



6042 Cornerstone Ct. West, Ste. B  
San Diego, CA 92121  
Tel: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### ***TIGIT:CD112 Homogeneous Assay Kit***

**Catalog #72030**

**Size: 384 reactions**

**BACKGROUND:** Human T-cell immunoreceptor with Ig and ITIM domains (TIGIT) is a receptor that is expressed on the surface of human T cells and NK cells that binds to CD155 and CD112 on the surface of dendritic cells. Binding of TIGIT with CD155 or CD112 results in inhibition of T cell and NK cell activation. Antibodies and other agents that inhibit this signaling pathway have been shown to increase the immune response, especially in the case of certain cancers.

**DESCRIPTION:** The *TIGIT:CD112 Homogeneous Assay Kit* is designed to measure the inhibition of human T-cell immunoreceptor with Ig and ITIM domains (TIGIT) binding to CD112 (PVRL2/Nectin-2). The *TIGIT:CD112 Homogeneous Assay Kit* comes in a convenient AlphaLISA<sup>®</sup> format with purified biotinylated TIGIT, His-tagged CD112, 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 TIGIT and an inhibitor of choice is incubated with the CD112 for 60 minutes. Next, acceptor beads are added, then donor beads, followed by reading the Alpha-counts.

#### **COMPONENTS:**

Catalog #	Component	Amount	Storage	
71251	TIGIT-Fc-biotin	2x 5 µg	-80°C	<b>(Avoid freeze/ thaw cycles!)</b>
71197	CD112-His	2x 6 µg	-80°C	
79311	3x Immuno Buffer 1	4 ml	-20°C	

#### **MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

AlphaLISA Ni Chelate Acceptor beads, 5 mg/ml (PerkinElmer #AL108C)  
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 TIGIT binding to CD112

**CONTRAINDICATIONS:** Only limited amounts of DMSO can be included, as it has been shown to disrupt TIGIT:CD112 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

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

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. Yu, X., *et al.*, *Nat. Immunol.* 2009; **10(1)**: 48-57.  
 2. Stanietzky, N., *et al.*, *Proc. Natl. Acad. Sci.* 2009; **106(42)**: 17858-17863.

**ASSAY PROTOCOL:**

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

**Step 1:**

- 1) Thaw **CD112-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: CD112-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 **CD112-His** in **1x Immuno Buffer 1** to 5 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 **CD112-His** + 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
CD112-His (5 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		
TIGIT-biotin (4 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%.*

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](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: 1.858.829.3082  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

- 6) Add 2  $\mu$ l of **1x Immuno Buffer 1** to the well designated "Blank".
- 7) Thaw **TIGIT-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  $-80^{\circ}\text{C}$  immediately. *Note: **TIGIT-biotin** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 8) Dilute **TIGIT-biotin** in **1x Immuno Buffer 1** to 4 ng/ $\mu$ l. Keep diluted proteins on ice until use. Discard any remaining unused diluted protein after use.
- 9) Initiate reaction by adding 2  $\mu$ l of diluted **TIGIT-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 Ni Chelate Acceptor beads (PerkinElmer #AL108C) 250-fold with **1x Immuno Buffer 1**. Add 10  $\mu$ l per well. Shake plate briefly. Incubate at room temperature for 30 minutes.

#### **Step 3:**

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

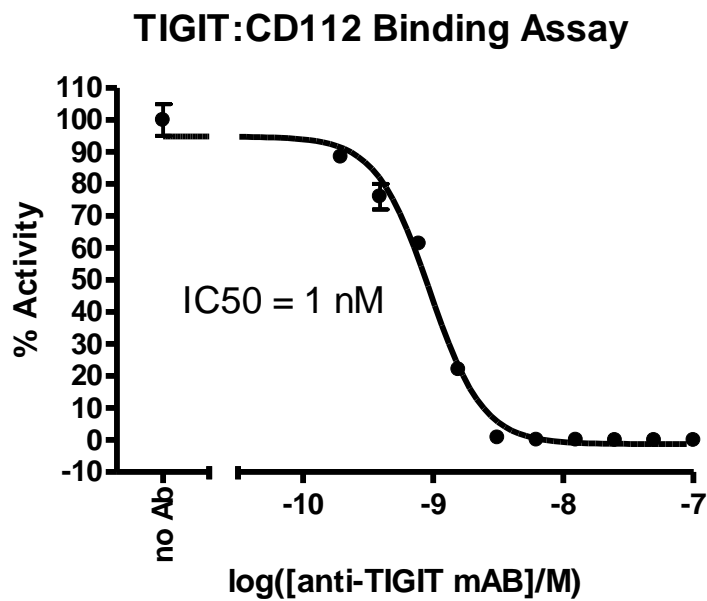
- 1) Dilute Streptavidin-conjugated donor beads (PE #6760002S) 125-fold with **1x Immuno Buffer 1**. Add 10  $\mu$ l per well. Incubate at room temperature for 30 minutes.
- 2) Read Alpha-counts.

*Due to lot to lot variability in AlphaScreen<sup>®</sup> bead performance, it may be necessary to optimize assay conditions. For example, slight adjustments to TIGIT-biotin or CD112-His 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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

**Example of Assay Results:**



TIGIT:CD112 inhibition, measured using the TIGIT:CD112 Inhibitor Screening Assay Kit, BPS Bioscience, Catalog #72030 and an anti-TIGIT antibody (BPS Cat. #71218). *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>
Human TIGIT	#71218	100 µg
Human TIGIT, Fc fusion, Biotin-labeled	#71251	50 µg
Human TIGIT, Fc fusion	#71186	100 µg
Human CD112, His-tag	#71197	100 µg
Human CD112, His-tag, Biotin-labeled	#71234	50 µg
Human CD155 (PVR), His-tag	#71181	100 µg
Mouse CD155 (PVR), His-tag	#71167	100 µg
Mouse CD155 (PVR), His-tag, Biotin-labeled	#71168	50 µg
TIGIT:CD155 Homogeneous Assay Kit	#72029	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 **1.858.829.3082**, 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)