



6042 Cornerstone Ct. West, Ste. B  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet ***PD-1:PD-L2 Homogeneous Assay Kit*** Catalog #72015

**DESCRIPTION:** The *PD-1:PD-L2 Homogeneous Assay Kit* is designed to measure the inhibition of PD-1 binding to PD-L2. The *PD-1:PD-L2 Homogeneous Assay Kit* comes in a convenient AlphaLISA<sup>®</sup> format with purified biotinylated PD-L2, FLAG-tagged PD-1, and assay buffer to perform a total of 384 reactions. With this kit, only three simple steps on a microtiter plate are required. First, a sample containing PD-1 and an inhibitor of choice is incubated with the biotinylated PD-L2 for 60 minutes. Next, acceptor beads are added, then donor beads, followed by reading the Alpha-counts.

### COMPONENTS:

Catalog #	Component	Amount	Storage	
71198	PD-1-FLAG-Avi-His	10 µg	-80°C	<b>(Avoid freeze/ thaw cycles!)</b>
71108	PD-L2-biotin	5 µg	-80°C	
79311	3x Immuno Buffer 1	4 ml	-20°C	

### MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

AlphaLISA FLAG acceptor beads, 5 mg/ml (PerkinElmer #AL112C)  
AlphaScreen Streptavidin-conjugated donor beads, 5 mg/ml (PerkinElmer #6760002S)  
Optiplate-384 (PerkinElmer #6007290)  
AlphaScreen microplate reader  
Adjustable micropipettor and sterile tips

**APPLICATIONS:** Useful for screening for inhibitors of PD-1 binding to PD-L2

**CONTRAINDICATIONS:** Only limited amounts of DMSO can be included, as it has been shown to disrupt PD-1:PD-L2 interaction. Avoid green and blue dyes that absorb light in the AlphaScreen signal emission range (520-620 nm), such as Trypan Blue. Avoid the use of the potent singlet oxygen quenchers such as sodium azide (NaN<sub>3</sub>) or metal ions (Fe<sup>2+</sup>, Fe<sup>3+</sup>, Cu<sup>2+</sup>, Zn<sup>2+</sup> and Ni<sup>2+</sup>). The presence of >1% RPMI 1640 culture medium leads to a signal reduction due to the presence of excess biotin and iron in this medium. MEM, which lacks these components, does not affect AlphaScreen assays.

**STABILITY:** At least one year from date of receipt when stored as directed.

**REFERENCES:** 1. Molnar, E. *et al. Proc. Natl. Acad. Sci. U.S.A.* 2008; **105**: 10483-10488  
2. Keir, M.E. *et al. Annu. Rev. Immunol.* 2008, **26**: 677-704.

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)

**ASSAY PROTOCOL:**

All samples and controls should be tested in duplicate. Use slow shaking for all incubations.

**Step 1:**

- 1) Thaw **PD-1-FLAG-Avi-His** on ice. Upon first thaw, briefly spin tube containing protein to recover full contents of the tube. Aliquot the protein into single use aliquots. Store remaining undiluted protein in aliquots at -80°C immediately. *Note: PD-1-FLAG-Avi-His is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 2) Dilute one part **3x Immuno Buffer 1** with 2 parts of distilled water (3-fold dilution) to make **1x Immuno Buffer 1**. Make only a sufficient quantity needed for the assay; store remaining stock solution in aliquots at -20°C.
- 3) Dilute **PD-1-FLAG-Avi-His** in **1x Immuno Buffer 1** to 10 ng/μl. Keep diluted protein on ice until ready to use. Discard any remaining unused diluted protein after use.
- 4) Prepare the master mixture: N wells × (2 μl **3x Immuno Buffer 1** + 2 μl diluted **PD-1-FLAG** + 2 μl distilled water). Add 6 μl of master mixture to every well.

