



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
CD112R:CD112[Biotinylated] Inhibitor Screening Assay Kit
Catalog #79732
Size: 96 reactions

BACKGROUND: CD112 (Poliovirus receptor-related 2, PVRL2), widely expressed on antigen-presenting cells and tumor cells, is the high affinity ligand of CD112R (Poliovirus receptor related immunoglobulin domain containing, PVRIG). CD112-CD112R interaction is a positive immune checkpoint that enhances human T cell response and it has emerged as an attractive therapeutic target for oncology. TIGIT (T cell immunoreceptor with Ig and ITIM domains) and CD226 (also called DNAM-1, DNAX Accessory Molecule-1) are other ligands of CD112.

DESCRIPTION: The *CD112R:CD112[Biotinylated] Inhibitor Screening Assay Kit* is designed for screening and profiling inhibitors of CD112R-CD112. This kit comes in a convenient 96-well format, with biotin-labeled CD112, purified CD112R, streptavidin-labeled HRP, and assay buffer for 100 binding reactions. The key to this kit is the high sensitivity of detection of biotin-labeled CD112 by streptavidin-HRP. Only a few simple steps on a microtiter plate are required for the assay. First, CD112R is coated on a 96-well plate. Next, CD112 is incubated with CD112R 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	
79116	CD112R (PVRIG), Fc-Fusion, Avi-Tag (Human) HiP™	10 µg	-80°C	Avoid multiple freeze/thaw cycles!
72231	CD112, Fc-Fusion, Avi-Tag, Biotin-labeled (Human)	5 µg	-80°C	
79311	3x Immuno Buffer 1	50 ml	-20°C	
	Blocking Buffer	50 ml	+4°C	
80611	Streptavidin-HRP	15 µl	-20°C	
79670	ELISA ECL Substrate A	6 ml	+4°C	
79670	ELISA ECL Substrate B	6 ml	+4°C	
79699	96-well white microplate	1	+4°C	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

PBS (Phosphate buffered saline)
Luminometer or fluorescent microplate reader capable of reading chemiluminescence
Adjustable micropipettor and sterile tips

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 screening for inhibitors of CD112 (PVRL2) binding to CD112R (PVRIG).

STABILITY: Up to 6 months from date of receipt, when stored as recommended.

REFERENCES:

Zhu, Y., *et al. J. Exp. Med.* 2016, **213(2)**:167-76
Torphy, R., *et al. Int. J. Mol. Sci.* 2017, **18(12)**: E2642

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Coating the plate with CD112R:

- 1) Thaw **CD112R** on ice. Upon first thaw, briefly spin tube containing **CD112R** to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining **CD112R** in aliquots at -80°C. Note: **CD112R** is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.
- 2) Dilute **CD112R** to 2 µg/ml in PBS.
- 3) Add 50 µl of diluted **CD112R** 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. Dilute only enough as required for this step of the assay.
- 5) Decant to remove supernatant. Wash the plate three 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 with slow shaking. Remove supernatant as described in step 5.

Step 1:

- 1) Prepare the master mixture: N wells × (10 µl **3x Immuno Buffer 1** + 15 µl distilled water)
- 2) Add 25 µl of master mixture to each well. Use uncoated wells for the "Ligand Control".
- 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 with slow shaking.
- 4) Thaw **CD112-biotin** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **CD112-biotin** into single use aliquots. Immediately store remaining undiluted enzyme in aliquots at - 80 °C. Note: **CD112-biotin** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
- 5) Dilute **CD112-biotin** to 1.6 ng/µl (25 nM) in **1x Immuno Buffer 1**. Keep diluted protein on ice until use. Discard any unused diluted protein after use.

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

- 6) Add 20 μ l of **1x Immuno Buffer 1** to the well designated “Blank”.

