



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
PRMT3 Chemiluminescent Activity Assay Kit
Catalog # 52005L
Size: 96 reactions

DESCRIPTION: The *PRMT3 Chemiluminescent Activity Assay kit* is designed to measure PRMT3 activity for screening and profiling applications. The *PRMT3 Chemiluminescent Activity Assay Kit* comes in a convenient format, with a 96-well plate precoated with specific histone substrate, the antibody against methylated substrate, the secondary HRP-labeled antibody, S-adenosylmethionine, methyltransferase assay buffer, and purified PRMT3 enzyme for 96 enzyme reactions. The key to the *PRMT3 Chemiluminescent Activity Assay Kit* is a highly specific antibody that recognizes methylated histone substrate. With this kit, only three simple steps on a microtiter plate are required for methyltransferase detection. First, S-adenosylmethionine is incubated with a sample containing assay buffer and methyltransferase enzyme. Next, primary antibody is added. Finally, the strip plates are treated with an HRP-labeled secondary antibody followed by addition of the HRP substrate to produce chemiluminescence that can then be measured using a chemiluminescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
51043	PRMT3 human recombinant enzyme	10 µg	-80°C	(Avoid freeze/thaw cycles!)
52120	20 µM S-adenosylmethionine*	250 µl	-80°C	
52150	Primary antibody 4	100 µl	-80°C	
52131H	Secondary HRP-labeled antibody 2	10 µl	-80°C	
52170	4x HMT assay buffer 2	3 ml	-20°C	
52100	Blocking buffer 4	50 ml	+4°C	
	HRP chemiluminescent A (transparent bottle)	6 ml	+4°C	
	HRP chemiluminescent B (brown bottle)	6 ml	+4°C	
	96-well plate precoated with histone substrate	1 plate	+4°C	

* Decreasing S-adenosylmethionine concentration will make the assay more sensitive to inhibitors.

MATERIALS REQUIRED BUT NOT SUPPLIED:

TBST buffer (1 x TBS, pH 8.0, containing 0.05% Tween20)
Luminometer or microplate reader capable of reading chemiluminescence
Adjustable micropipettor and sterile tips
Rotating or rocker platform

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

APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

CONTRAINDICATIONS: DMSO >1%, strong acids or bases, ionic detergents, high salt

STABILITY: One year from date of receipt when stored as directed.

REFERENCE: Dillon SC, Zhang X, Trievel RC, Cheng X. *Genome Biology* 2005; **6**:227.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Step 1:

- 1) Rehydrate the microwells by adding 150 μ l of TBST buffer (1x TBS, pH 8.0, containing 0.05% Tween-20) to every well. Incubate 15 minutes at room temperature. Tap the plate onto clean paper towels to remove liquid.
- 2) Thaw **S-adenosylmethionine** on ice. Upon first thaw, briefly spin tube containing **S-adenosylmethionine** to recover full contents of the tube. Aliquot **S-adenosylmethionine** into single-use aliquots and store at -80°C . *Note: S-adenosylmethionine is very sensitive to freeze/thaw cycles. Avoid multiple freeze/thaw cycles.*
- 3) Prepare the master mixture: N wells \times (7.5 μ l **4x HMT assay buffer 2** + 2.5 μ l **20 μ M S-adenosylmethionine** + 15 μ l water). Add 25 μ l of master mixture to all wells labeled "Positive Control", "Test Sample" and "Blank". For wells labeled "Substrate control", add 7.5 μ l **4x HMT assay buffer 2** + 17.5 μ l water.

	Blank	Substrate Control	Positive Control	Test Sample
4x HMT assay buffer 2	7.5 μ l	7.5 μ l	7.5 μ l	7.5 μ l
20 μ M S-adenosylmethionine	2.5 μ l	-	2.5 μ l	2.5 μ l
H ₂ O	15 μ l	17.5 μ l	15 μ l	15 μ l
Test Inhibitor/Activator	-	-	-	5 μ l
Inhibitor buffer (no inhibitor)	5 μ l	5 μ l	5 μ l	
1X HMT assay buffer 2	20 μ l	-	-	-
Diluted PRMT3 (0.25-1 ng/ μ l)		20 μ l	20 μ l	20 μ l
Total	50 μl	50 μl	50 μl	50 μl

