



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

Data sheet
CD27:CD70[Biotinylated] Inhibitor Screening Assay Kit
Catalog #79695
Size: 96 reactions

BACKGROUND:

The tumor necrosis factor (TNF) receptor family member CD27 (TNFRSF7) and its TNF-like ligand CD70 (CD27L, TNFSF7) are a key costimulatory immune checkpoint signal for lymphocyte activation. Among other effects, this interaction promotes the differentiation of T and B cells and participates in immunoglobulin synthesis. CD27:CD70 interaction has been demonstrated to have important roles in autoimmunity, inflammation, infection and cancer, and antibodies targeting either CD27 or CD70 are in clinical trials for oncology.

DESCRIPTION: The *CD27:CD70[Biotinylated] Inhibitor Screening Assay Kit* is designed for screening and profiling inhibitors of CD27:CD70 signaling. This kit comes in a convenient 96- well format, with biotin-labeled CD70, purified CD27, streptavidin-labeled HRP, and assay buffer for 100 binding reactions. The key to this kit is the high sensitivity of detection of biotin-labeled CD70 by streptavidin-HRP. Only a few simple steps on a microtiter plate are required for the assay. First, CD27 is coated on a 96-well plate. Next, CD70 is incubated with CD27 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	
71176	CD27, Fc fusion (Human)	10 µg	-80 °C	Avoid multiple freeze/thaw cycles!
100228	CD70 His-Avi-Tag, Biotin-Labeled	3 µg	-80 °C	
79311	3x Immuno Buffer 1	50 ml	-20 °C	
	Blocking Buffer	50 ml	+4 °C	
	Streptavidin-HRP	15 µl	-20 °C	
79670	ECL Substrate A (transparent bottle)	6 ml	+4 °C	
79670	ECL Substrate B (transparent bottle)	6 ml	+4 °C	
	96-well white microplate	1	+4 °C	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

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

APPLICATIONS: This kit is useful for screening for inhibitors of CD70 binding to CD27.

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

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

REFERENCES:

Granger, S.W., *et al. Cytokine Growth Factor Rev.* 2003, **14(3-4)**: 289-296

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Coating the plate with CD27:

- 1) Thaw **CD27** on ice. Upon first thaw, briefly spin tube containing **CD27** to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining **CD27** in aliquots at -80°C. Note: **CD27** is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.
- 2) Dilute **CD27** to 2 µg/ml in PBS.
- 3) Add 50 µl of dilute **CD27** 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.
- 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 **CD70-biotin** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **CD70-biotin** into single use aliquots. Immediately store remaining undiluted enzyme in aliquots at - 80 °C. Note: **CD70-biotin** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.

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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

	Blank	Ligand Control	Positive Control	Test Inhibitor
3x Immuno Buffer	10 µl	10 µl	10 µl	10 µl
Distilled water	15 µl	15 µl	15 µl	15 µl
Test Inhibitor	-	-	-	5 µl
Inhibitor buffer (no inhibitor)	5 µl	5 µl	5 µl	-
1x Immuno Buffer 1	20 µl	-	-	-
CD70-biotin (0.25 ng/µl)	-	20 µl	20 µl	20 µl
Total	50 µl	50 µl	50 µl	50 µl

- 5) Dilute **CD70-biotin** to 0.25 ng/µl (12.5 nM) in **1x Immuno Buffer 1**. Keep diluted protein on ice until use. Discard any unused diluted protein after use.
- 6) Add 20 µl of **1x Immuno Buffer 1** to the well designated "Blank".
- 7) Initiate reaction by adding 20 µl of diluted **CD70-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 µl/well **1x Immuno Buffer 1**. Tap plate onto clean paper towels to remove liquid.
- 9) Block wells by adding 100 µl 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 µl 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 µl 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 µl **ECL Substrate A** and 50 µl **ECL Substrate B**, then add 100 µl 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.

