



6042 Cornerstone Court W, 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**

### ***DNMT1 Direct Activity Assay Kit***

**Catalog # 52050L**

**DESCRIPTION:** The *DNMT1 Direct Activity Assay Kit* is designed to measure DNMT1 activity using purified DNMT1. The *DNMT1 Direct Activity Assay Kit* comes in a convenient format, with a 96-well plate precoated with DNA substrate, the antibody against 5-methylcytosine, the secondary HRP-labeled antibody, S-adenosylmethionine, DNMT1 assay buffer, and purified DNMT1 for 100 enzyme reactions. The key to the *DNMT1 Direct Activity Assay Kit* is a highly specific antibody that recognizes 5-methylcytosine of the substrate. With this kit, only three simple steps on a microtiter plate are required for DNMT1 detection. First, S-adenosylmethionine is incubated with a sample containing assay buffer and DNMT1 for one hour. 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 #	Component	Amount	Storage	
51101	DNMT1	2 x 10 µg	-80°C	<b>(Avoid freeze/thaw cycles!)</b>
52120	400 µM S-adenosylmethionine	2 x 250 µl	-80°C	
	Anti-5-methylcytosine antibody	25 µl	-80°C	
52130H	Secondary HRP-labeled antibody 1	10 µl	-80°C	
52200	4x DNMT assay buffer 1*	5 ml	-20°C	
52100	Blocking buffer 4	50 ml	+4°C	
	HRP chemiluminescent substrate A	6 ml	+4°C	
	HRP chemiluminescent substrate B (brown bottle)	6 ml	+4°C	
	Plate precoated with DNA substrate	1	+4°C	

\*Add 260 µl of 0.5 M DTT before use.

#### **MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

TBST buffer (1 x TBS, pH 8.0, containing 0.05% Tween20)  
Luminometer or fluorescent microplate reader capable of reading chemiluminescence  
Adjustable micropipettor and sterile tips  
Rotating or rocker platform  
0.5M DTT

**APPLICATIONS:** Useful for screening enzyme inhibitors.

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

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

**REFERENCE(S):** Svedruzic ZM. *Curr. Med. Chem.* 2008; **15**(1):92-106

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)

## ASSAY PROTOCOL:

*All samples and controls should be tested in duplicate.*

### Step 1:

- 1) Rehydrate the microwells by adding 150 µl of TBST buffer (1 x 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 content of the tube. Aliquot S-adenosylmethionine into single use aliquots. Store remaining S-adenosylmethionine in aliquots at -80°C immediately. *Note: S-adenosylmethionine is very sensitive to freeze/thaw cycles. Avoid multiple freeze-thaw cycles.*
- 3) Add 260 µl of 0.5 M DTT to 4x DNMT assay buffer 1. Prepare the master mixture: N wells × (7.5 µl **4x DNMT assay buffer 1** + 5 µl **400 µM S-adenosylmethionine** + 12.5 µl **H<sub>2</sub>O** )
- 4) Thaw **DNMT1** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **DNMT1** into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. *Note: DNMT1 enzyme is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 5) Dilute **DNMT1** in **1x DNMT** assay buffer 1 at 10 ng/µl (200 ng/20 µl). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.
- 6) Add 25 µl of master mixture to each well designated for the “Positive Control” and “Test Inhibitor”. For the “Substrate Control”, add 7.5 µl **4x DNMT assay buffer 1** + 17.5 µl **H<sub>2</sub>O**. For the “Blank”, add 12.5 µl **4x DNMT assay buffer 1** + 5 µl **400 µM S-adenosylmethionine** + 27.5 µl **H<sub>2</sub>O**. (Refer to chart)

Component	Positive Control	Test Sample	Substrate Control	Blank
4x DNMT assay buffer 1	7.5 µl	7.5 µl	7.5 µl	12.5 µl
400 µM S-adenosylmethionine	5 µl	5 µl	–	5 µl
H <sub>2</sub> O	12.5 µl	12.5 µl	17.5 µl	27.5 µl
DNMT1 (10 ng/µl)	20 µl	20 µl	20 µl	–
Inhibitor	–	5 µl	–	–
Inhibitor Buffer (no inhibitor)	5 µl	–	5 µl	5 µl
<b>Total</b>	<b>50 µl</b>	<b>50 µl</b>	<b>50 µl</b>	<b>50 µl</b>

