



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

### ***TIGIT:CD155 Homogeneous Assay Kit***

**Catalog #72029**

**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:CD155 Homogeneous Assay Kit* is designed to measure the inhibition of TIGIT binding to CD155. The *TIGIT:CD155 Homogeneous Assay Kit* comes in a convenient AlphaLISA<sup>®</sup> format with purified biotinylated TIGIT, His-tagged CD155, 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 CD155 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 3 µg	-80°C	<b>(Avoid freeze/ thaw cycles!)</b>
71181	CD155-His	2x 5 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 CD155

**CONTRAINDICATIONS:** Only limited amounts of DMSO can be included, as it has been shown to disrupt TIGIT:CD155 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.

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)

**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 **CD155-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: CD155-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 **CD155-His** in **1x Immuno Buffer 1** to 4 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 **CD155-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
CD155-His (4 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 (2 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”.

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)

- 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°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 2 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 **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 μl per well. Shake plate briefly. Incubate at room temperature for 30 minutes.

**Step 3:**

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

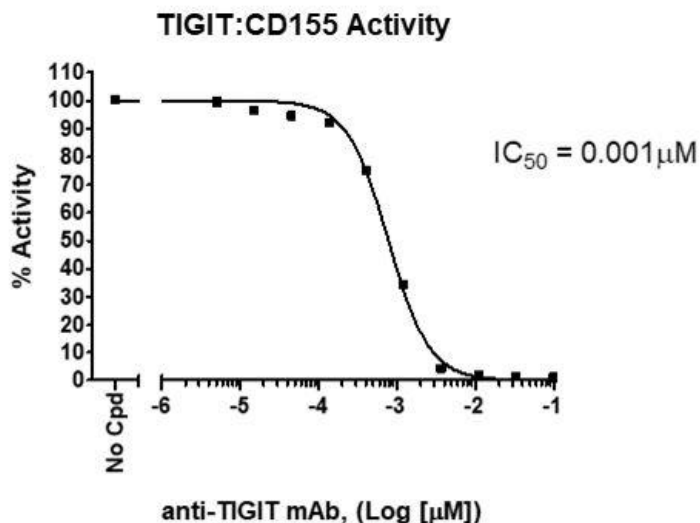
- 1) Dilute Streptavidin-conjugated donor beads (PE #6760002S) 125-fold with **1x Immuno Buffer 1**. Add 10 μl per well. Incubate at room temperature for 30 minutes.
- 2) Read Alpha-counts.

*Due to lot to lot variability in AlphaScreen® bead performance, it may be necessary to optimize assay conditions. For example, slight adjustments to TIGIT-biotin or CD155-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 **1.858.202.1401** 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:**



TIGIT:CD155 inhibition, measured using the TIGIT:CD155 Inhibitor Screening Assay Kit, BPS Bioscience, Catalog #72029 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 $\mu g$
Human TIGIT, Fc fusion, Biotin-labeled	#71251	50 $\mu g$
Human TIGIT, Fc fusion	#71186	100 $\mu g$
Human CD112, His-tag	#71197	100 $\mu g$
Human CD112, His-tag, Biotin-labeled	#71234	50 $\mu g$
Human CD155 (PVR), His-tag	#71181	100 $\mu g$
Mouse CD155 (PVR), His-tag	#71167	100 $\mu g$
Mouse CD155 (PVR), His-tag, Biotin-labeled	#71168	50 $\mu g$
TIGIT:CD112 Homogeneous Assay Kit	#72030	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)