



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

PRMT8 Chemiluminescent Assay Kit

Catalog # 52062

Size: 96 reactions

DESCRIPTION: The *PRMT8 Chemiluminescent Assay Kit* is designed to measure PRMT8 activity for screening and profiling applications. The *PRMT8 Chemiluminescent Assay Kit* comes in a convenient format with a 96-well plate precoated with histone H4 peptide substrate, the antibody against methylated arginine 3 (R3) residue of Histone H4, the secondary HRP-labeled antibody, S-adenosylmethionine, methyltransferase assay buffer, and purified PRMT8 enzyme for 100 enzyme reactions. The key to the *PRMT8 Chemiluminescent Assay Kit* is a highly specific antibody that recognizes methylated R3 residue of Histone H4. 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 plate is 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 #	Components	Amount	Storage	
51052	PRMT8 His-FLAG-tags	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	
79670	ELISA ECL chemiluminescent substrate A (transparent bottle)	6 ml	Room Temp	
	ELISA ECL chemiluminescent substrate B (brown bottle)	6 ml	Room Temp	
	96-well plate precoated with histone substrate	1	+4°C	

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

MATERIALS REQUIRED BUT NOT SUPPLIED:

TBST buffer (1x TBS, pH 8.0, containing 0.05% Tween-20)

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



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: Yang, Y., Bedford, M.T. 2013. Nat Rev Cancer. 13(1):37-50.

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 **20 μ M S-adenosylmethionine** on ice. Upon first thaw, briefly spin tube containing **20 μ M S-adenosylmethionine** to recover full contents of the tube. Aliquot **20 μ M S-adenosylmethionine** into single use aliquots and store at -80°C . *Note: 20 μ M 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	–	–	–	5 μ l
10% DMSO in water (Inhibitor buffer)	5 μ l	5 μ l	5 μ l	–
1X HMT assay buffer 2	20 μ l	–	–	–
Diluted PRMT8 (2.5-5 ng/ μ l)	–	20 μ l	20 μ l	20 μ l
Total	50 μl	50 μl	50 μl	50 μl

- 4) Add 5 μ l of inhibitor solution of each well designated "Test Inhibitor."

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



6405 Mira Mesa Blvd Ste 100

San Diego CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: support@bpsbioscience.com

- 5) For the "Positive Control," "Substrate Control," and "Blank," add 5 µl of 10% DMSO in water (inhibitor buffer).
- 6) Dilute one part **4x HMT assay buffer 2** with three parts water to make **1x HMT assay buffer 2**. Dilute only enough for the assay; store any remaining **4x HMT assay buffer 2** at -20°C.
- 7) Thaw **PRMT8 enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **PRMT8** into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: **PRMT8** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 8) Dilute **PRMT8** in **1x HMT assay buffer 2** at 2.5-5 ng/µl (50-100 ng/20 µl). 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.*
- 9) Add 20 µl of **1x HMT buffer 2** to the wells designated "Blank".
- 10) Initiate reaction by adding 20 µl of diluted **PRMT8** to the wells designated "Positive Control," "Substrate Control," and "Test Sample." Incubate at room temperature for 1 hour.
- 11) Remove the supernatant from the wells and wash the strip three times with 200 µl TBST buffer. Blot dry onto clean paper towels.
- 12) Add 100 µl 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 µl per well. Incubate 1 hour at room temperature with slow shaking.
- 3) Wash plate three times with 200 µl TBST buffer and incubate in **Blocking buffer 4** as in steps 1-11 and 1-12.

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



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 µl per well. Incubate for 30 min. at room temperature with slow shaking.
- 3) Wash plate three times with TBST buffer and incubate in **Blocking buffer 4** as in steps 1-11 and 1-12.
- 4) Just before use, mix on ice 50 µl **ELISA ECL chemiluminescent substrate A** and 50 µl **ELISA chemiluminescent substrate B** and add 100 µl 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 readings.

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.

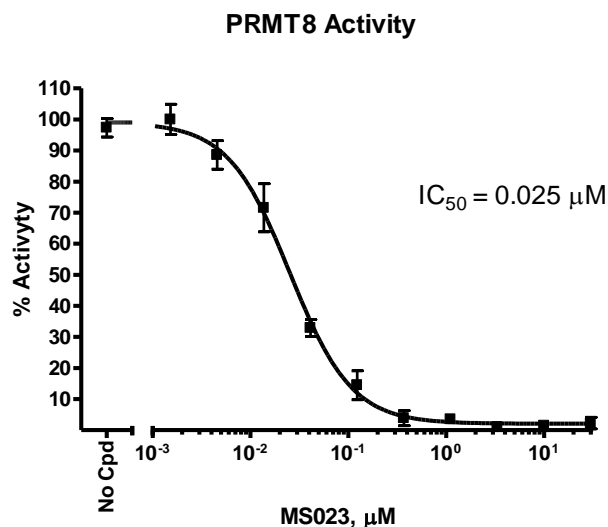
To properly read chemiluminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second, delay after plate movement is 100 msec. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader are: use the "hole" position 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

Example of Assay Results:



PRMT8 enzyme activity, measured using the *PRMT8 Chemiluminescent Assay Kit*, BPS Bioscience Catalog #52062. 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>
PRMT8 (expressed in Sf9)	51052	20 μg
PRMT1 (expressed in Sf9 cells)	51041	20 μg
PRMT3 (expressed in <i>E. coli</i>)	51043	50 μg
PRMT4/CARM1 (expressed in HEK293)	51047	20 μg
PRMT5/MEP50 (expressed in HEK293)	51045	20 μg
PRMT5/MEP50 (expressed in Sf9 cells)	51048	20 μg
PRMT6 (expressed in HEK293)	51049	20 μg
PRMT7 (expressed in Sf9)	51054	20 μg
PRMT9 (expressed in Sf9)	51053	20 μg
PRMT1 Chemiluminescent Assay Kit	52004L	96 reactions
PRMT3 Chemiluminescent Assay Kit	52005L	96 reactions
PRMT4 Chemiluminescent Assay Kit	52041L	96 reactions
PRMT6 Chemiluminescent Assay Kit	52046	96 reactions
PRMT8 Homogeneous Assay Kit	52058	384 reactions
PRMT1 Homogeneous Assay Kit	52054	384 reactions
PRMT3 Homogeneous Assay Kit	52055	384 reactions
PRMT5 Homogeneous Assay Kit	52052	384 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



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	PRMT8 enzyme has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh enzyme (PRMT8, BPS Bioscience #51052). 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	Record light signals at 5 second intervals. Refer to instrument instructions for settings to increase sensitivity of light detection. See "Reading Chemiluminescence" section.
	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 (PRMT8, BPS Bioscience #51052) 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