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

- 4) Add 5 μ l of inhibitor solution of each well designated "Test Inhibitor".
- 5) For the "Positive Control", "Substrate Control" and "Blank", add 5 μ l of the same solution without inhibitor (inhibitor buffer).
- 6) Thaw **PRMT3 enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **PRMT3 enzyme** into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C . *Note: **PRMT3 enzyme** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 7) Dilute **PRMT3 enzyme** in **1x HMT assay buffer 2** to 0.25-1 ng/ μ l (5-20 ng/20 μ l). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use. *Note: Diluted enzyme may not be stable. Dilute the enzyme immediately before use.*
- 8) Add 20 μ l of **1x HMT assay buffer 2** to the wells designated "Blank".
- 9) Initiate reaction by adding 20 μ l of diluted **PRMT3 enzyme** to the wells designated "Positive Control", "Substrate Control", and "Test Sample ". Incubate at room temperature for 1 hour.
- 10) Remove the supernatant from the wells and wash the strip three times with 200 μ l TBST buffer. Blot dry onto clean paper towels.
- 11) Add 100 μ l of **Blocking buffer 4** to every well. Shake on a rotating platform for 10 min. Remove supernatant as above.

Step 2:

- 1) Dilute "**Primary antibody 4**" 100-fold with **Blocking buffer 4**.
- 2) Add 100 μ l per well. Incubate 1 hour at room temperature with slow shaking.
- 3) Remove the supernatant from the wells and wash the strip three times with 200 μ l TBST buffer and incubate in **Blocking buffer 4** as described in steps 1-10 and 1-11.

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

Step 3:

- 1) Dilute "**Secondary HRP-labeled antibody 2**" 1,000-fold with **Blocking buffer 4**.
- 2) Add 100 μ l per well. Incubate for 30 min. at room temperature with slow shaking.
- 3) Wash plate with TBST buffer and **Blocking buffer 4** as in steps 1-10 and 1-11.
- 4) Just before use, mix on ice 50 μ l **HRP chemiluminescent substrate A** and 50 μ l **HRP chemiluminescent substrate B** and add 100 μ l per well. Discard any unused chemiluminescent reagent after use.
- 5) Immediately read sample in a luminometer or microtiter-plate capable of reading chemiluminescence. "Blank" value is subtracted from all other values.

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 msec. 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).

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**

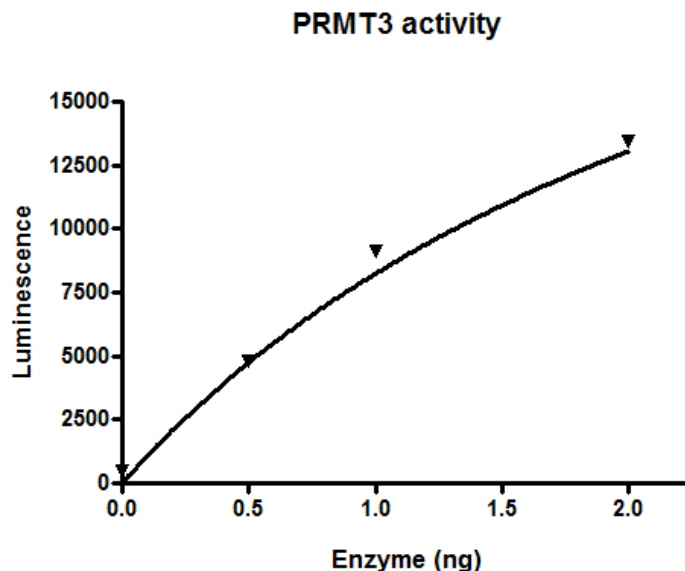
Or you can Email us at: support@bpsbioscience.com

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

Example of Assay Results:



PRMT3 enzyme activity, measured using the PRMT3 Chemiluminescent Assay Kit, BPS Bioscience #52005L. 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</u>	<u>Catalog</u>	<u>Size</u>
PRMT1 (expressed in E. coli)	#51040	50 µg
PRMT1 (expressed in Sf9 cells)	#51041	20 µg
PRMT3 (expressed in E. coli)	#51043	50 µg
PRMT4 (expressed in HEK293)	#51047	20 µg
PRMT4 (expressed in Sf9 cells)	#51044	20 µg
PRMT5 (expressed in HEK293)	#51045	20 µg
PRMT5 (expressed in Sf9 cells)	#51048	20 µg
PRMT6 (expressed in HEK293)	#51046	20 µg
PRMT8 (expressed in Sf9 cells)	#51052	20 µg
PRMT1 Chemiluminescent Assay Kit	#52004L	96 reactions
PRMT4 Chemiluminescent Assay Kit	#52041L	96 reactions
PRMT5 Chemiluminescent Assay Kit	#52002L	96 reactions
PRMT6 Chemiluminescent Assay Kit	#52046	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	PRMT3 enzyme has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh enzyme (PRMT3, BPS Bioscience #51043). Store enzyme 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 TBST.
	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 enzyme (PRMT3, BPS Bioscience #51043) 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