



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

### ***CD28:B7-2[Biotinylated] Inhibitor Screening Assay Kit***

**Catalog # 72062**  
**Size: 96 reactions**

**BACKGROUND:** The activation of naïve T cells requires two signals; the specific T cell receptor recognition of MHC/Antigen on the surface of the antigen-presenting cell (APC), and the binding of B7-2 (CD86) ligand on the APC with the CD28 receptor on the T cell surface. CD28:B7-2 interaction is an important drug target for the regulation of T cells involved in autoimmunity, inflammation, tumor recognition, and immune tolerance.

**DESCRIPTION:** The *CD28:B7-2[Biotinylated] Inhibitor Screening Assay Kit* is designed for screening and profiling inhibitors of CD28:B7-2 signaling. This kit comes in a convenient 96-well format, with biotin-labeled B7-2, purified CD28, streptavidin-labeled HRP, and assay buffer for 100 binding reactions. The key to this kit is the high sensitivity of detection of biotin-labeled B7-2 by streptavidin-HRP. Only a few simple steps on a microtiter plate are required for the assay. First, CD28 is coated on a 96-well plate. Next, B7-2 is incubated with CD28 on the plate. Finally, the plate is treated with streptavidin-HRP followed by addition of an HRP substrate to produce chemiluminescence, which can be measured using a chemiluminescence reader.

#### **COMPONENTS:**

Catalog #	Component	Amount	Storage	
71159	B7-2, Biotin-labeled	20 µg	-80°C	<b>(Avoid freeze/thaw cycles!)</b>
71113	CD28	20 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](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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 B7-2 binding to CD28.

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

**REFERENCES:**

1. Keir, M.E., *et al. Immunol. Rev.* 2005, **204**: 128-143.
2. Yao, S., *et al. Immunity* 2011, **34(5)**:729-40.

**ASSAY PROTOCOL:**

All samples and controls should be tested in duplicate.

**Coating the plate with CD28:**

- 1) Thaw **CD28** on ice. Upon first thaw, briefly spin tube containing **CD28** to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining **CD28** in aliquots at -80°C. *Note: CD28 is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.*
- 2) Dilute **CD28** to 4 ng/μl in PBS.
- 3) Add 50 μl of diluted **CD28** 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** with 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 to step 5.

**Step 1:**

- 1) Prepare the master mixture: N wells × (10 μl **3x Immuno Buffer 1** + 15 μl H<sub>2</sub>O).

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

- 2) Add 25  $\mu$ l of master mixture to each well. Use uncoated wells for the “Ligand Control”.
- 3) Add 5  $\mu$ l of inhibitor solution to each well designated “Test Inhibitor”. For the “Positive Control”, “Ligand Control” and “Blank”, add 5  $\mu$ l of the same solution without inhibitor (inhibitor buffer). Incubate at room temperature for one hour.
- 4) Thaw **B7-2-biotin** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **B7-2-biotin** into single use aliquots. Immediately store remaining undiluted enzyme in aliquots at  $-80^{\circ}\text{C}$ . *Note: B7-2-biotin is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*

	Blank	Ligand Control	Positive Control	Test Inhibitor
3x Immuno Buffer 1	10 $\mu$ l	10 $\mu$ l	10 $\mu$ l	10 $\mu$ l
H <sub>2</sub> O	15 $\mu$ l	15 $\mu$ l	15 $\mu$ l	15 $\mu$ l
Test Inhibitor/Activator	-	-	-	5 $\mu$ l
Inhibitor buffer (no inhibitor)	5 $\mu$ l	5 $\mu$ l	5 $\mu$ l	-
1x Immuno Buffer 1	20 $\mu$ l	-	-	-
B7-2-biotin (10 ng/ $\mu$ l)	-	20 $\mu$ l	20 $\mu$ l	20 $\mu$ l
<b>Total</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>

- 5) Dilute **B7-2-biotin** to 10 ng/ $\mu$ l in **1x Immuno Buffer 1**. Keep diluted protein on ice until use. Discard any unused diluted protein after use.
- 6) Add 20  $\mu$ l of **1x Immuno Buffer 1** to the well designated “Blank”.
- 7) Initiate reaction by adding 20  $\mu$ l of diluted **B7-2-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 three times with 100  $\mu$ l/well **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 9) Block wells by adding 100  $\mu$ l 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](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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  $\mu$ 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 plate onto clean paper towels to remove liquid.
- 4) Block wells by adding 100  $\mu$ 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  $\mu$ l **HRP Chemiluminescent Substrate A** and 50  $\mu$ l **HRP Chemiluminescent Substrate B**, then add 100  $\mu$ 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 uses 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 enzyme (typically we set this value as 100).

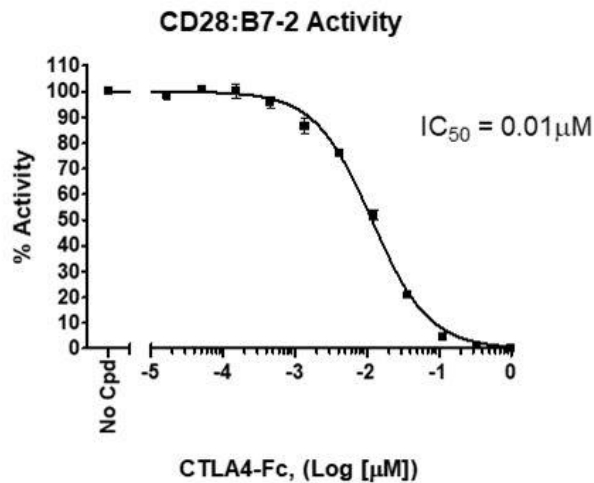
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

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



Inhibition of CD28:B7-2 interaction by CTLA4-Fc, Cat. #71149, using the *CD28:B7-2[Biotinylated] Inhibitor Screening Assay Kit*, Cat. #72062. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [support@bpsbioscience.com](mailto:support@bpsbioscience.com).

**RELATED PRODUCTS:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
CD28	71113	200 µg
B7-1	71125	100 µg
B7-1, Biotin labeled	71114	50 µg
B7-2	71150	100 µg
B7-2, Biotin labeled	71159	50 µg
CTLA-4	71149	100 µg
CTLA-4, Biotin labeled	71152	50 µg
CTLA4[Biotinylated]:B7-1 Inhibitor Screening Assay Kit	72009	96 rxns
CTLA4[Biotinylated]:B7-2 Inhibitor Screening Assay Kit	72024	96 rxns
CD28:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72007	96 rxns
PD-L1:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72026	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-L1, Biotin-labeled	71105	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](mailto:support@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://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	CD28 or B7-2 has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh B7-2-biotin, (BPS Bioscience #71159) and fresh CD28 (BPS Bioscience #71113). Store proteins in single-use aliquots. Increase time of enzyme incubation. Increase enzyme 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.
Luminescent 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.
Background (signal to noise ratio) is high	Insufficient washes	Increase number of washes. Increase wash volume. Increase Tween-20 concentration to 0.1% in PBST.
	Sample solvent is inhibiting the enzyme	Run negative control assay including solvent. Maintain DMSO level at <1% Increase time of enzyme incubation.
	Results are outside the linear range of the assay	Use different concentrations of B7-2-biotin (BPS Bioscience #71159) 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](mailto:support@bpsbioscience.com)

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