | 100 μg |
| PD-L2, Biotin-labeled | 71108 | 50 μg |

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
 San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

TROUBLESHOOTING GUIDE

| Problem | Possible Cause | Solution |
|--|---|--|
| Luminescence signal of positive control reaction is weak | CTLA4 or B7-2 has lost binding capacity | Protein loses activity upon repeated freeze/thaw cycles. Use fresh B7-2-biotin, (BPS Bioscience #71150) and fresh CTLA4 (BPS Bioscience #71152). Store proteins in single-use aliquots. Increase time of protein incubation. Increase protein concentration. |
| | Incorrect settings on instruments | Refer to instrument instructions for settings to increase sensitivity of light detection. |
| | Chemiluminescent reagents mixed too soon | Chemiluminescent solution should be used within 15 minutes of mixing. Ensure both reagents are properly mixed. |
| | Inaccurate pipetting/technique | Run duplicates of all reactions. Use a multichannel pipettor. Use master mixes to minimize errors. |
| Luminescent signal is erratic or varies widely among wells | Bubbles in wells | Pipette slowly to avoid bubble formation. Tap plate lightly to disperse bubbles; be careful not to splash between wells. |
| | Insufficient washes | Increase number of washes. Increase wash volume. Add Tween-20 to 0.1% in washing buffer. |
| Background (signal to noise ratio) is high | 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 CTLA-4-biotin (BPS Bioscience #71152) 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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com