



6044 Cornerstone Ct. West, Ste. E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

PD-1:PD-L1 Homogeneous Assay Kit

Catalog # 72014

DESCRIPTION: The *PD-1:PD-L1 Homogeneous Assay Kit* is designed to measure the inhibition of PD-1 binding to PD-L1. The *PD-1:PD-L1 Homogeneous Assay Kit* comes in a convenient AlphaLISA[®] format with purified biotinylated PD-L1, 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-L1 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	30 µg	-80°C	(Avoid freeze/ thaw cycles!)
71105	PD-L1-biotin	5 µg	-80°C	
79311	3x Immunobuffer 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-L1

CONTRAINDICATIONS: Only limited amounts of DMSO can be included, as it has been shown to disrupt PD-1-PD-L1 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 (NaN₃) or metal ions (Fe²⁺, Fe³⁺, Cu²⁺, Zn²⁺ and Ni²⁺). 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. Lin, D., *et al. Proc Natl Acad Sci U.S.A.* 2008, **105**: 3011-3016.
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 **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Ct. West, Ste. E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@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 Immunobuffer 1** with 2 parts of distilled water (3-fold dilution) to make **1x Immunobuffer 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 Immunobuffer 1** to 25 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 Immunobuffer 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 Immunobuffer 1	2 μl	2 μl	2 μl
PD-1-FLAG-Avi-His (25 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 Immunobuffer 1	2 μl		
PD-L1-biotin (3 ng/μl)	-	2 μl	2 μl
Total	10 μl	10 μl	10 μl

- 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 Immunobuffer 1** to the well designated “Blank”.
- 7) Thaw **PD-L1-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-L1-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 **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Ct. West, Ste. E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 8) Dilute **PD-L1-biotin** in **1x Immunobuffer 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-L1-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 Immunobuffer 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!

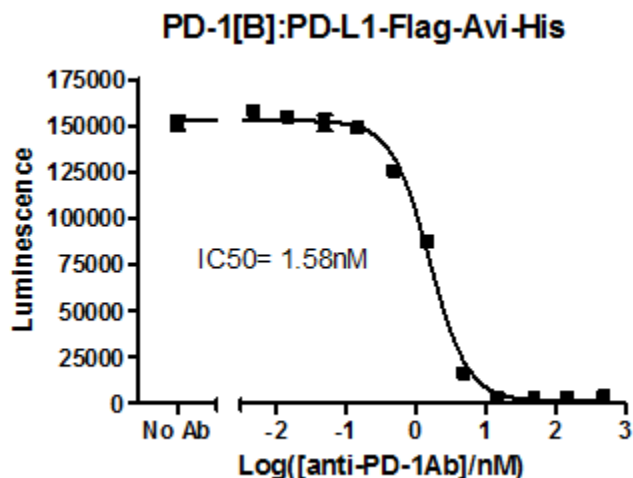
- 1) Dilute Streptavidin-conjugated donor beads (PE #6760002S) 125-fold with **1x Immunobuffer 1**. Add 10 μl per well. Incubate at room temperature for 30 minutes.
- 2) Read Alpha-counts.

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 **1.858.829.3082**, Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com

Example of Assay Results:



PD-1:PD-L1 inhibition, measured using the PD-1:PD-L1 Inhibitor Screening Assay Kit, BPS Bioscience, Catalog #72014 and PD-1 neutralizing antibody, Catalog #71120. *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>
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

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 **1.858.829.3082**, Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com