



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
BCMA:APRIL[Biotinylated] Inhibitor Screening Assay Kit
Catalog #79722
Size: 96 reactions

BACKGROUND: A proliferation-inducing ligand (APRIL, TNSF13 or CD256) and B-cell maturation antigen (BCMA, also known as TNFRSF17 or CD269) create a signaling pathway that participates in the regulation of B-cell development, autoimmunity and long-term plasma cell survival. APRIL expression is significantly elevated in multiple myeloma (MM) and it has been demonstrated that activation of BCMA by APRIL promotes tumor growth, chemoresistance and immunosuppression in the bone marrow microenvironment. Anti-APRIL monoclonal antibodies have been recently progressed to clinical trials for MM.

DESCRIPTION: The *BCMA:APRIL[Biotinylated] Inhibitor Screening Assay Kit* is designed for screening and profiling inhibitors of BCMA:APRIL. This kit comes in a convenient 96-well format, with biotin-labeled APRIL, purified BCMA, streptavidin-labeled HRP, and assay buffer for 100 binding reactions. The key to this kit is the high sensitivity of detection of biotin-labeled APRIL by streptavidin-HRP. Only a few simple steps on a microtiter plate are required for the assay. First, BCMA is coated on a 96-well plate. Next, biotinylated APRIL is incubated with BCMA 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 | |
|-----------|------------------------------------|--------|------------|------------------------------------|
| 79465 | BCMA, Fc-Fusion, Avi-Tag HiP™ | 10 µg | -80 °C | Avoid multiple freeze/thaw cycles! |
| 100262 | APRIL, His-Avi-Tag, Biotin-Labeled | 5 µg | -80 °C | |
| 79311 | 3x Immuno Buffer 1 | 50 ml | -20 °C | |
| 79728 | Blocking Buffer 2 | 50 ml | +4 °C | |
| 79742 | Streptavidin-HRP | 15 µl | +4 °C | |
| 79670 | ELISA ECL Substrate A | 6 ml | Room Temp. | |
| | ELISA ECL Substrate B | 6 ml | Room Temp. | |
| 79699 | 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 BCMA binding to APRIL.

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

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

REFERENCES:

1. Yu, G., *et al. Nature Immunol.* 2000, **1(3)**: 252
2. Tai Y, T., *et al. Blood* 2016, **127**: 3225-3236.
3. Cho, S.-F., *et al. Front. Immunol.* 2018, **9**:1821

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Coating the plate with BCMA:

- 1) Thaw **BCMA** on ice. Upon first thaw, briefly spin tube containing **BCMA** to recover the full contents of the tube. Aliquot into single use aliquots. Immediately store remaining **BCMA** in aliquots at -80°C. Note: **BCMA** is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.
- 2) Dilute **BCMA** to 2 µg/ml in PBS.
- 3) Add 50 µl of diluted **BCMA** 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 is required for the washing steps below.
- 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 2** 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 **APRIL-biotin** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **APRIL-biotin** into single use aliquots. Immediately store remaining undiluted enzyme in aliquots at -80°C. Note: **APRIL-biotin** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
- 5) Dilute **APRIL-biotin** to 0.25 ng/µl (5 ng/20 µl) 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: 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

- 6) Add 20 µl of 1x Immuno Buffer 1 to the well designated "Blank".

| | 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 | - | - | - |
| APRIL-biotin (0.25 ng/µl) | - | 20 µl | 20 µl | 20 µl |
| Total | 50 µl | 50 µl | 50 µl | 50 µl |

- 7) Initiate reaction by adding 20 µl of diluted **APRIL-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 2** 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**.
- 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 2** 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 **ELISA ECL Substrate A** and 50 µl **ELISA 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



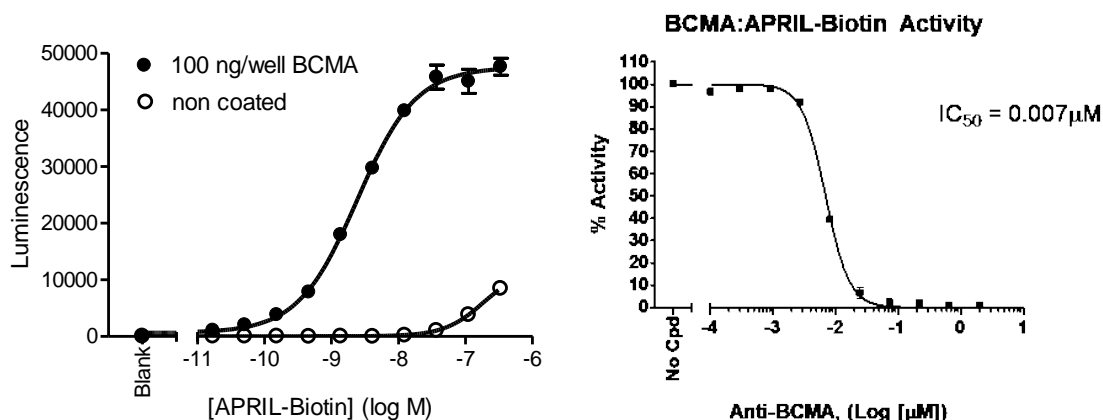
6405 Mira Mesa Blvd Ste. 100
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: support@bpsbioscience.com

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:



BCMA:APRIL binding activity, measured using the using the *BCMA:APRIL[Biotinylated] Inhibitor Screening Assay Kit*, BPS Bioscience #79722 (left). Inhibition of BCMA:APRIL binding using the Anti-BCMA Antibody (scFv), His-Tag, BPS Bioscience #100173 in the *BCMA:APRIL[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 @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> |
|---|------------------------|--------------------|
| BCMA, Fc-Fusion, Avi-Tag HiP™ | 79465 | 100 µg |
| BCMA, Fc-fusion (IgG1), Avi-Tag, Biotin-Labeled HiP™ | 79467 | 50 µg |
| APRIL, His-Avi-Tag, Biotin-Labeled | 100262 | 50 µg |
| Anti-BCMA Antibody (scFv), His-Tag | 100173-1 | 50 µg |
| BAFF:BCMA[Biotinylated] Inhibitor Screening Assay Kit | 79667 | 96 reactions |
| BCMA CHO Recombinant Cell Line (High Expression) | 79500-H | 2 vials |
| BCMA CHO Recombinant Cell Line (Medium Expression) | 79500-M | 2 vials |
| BCMA CHO Recombinant Cell Line (Low Expression) | 79500-L | 2 vials |
| ELISA ECL Substrate | 79670-1 | 200 ml |
| Immuno Buffer 1 | 79311 | 50 ml |

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 | BCMA or APRIL has lost activity | Proteins lose activity upon repeated freeze/thaw cycles. Use fresh APRIL-biotin, (BPS Bioscience #100262) and fresh BCMA (BPS Bioscience #79465). 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 APRIL-Biotin (BPS Bioscience #100262) 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