	Blank	Ligand Control	Positive Control	Test Inhibitor
3x Immuno Buffer	10 μ l	10 μ l	10 μ l	10 μ l
H ₂ O	15 μ l	15 μ l	15 μ l	15 μ l
Test Inhibitor	-	-	-	5 μ l
Inhibitor buffer (no inhibitor)	5 μ l	5 μ l	5 μ l	-
1x Immuno Buffer 1	20 μ l	-	-	-
CD112-biotin (1.6 ng/ μ l)	-	20 μ l	20 μ l	20 μ l
Total	50 μl	50 μl	50 μl	50 μl

- 7) Initiate reaction by adding 20 μ l of diluted **CD112-biotin** (see Step 1-5) to wells labeled “Positive Control”, “Ligand Control” and “Test Inhibitor”. Incubate at room temperature for two hours with slow shaking.
- 8) Decant to remove supernatant. Wash the plate 3 times with 100 μ l/well **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 9) Block wells by adding 100 μ l of **Blocking Buffer** to each well. Incubate for 10 minutes at room temperature. Remove supernatant as in Step 1-8.

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 plate onto clean paper towel 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 **ELISA ECL Substrate A** and 50 μ l **ELISA ECL Substrate B**, 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.

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

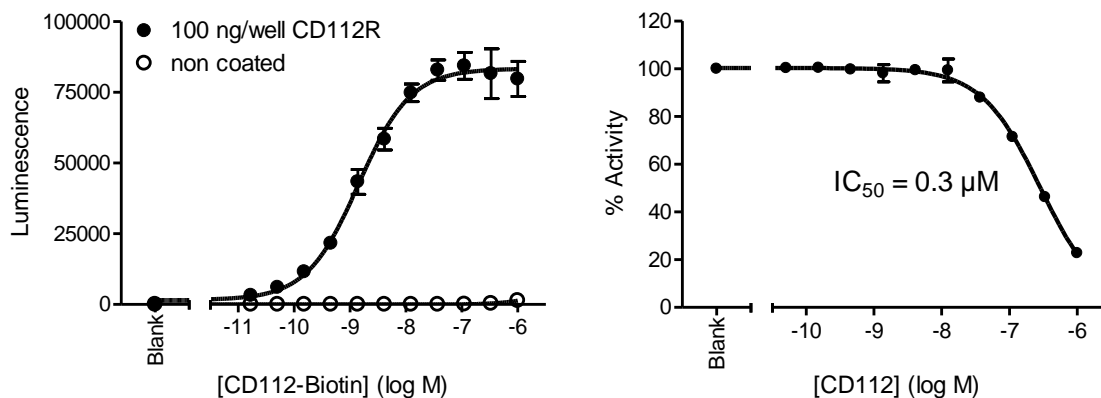
Please visit our website at: www.bpsbioscience.com

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 milliseconds. 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).

Example of assay results:



CD112R:CD112 binding activity, measured using the using the *CD112R:CD112[Biotinylated] Inhibitor Screening Assay Kit*, BPS Bioscience, #79732 (left). Inhibition of CD112R:CD112 binding using the non-biotinylated CD112, BPS Bioscience, #11079 in the *CD112R:CD112[Biotinylated] Inhibitor Screening Assay Kit* (right). 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>
CD112, His-tag (Human) HiP™	71197	100 μg
CD112, His-tag, Biotin-labeled (Human) HiP™	71234	50 μg
CD112, Fc-Fusion, Avi-Tag, Biotin-labeled (Human)	72231	50 μg
CD112R (PVRIG), Fc-Fusion, Avi-Tag (Human) HiP™	79116	100 μg
CD112R (PVRIG), Fc-Fusion, Avi-Tag, Biotin (Human) HiP™	79270	50 μg
TIGIT:CD112 Homogeneous Assay Kit	72030	384 reactions
CD226:CD112 Homogeneous Assay Kit	72051	384 reactions

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

TROUBLESHOOTING GUIDE

Problem	Possible cause	Solution
Luminescence signal of positive control reaction is weak	CD112R or CD112 has lost activity	Proteins lose activity upon repeated freeze/thaw cycles. Use fresh CD112-biotin, (BPS Bioscience #72231) and fresh CD112R (BPS Bioscience #79116). 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 CD112-Biotin (BPS Bioscience #72231) 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