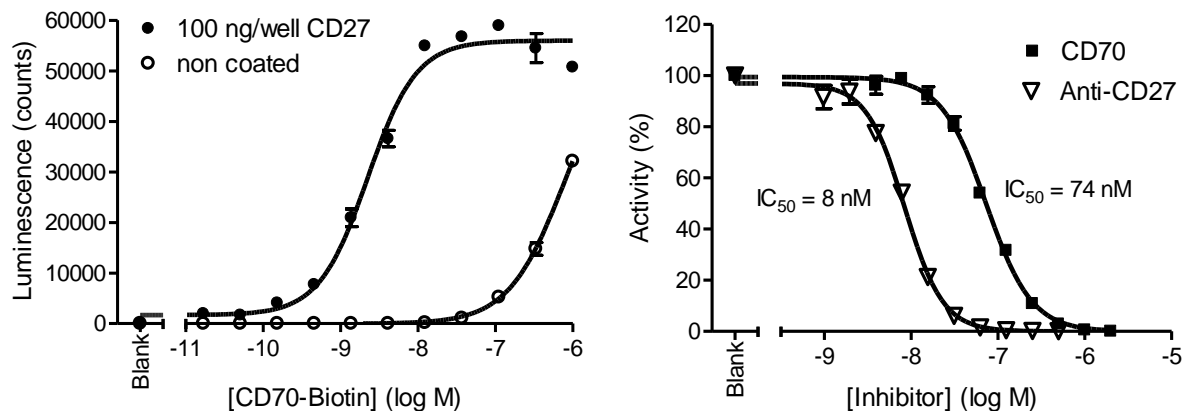
To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: support@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com

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 mseconds. 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:



CD27:CD70 binding activity, measured using the using the *CD27:CD70[Biotinylated] Inhibitor Screening Assay Kit*, BPS Bioscience #79695 (left). Inhibition of CD27:CD70 binding using the non-biotinylated CD70, BPS Bioscience #71178, and the Anti-CD27 Agonist Antibody, BPS Bioscience #100111, in the *CD27:CD70 [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.

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: support@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog#</u>	<u>Size</u>
CD27, Fc fusion (Human)	71176	100 µg
CD27, Fc Fusion (IgG1) Avi-Tag, Biotin Labeled	100250	50 µg
CD27, Fc-Fusion (IgG1), Avi-Tag (Mouse) HiP™	79067	100 µg
CD27 (Mouse), Fc-Fusion (Human), Avi-Tag, Biotin	79107	50 µg
Anti-CD27 Agonist Antibody	100111	100 µg
CD27 CHO-K1 Stable Recombinant Cell Line	60624	2 vials
CD70, His-tag (Human)	71178	100 µg
CD70, His-Avi-Tag, Biotin-Labeled	100228	50 µg
CD70, His-Tag (Mouse)	79066	100 µg
CD70-CHO Recombinant Cell line	79510	2 vials
PD-1:PD-L1[Biotinylated] Inhibitor Screening Assay Kit	72003	96 reactions
PD-1:PD-L2[Biotinylated] Inhibitor Screening Assay Kit	72004	96 reactions
PD-L1 Inhibitor Screening Assay Kit	72005	96 reactions
PD-L2 Inhibitor Screening Assay Kit	72006	96 reactions
CD28:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72007	96 reactions
BTLA:HVEM[Biotinylated] Inhibitor Screening Assay Kit	72008	96 reactions
CTLA4:B7-1[Biotinylated] Inhibitor Screening Assay Kit	72009	96 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: support@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
 San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

TROUBLESHOOTING GUIDE

Problem	Possible cause	Solution
Luminescence signal of positive control reaction is weak	CD27 or CD70 has lost activity	Proteins lose activity upon repeated freeze/thaw cycles. Use fresh CD70-biotin, (BPS Bioscience #100228) and fresh CD27 (BPS Bioscience #71176). Store proteins in single-use aliquots. Increase time of enzyme incubation. Increase enzyme concentration.
	Antibody reaction is insufficient	Increase time for primary antibody incubation. Avoid freeze/thaw cycles of antibodies.
	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 CD70-Biotin (BPS Bioscience #100228) 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: support@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com