



6405 Mira Mesa Blvd Ste 100

San Diego CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

**Data Sheet**  
***NSD2 Chemiluminescent Assay Kit***  
**Catalog # 53009**  
**Size: 96 reactions**

**DESCRIPTION:** The *NSD2 Chemiluminescent Assay Kit* is designed to measure NSD2 activity for screening and profiling applications. The *NSD2 Chemiluminescent Assay Kit* comes in a convenient format, with 96-well plate precoated with histone H3 substrate, the antibody against methylated lysine residue of Histone H3, a secondary HRP-labeled antibody, S-adenosylmethionine, methyltransferase assay buffer, and purified NSD2 enzyme for 96 enzyme reactions. The key to the *NSD2 Chemiluminescent Activity Assay Kit* is a highly specific antibody that recognizes methylated 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 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	
51026	NSD2 human enzyme	50 µg	-80°C	<b>(Avoid freeze/thaw cycles!)</b>
52120	100 µM S-adenosylmethionine	250 µl	-80°C	
52140P2	Primary antibody 16-2	12.5 µl	-80°C	
52131H	Secondary HRP-labeled antibody 2	10 µl	-80°C	
52193Z	4x HMT assay buffer 7	3 ml	-20°C	
79556	Blocking buffer 1	50 ml	+4°C	
	HRP chemiluminescent substrate A (translucent bottle)	6 ml	+4°C	
	HRP chemiluminescent substrate B (brown bottle)	6 ml	+4°C	
	White microplate precoated with histone substrate	1	+4°C	

**Note:**

Since May of 2018, this kit has been improved with reformulated buffers. The previous version of the NSD2 kit #53009 can still be purchased upon special request.

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](mailto:support@bpsbioscience.com)

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

**MATERIALS REQUIRED BUT NOT SUPPLIED:**

TBST buffer (1 x Tris-buffered saline (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

**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(S):**

1. 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  $\mu$ 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 **100  $\mu$ M 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: 100  $\mu$ M S-adenosylmethionine is sensitive to freeze/thaw cycles. Avoid multiple freeze-thaw cycles.*
- 3) Prepare the master mixture: N wells  $\times$  (7.5  $\mu$ l **4x HMT assay buffer 7** + 2.5  $\mu$ l **100  $\mu$ M S-adenosylmethionine** + 15  $\mu$ l **H<sub>2</sub>O**)
- 4) Add 25  $\mu$ l of master mixture to each well designated for the "Positive Control," "Test Inhibitor," and "Blank." For the "Substrate Control," add 7.5  $\mu$ l **4x HMT assay buffer 7** + 17.5  $\mu$ l **H<sub>2</sub>O**.

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](mailto:support@bpsbioscience.com)

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

- 5) Thaw **NSD2 human enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **NSD2 human enzyme** into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: **NSD2 human enzyme** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 6) Dilute **NSD2 human enzyme** in **1x HMT assay buffer 7** to 10-25 ng/μl (200-500 ng/20 μl). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.

	Blank	Substrate Control	Positive Control	Test Inhibitor
4x HMT assay buffer 7	7.5 μl	7.5 μl	7.5 μl	7.5 μl
100 μM S-adenosylmethionine	2.5 μl	–	2.5 μl	2.5 μl
H <sub>2</sub> O	15 μl	17.5 μl	15 μl	15 μl
Test Inhibitor/Activator	–	–	–	5 μl
Inhibitor buffer (no inhibitor)	5 μl	5 μl	5 μl	–
1x HMT assay buffer 7	20 μl	–	–	–
NSD2 (10-25 ng/μl)	–	20 μl	20 μl	20 μl
<b>Total</b>	<b>50 μl</b>	<b>50 μl</b>	<b>50 μl</b>	<b>50 μl</b>

- 7) Add 5 μl of inhibitor solution of each well designated “Test Inhibitor.” For the “Positive Control,” “Substrate Control,” and “Blank,” add 5 μl of the same solution without inhibitor (**inhibitor buffer**). *Note: The NSD2 Chemiluminescent Assay Kit is compatible with up to 1% final DMSO concentration. We recommend preparing the inhibitor in 10% DMSO aqueous solution and using 5 μl per NSD2 reaction.*
- 8) Add 20 μl of **1 x HMT buffer 7** to the well designated “Blank.”
- 9) Initiate reaction by adding 20 μl of diluted **NSD2 Human Enzyme** (prepared as described above) to the wells labeled “Test Inhibitor,” “Positive Control,” and “Substrate Control.” Incubate two hours at room temperature on a rotating platform.
- 10) Wash the wells three times with 200 μl TBST buffer. Blot dry onto clean paper towels.
- 11) Add 100 μl of **Blocking buffer 1** to every well. Shake on a rotating platform for 10 min. Remove supernatant as above.

