



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet
EZH2 Chemiluminescent Assay Kit
Catalog # 52085
Size: 384 reactions

DESCRIPTION: The *f WT (Wild Type) Chemiluminescent Assay Kit* is designed to measure activity of the EZH2 complex (EZH2/EED/SUZ12/RbAp48/ AEBP) for screening and profiling purposes. The *EZH2 Chemiluminescent Assay Kit* comes in a convenient format, with wells precoated with histone H3 peptide substrate, an antibody against methylated K27 residue of Histone H3, the secondary HRP-labeled antibody, S-adenosylmethionine, methyltransferase assay buffer, and purified EZH2 WT complex for 384 enzyme reactions. The key to the EZH2 Assay Kit is a highly specific antibody that recognizes methylated Histone H3K27. 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 for one hour. 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	
51004	EZH2/EED/SUZ12/RbAp48/AEBP2	4 x 25 µg	-80°C	Avoid freeze/ thaw cycles!
52120	100 µM S-adenosylmethionine	4 x 250 µl	-80°C	
52140F	Primary Antibody 6	2 x 12.5 µl	-80°C	
52131H	Secondary HRP-labeled Antibody 2	2 x 10 µl	-80°C	
52170	4x HMT Assay buffer 2	2 x 3 ml	-20°C	
79556	Blocking buffer 1	2 x 50 ml	+4°C	
	HRP chemiluminescent substrate A (translucent bottle)	2 x 6 ml	+4°C	
	HRP chemiluminescent substrate B (brown bottle)	2 x 6 ml	+4°C	
	384-well plate precoated with histone substrate	1 plate	+4°C	

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

TBST buffer (1x Tris-buffered saline, pH 8.0, containing 0.05% Tween-20)
Luminometer or fluorescent microplate reader capable of reading chemiluminescence
Rotating or rocker platform

APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

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

REFERENCE(S):

1. Dillon, S.C., *et al. Genome Biology* 2005; **6**:227.
2. Morin, R.D., *et al. Nat Genet.* 2010, **42**(2):181.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Step 1:

- 1) Rehydrate the microwells by adding 90 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 **100 μ M S-adenosylmethionine** on ice. Upon first thaw, briefly spin tube containing **100 μ M S-adenosylmethionine** to recover full content of the tube. Aliquot **100 μ M S-adenosyl-methionine** into single use aliquots. Store remaining **100 μ M S-adenosylmethionine** in aliquots at -80°C immediately. *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 **4 \times HMT Assay Buffer 2** + 2.5 μ l **100 μ M S-adenosylmethionine** + 15 μ l **H₂O**)

	Blank	Substrate Control	Positive Control	Test Inhibitor
4 \times HMT assay Buffer 2	7.5 μ l	7.5 μ l	7.5 μ l	7.5 μ l
100 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	-
1 \times HMT assay Buffer 2	20 μ l	-	-	-
EZH2 (12.5 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 4) Add 25 μ l of master mixture to each well designated for the "Positive Control", "Test Inhibitor", and "Blank". For the "Substrate Control", add 7.5 μ l **4 \times HMT Assay Buffer 2** + 17.5 μ l **H₂O**
- 5) Add 5 μ l of inhibitor solution of each well designated "Test Inhibitor". For the "Positive Control", "Substrate Control" and "Blank", add 5 μ l of the same solution without inhibitor (inhibitor buffer).
- 6) Add 20 μ l of **1 \times HMT assay Buffer 2** to the well designated "Blank".
- 7) Thaw **EZH2 enzyme** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **EZH2 enzyme** into single use aliquots. Store remaining undiluted enzyme in aliquots at -70°C immediately. *Note: **EZH2 enzyme** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 8) Dilute **EZH2 enzyme** in **1 \times HMT assay Buffer 2** at 12.5 ng/ μ l (250 ng/20 μ l). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.
- 9) Initiate reaction by adding 20 μ l of diluted **EZH2 enzyme** prepared as described above. Incubate at room temperature for one hour.
- 10) Wash the strips three times with 90 μ l TBST buffer. Blot dry onto clean paper towels.
- 11) Add 50 μ l of **Blocking Buffer 1** to every well. Shake on a rotating platform for 10 min. Remove supernatant as described above.

Step 2:

- 1) Dilute **Primary Antibody 6** 800-fold with **Blocking Buffer 1**.
- 2) Add 50 μ l per well. Incubate 1 hour at room temperature with slow shaking.
- 3) Wash the strips three times with TBST buffer and incubate in **Blocking Buffer 1** as described 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 50 μ l per well. Incubate for 30 min. at room temperature with slow shaking.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 3) Wash the strips with TBST buffer and incubate in **Blocking Buffer 1** as described in step 1-10 and 1-11.
- 4) Just before use, mix on ice 25 μ l **HRP chemiluminescent substrate A** and 25 μ l **HRP chemiluminescent substrate B** and add 50 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 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 **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: [**info@bpsbioscience.com**](mailto:info@bpsbioscience.com)

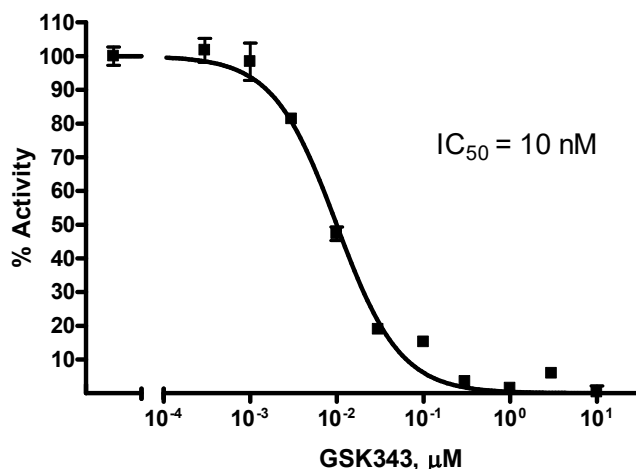
Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd Ste 100
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Example of Assay Results:

EZH2 Activity



EZH2 WT enzyme activity, measured using the *EZH2 Chemiluminescent Assay Kit*, BPS Bioscience #52085. 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 Name</u>	<u>Catalog #</u>	<u>Size</u>
EZH2 (Y641F)/EED/SUZ12/RbAp48/AEBP2	51017	20 µg
EZH2 (Y641C)/EED/SUZ12/RbAp48/AEBP2	51029	20 µg
EZH2 (Y641N)/EED/SUZ12/RbAp48/AEBP2	51028	20 µg
EZH2 (Y641S)/EED/SUZ12/RbAp48/AEBP2	51013	20 µg
EZH2 (Y641H)/EED/SUZ12/RbAp48/AEBP2	51011	20 µg
EZH2/EED/SUZ12/RbAp48/AEBP2	51004	50 µg
EZH2 Chemiluminescent Assay Kit	52009L	96 rxns.
EZH2 (Y641F) Chemiluminescent Assay Kit	52075	96 rxns.
EZH2 (Y641N) Chemiluminescent Assay Kit	52076	96 rxns.
EZH2 WT Chemiluminescent Assay Kit	52067	96 rxns.
EZH2 Homogeneous Assay Kit	52059	384 rxns.
EZH2 (Y641N) TR-FRET Assay Kit	52078	384 rxns.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6405 Mira Mesa Blvd Ste 100
 San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Luminescence signal of positive control reaction is weak	EZH2 Complex has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh EZH2 WT Complex, BPS Bioscience #51004. 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 EZH2 Complex, BPS #51004 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com