	Blank	Positive Control	Test Inhibitor
3x Immuno Buffer 1	2 μl	2 μl	2 μl
PD-1-FLAG-Avi-His (10 ng/μl)	2 μl	2 μl	2 μl
Distilled water	2 μl	2 μl	2 μl
Test Inhibitor	–	–	2 μl
Inhibitor buffer (no inhibitor)	2 μl	2 μl	–
1x Immuno Buffer 1	2 μl		
PD-L2-biotin (3 ng/μl)	–	2 μl	2 μl
<b>Total</b>	<b>10 μl</b>	<b>10 μl</b>	<b>10 μl</b>

- 5) Add 2 μl of inhibitor solution to each well designated “Test Inhibitor”. For the “Positive Control” and “Blank”, add 2 μl of the same solution without inhibitor (inhibitor buffer). *Note: If possible, keep final DMSO concentration below 0.5%.*
- 6) Add 2 μl of **1x Immuno Buffer 1** to the well designated “Blank”.
- 7) Thaw **PD-L2-biotin** on ice. Upon first thaw, briefly spin tube containing protein to recover full contents of the tube. Aliquot the protein into single use aliquots. Store remaining undiluted protein in aliquots at -80°C immediately. *Note: PD-L2-biotin is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*

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 Ct. West, Ste. B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

- 8) Dilute **PD-L2-biotin** in **1x Immuno Buffer 1** to 3 ng/μl. Keep diluted proteins on ice until use. Discard any remaining unused diluted protein after use.
- 9) Initiate reaction by adding 2 μl of diluted **PD-L2-biotin** prepared as described above to each well designated "Positive Control" and "Test Inhibitor". Incubate at room temperature for 60 minutes.

**Step 2:**

**Note: Protect your samples from direct exposure to light!**

- 1) Dilute FLAG Acceptor beads (PerkinElmer #AL112C) 250-fold with **1x Immuno Buffer 1**. Add 10 μl per well. Shake plate briefly. Incubate at room temperature for 30 minutes.

**Step 3:**

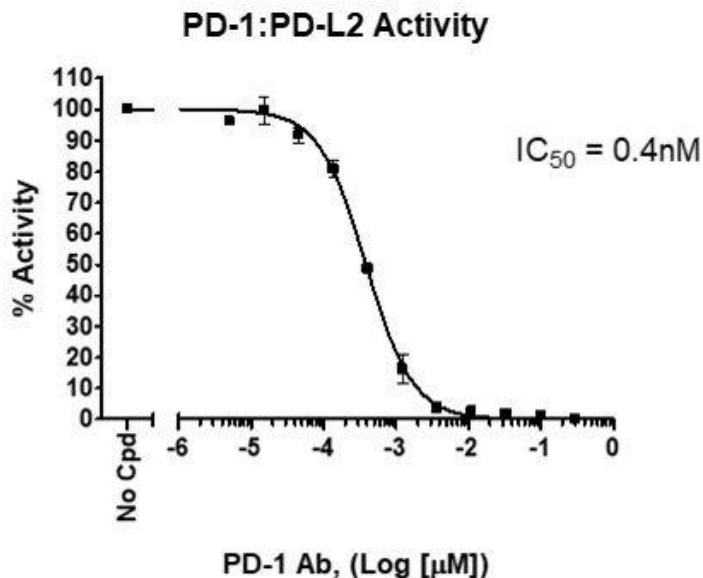
**Note: Protect your samples from direct exposure to light!**

- 1) Dilute Streptavidin-conjugated donor beads (PE #6760002S) 125-fold with **1x Immuno Buffer 1**. Add 10 μl per well. Incubate at room temperature for 30 minutes.
- 2) Read Alpha-counts.

*Due to lot to lot variability in AlphaScreen® bead performance, it may be necessary to optimize assay conditions. For example, slight adjustments to PD-1 or PD-1L concentrations may improve signal-to-noise ratio.*

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)

**Example of Assay Results:**


PD-1:PD-L2 inhibition, measured using the PD-1:PD-L2 Inhibitor Screening Assay Kit, BPS Bioscience, Catalog #72015 and PD-1 neutralizing antibody, Catalog #71213. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com).*

**RELATED PRODUCTS:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
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
PD-1:PD-L2[Biotinylated] Inhibitor Screening Kit	72004	96 rxns
PD-1:PD-L1[Biotinylated] Inhibitor Screening Kit	72003	96 rxns
PD-L1 Inhibitor Screening Kit	72005	96 rxns
PD-L2 Inhibitor Screening Kit	72006	96 rxns
PD-1 Neutralizing Antibody	71120	50 µg
PD-L1 Neutralizing Antibody	71213	50 µg
PD-1:PD-L1 Homogenous Assay Kit	72014	384 rxns

AlphaScreen® and AlphaLISA® are registered trademarks of PerkinElmer, Inc.

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)