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 Court W, 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) Add 5  $\mu$ l of inhibitor solution of each well designated "Test Inhibitor". For the "Positive Control", "Substrate Control" and "Blank", add 5  $\mu$ l of the same solution without inhibitor (inhibitor buffer).
- 8) Initiate reaction by adding 20  $\mu$ l of diluted **DNMT1** (prepared as described above). Incubate at 37°C for 1 hour.
- 9) Remove the supernatant from the wells and wash the plate three times with 200  $\mu$ l TBST buffer. Blot dry onto clean paper towels.
- 10) Add 100  $\mu$ l of **Blocking buffer 4** to every well. Shake on a rotating platform for 10 min. Remove supernatant as described above.

#### Step 2:

- 1) Dilute "**Anti-5-methylcytosine antibody**" 400-fold with **Blocking buffer 4**.
- 2) Add 100  $\mu$ l per well. Incubate 1 hour at room temperature with slow shaking.
- 3) Remove the supernatant from the wells and wash plate three times with 200  $\mu$ l TBST buffer and incubate in **Blocking buffer 4** as described in steps 1-9 and 1-10.

#### Step 3:

- 1) Dilute "**Secondary HRP-labeled antibody 1**" 1,000-fold with **Blocking buffer 4**.
- 2) Add 100  $\mu$ l per well. Incubate for 30 min. at room temperature with slow shaking.
- 3) Remove the supernatant from the wells and wash plate three times with TBST buffer and incubate in **Blocking buffer 4** as described in steps 1-9 and 1-10.
- 4) Just before use, mix on ice 50  $\mu$ l **HRP chemiluminescent substrate A** and 50  $\mu$ l **HRP chemiluminescent substrate B** and add 100  $\mu$ l per well. Discard any unused chemiluminescent reagent after use.
- 5) Immediately read sample in a luminometer or microtiter-plate reader 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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



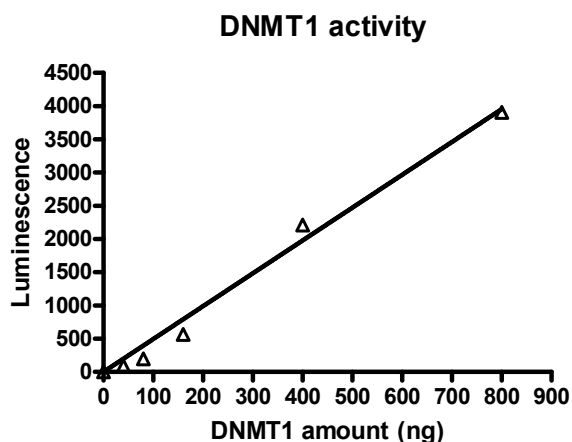
6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

### 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 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).

### Example of Assay Results:



DNMT1 enzyme activity, measured using the DNMT1 Assay Kit, BPS Bioscience #52050. 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](mailto: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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6042 Cornerstone Court W, Ste B  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

#### **RELATED PRODUCTS**

<b><u>Product Name</u></b>	<b><u>Catalog #</u></b>	<b><u>Size</u></b>
DNMT1	51101	10 µg
DNMT2	51102	10 µg
DNMT3a	51103	10 µg
4x DNMT Assay Buffer 1	52200	30 ml
EZH2 Assay Kit	52009	96 reactions
G9a Assay Kit	52001	96 reactions
MLL Assay Kit	52008	96 reactions
PRMT1 Assay Kit	52004	96 reactions
PRMT3 Assay Kit	52005	96 reactions
PRMT5 Assay Kit	52002	96 reactions
SET7/9 Assay Kit	52003	96 reactions
SUV39H1 Assay Kit	52006	96 reactions
SUV39H2 Assay Kit	52007	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: [\*\*info@bpsbioscience.com\*\*](mailto:info@bpsbioscience.com)

Please visit our website at: [\*\*www.bpsbioscience.com\*\*](http://www.bpsbioscience.com)

## TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Luminescence signal of positive control reaction is weak	DNMT1 enzyme has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh enzyme (DNMT1, BPS Bioscience #51101). 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. See section on "Reading Chemiluminescence" above.
	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	Be sure to include blocking steps after wash steps. 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 (DNMT1, BPS Bioscience #51101) 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: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)