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](mailto:support@bpsbioscience.com)

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

### Step 2:

- 1) Dilute **Primary antibody 16-2** 800-fold with **Blocking buffer 1**.
- 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 1** as in steps 1-10 and 1-11.

### Step 3:

- 1) Dilute **Secondary HRP-labeled antibody 2** 1,000-fold with **Blocking buffer 1**.
- 2) Add 100 µl per well. Incubate for 30 minutes at room temperature with slow shaking.
- 3) Wash plate three times with TBST buffer and incubate in **Blocking buffer 1** as in steps 1-10 and 1-11.
- 4) Just before use, mix on ice 50 µl **HRP chemiluminescent substrate A** and 50 µl **HRP 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 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, and 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](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd Ste 100

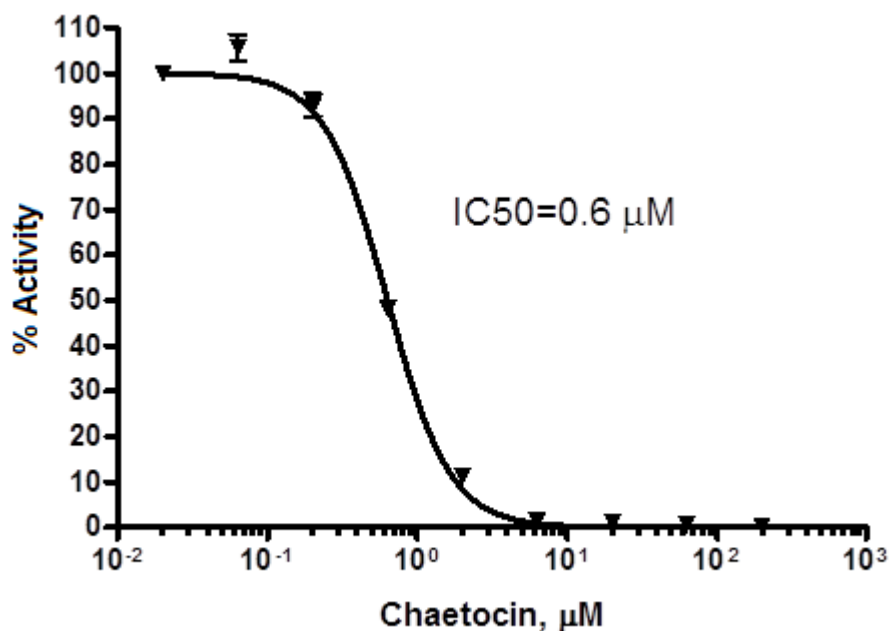
San Diego CA 92121

Tel: 1.858.202.1401

Fax: 1.858.481.8694

Email: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

#### Example of Assay Results:



NSD2 enzyme activity, measured using the *NSD2 Chemiluminescent Assay Kit*, BPS Bioscience Catalog #53009. 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).

#### RELATED PRODUCTS

Product	Catalog #	Size
NSD2 enzyme (catalytic)	#51026	50 $\mu$ g
NSD2 (782-end)/ReBPII enzyme	#51025	20 $\mu$ g
NSD1 enzyme	#51024	50 $\mu$ g
NSD3 (1021-1322) enzyme	#51036	50 $\mu$ g
SETD2 enzyme	#53019	50 $\mu$ g
SETD2 Chemiluminescent Assay Kit	#52060	96 rxns
Chaetocin	#27221	1 mg
4x HMT Assay Buffer 7	#52193B	30 mL
FBXL10(KDM2B, JHDM1B) enzyme	#50120	20 $\mu$ g
FBXL11(KDM2A) enzyme	#50102	20 $\mu$ g

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](mailto:support@bpsbioscience.com)

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

## TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Luminescence signal of positive control reaction is weak	NSD2 enzyme has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh enzyme (NSD2, BPS Bioscience #51026). 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 (NSD2, BPS Bioscience #51026) 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](mailto:support@bpsbioscience.com)

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