



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

SetDB1 Chemiluminescent Assay Kit

Catalog # 51056
Size: 96 reactions

DESCRIPTION: The *SetDB1 Chemiluminescent Assay Kit* is designed to measure SetDB1 activity for screening and profiling applications. The *SetDB1 Chemiluminescent Assay Kit* comes in a convenient format, with 8-well strips precoated with histone H3 peptide substrate, an antibody against methylated lysine residue of Histone H3, a secondary HRP-labeled antibody, S-adenosylmethionine, methyltransferase assay buffer, and purified SetDB1 enzyme for 96 enzyme reactions. The key to the *SetDB1 Direct Activity Assay Kit* is a highly specific antibody that recognizes the methylated K9 residue of Histone H3. With this kit, only three simple steps 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 strips 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	
51055	SetDB1 human enzyme*	5 µg	-80°C	(Avoid freeze/thaw cycles!)
52120	20 µM S-adenosylmethionine	250 µl	-80°C	
52140A	Primary antibody 1	100 µl	-80°C	
52130H	Secondary HRP-labeled antibody 1	10 µl	-80°C	
52160	4x HMT assay buffer 1**	3 ml	-20°C	
52100	Blocking buffer 4	50 ml	+4°C	
	HRP chemiluminescent substrate A (transparent bottle)	6 ml	+4°C	
	HRP chemiluminescent substrate B (brown bottle)	6 ml	+4°C	
	8-well strip plate precoated with histone substrate	1 plate (12 x 8-well strips)	+4°C	

*The concentration of SetDB1 is lot-specific and will be indicated on the tube containing the enzyme.

**Add 125 µl of 0.5M DTT before use.

MATERIALS REQUIRED BUT NOT SUPPLIED:

TBST buffer (1 x TBS, pH 8.0, containing 0.05% Tween-20)
Luminometer or fluorescent microplate reader capable of reading chemiluminescence
Adjustable micropipettor and sterile tips
Rotating or rocker platform

APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

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

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 strip 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. 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 125 μ l of 0.5 M DTT to **4x HMT assay buffer 1** before use.
- 4) Prepare the master mixture: N wells \times (7.5 μ l **4X HMT assay buffer 1** + 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 1** + 17.5 μ l water.

	Blank	Substrate Control	Positive Control	Test Sample
4X HMT assay buffer 1	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
Inhibitor buffer (no inhibitor)	5 μ l	5 μ l	5 μ l	-
1X HMT assay buffer 1	20 μ l	-	-	-
Diluted SetDB1 (1.25-2.5 ng/ μ l)		20 μ l	20 μ l	20 μ l
Total	50 μl	50 μl	50 μl	50 μl

- 5) 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
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) For "Positive Control", "Substrate Control" and "Blank", add 5 μ l of the same solution without inhibitor (inhibitor buffer).
- 7) Thaw **SetDB1 enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot **SetDB1 enzyme** into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: **SetDB1 enzyme** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 8) Dilute **4x HMT assay buffer 1** with water to prepare **1x HMT assay buffer 1**. Prepare only the amount required for the assay.
- 9) Dilute **SetDB1 enzyme** in **1x HMT assay buffer 1** to 1.25 – 2.5 ng/ μ l (25-50 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.***
- 10) Add 20 μ l of **1x HMT buffer 1** to the well designated "Blank".
- 11) Initiate reaction by adding 20 μ l of diluted **SetDB1 enzyme** to the wells designated "Positive Control", "Substrate Control", and "Test Sample". Incubate at room temperature for 1 hour.
- 12) Remove the supernatant from the wells and wash the strip three times with 200 μ l TBST buffer. Blot dry onto clean paper towels.
- 13) Add 100 μ l of **Blocking buffer 4** to every well. Shake on a rotating platform for 10 min. Blot dry onto clean paper towels.

Step 2:

- 1) Dilute **Primary antibody 1** 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. Add 100 μ l of **Blocking buffer 4** to every well and incubate in **Blocking buffer 4** on a rotating platform for 10 min. Blot dry onto clean paper towels.

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 1** 1,000-fold with **Blocking buffer 4**. Add 100 μ l per well. Incubate for 30 minutes at room temperature with slow shaking.
- 2) Wash strip three times with 200 μ l TBST buffer. Add 100 μ l of **Blocking buffer 4** to every well and incubate in **Blocking buffer 4** on a rotating platform for 10 min. Blot dry onto clean paper towels.
- 3) 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.
- 4) Immediately read sample in a luminometer or microtiter-plate reader capable of reading chemiluminescence. "Blank" value is subtracted from all readings.

Reading Chemiluminescence:

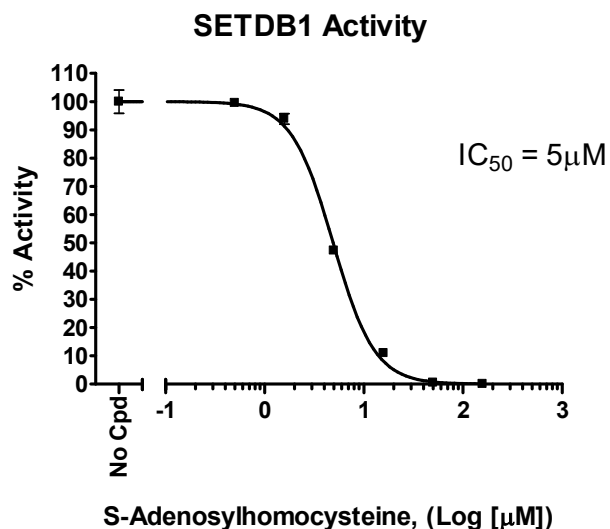
Chemiluminescence is the emission of light (luminescence) which results from a chemical reaction. The detection of chemiluminescence requires no wavelength selection because the method used is emission photometry and is not emission spectrophotometry.

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
Please visit our website at: www.bpsbioscience.com

Example of Assay Results:



SetDB1 enzyme activity, measured using the *SetDB1 Chemiluminescent Assay Kit*, BPS Bioscience Catalog #51056 in the presence of S-Adenosylhomocysteine. 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 support@bpsbioscience.com.*

RELATED PRODUCTS

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
SetDB1 enzyme	51055	50 µg
SetD2 enzyme	53019	50 µg
Set7/Set9 enzyme	51010	100 µg
Set8 enzyme	51008	50 µg
G9a enzyme (<i>E. coli</i>)	51000	50 µg
G9a enzyme (Sf9 cells)	51001	20 µg
G9a Homogeneous Assay Kit	52051	384 reactions
SETD2 Chemiluminescent Assay Kit	52060	96 reactions
H3(K9) Universal Methyltransferase Assay Kit	52072	96 reactions
SUV39H1 (82-end) enzyme	51070	50 µg
SUV39H1 (full length) enzyme	51071	5 µg
SUV39H2 enzyme	51080	50 µg
SUV39H1 Chemiluminescent Assay Kit	52045	96 reactions
SUV39H2 Chemiluminescent Assay Kit	52008	96 reactions
Anti-SETD1a Polyclonal Antibody	25314	100 µl
Anti-SETD1b Polyclonal Antibody	25313	100 µl
Anti-SETD8 Polyclonal Antibody	25315	100 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

TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Luminescence signal of positive control reaction is weak	SetDB1 enzyme has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh enzyme (SetDB1, BPS Bioscience #51055). Store enzyme in single-use aliquots. Increase time of enzyme incubation. Increase enzyme concentration.
	Antibody reaction is insufficient	Increase time for antibody incubation. Avoid freeze/thaw cycles of antibody.
	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 strip 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 (SetDB1, BPS Bioscience 51055) 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