



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet ***EZH2 Assay Kit*** **Catalog # 52009L**

DESCRIPTION: The *EZH2/EED/SUZ12/RbAp48/AEBP2 Direct Activity Assay Kit (EZH2 Assay Kit)* is designed to measure activity of the EZH2 complex (EZH2/EED/SUZ12/RbAp48/AEBP) for screening and profiling purposes. The *EZH2 Assay Kit* comes in a convenient format, with a 96-well plate 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 complex for 100 enzyme reactions. The key to the EZH2 Direct Activity 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 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 | |
|------------------|---|---------------|----------------|------------------------------------|
| 51004 | EZH2/EED/SUZ12/RbAp48/AEBP2 | 50 µg | -80 °C | (Avoid freeze/thaw cycles!) |
| 52120 | 400 µM S-adenosylmethionine | 250 µl | -80 °C | |
| 52140C | Primary antibody 3 | 100 µl | -80 °C | |
| 52130H | Secondary HRP-labeled antibody 1 | 10 µl | -80 °C | |
| 52170 | 4x HMT assay buffer 2 | 3 ml | -20 °C | |
| 52100 | Blocking buffer | 50 ml | +4 °C | |
| | HRP chemiluminescent substrate (2 components) | 6 ml each | +4 °C | |
| | Black microplate precoated with histone substrate | 1 | +4 °C | |

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
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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

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

REFERENCES:

Dillon SC, Zhang X, Trievel RC, Cheng X. Genome Biology 2005; **6**:227.
Morin, RD., 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 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 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) Prepare the master mixture: N wells \times (7.5 μ l **4 \times HMT assay buffer 2** + 1.25 μ l **400 μ M S-adenosylmethionine** + 16.25 μ l **H₂O**)
- 4) 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 -80°C immediately. *Note: EZH2 enzyme is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 5) Dilute **EZH2 enzyme** in 1 \times HMT assay buffer 2 at 25 ng/ μ l (500 ng/20 μ l). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.

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



6044 Cornerstone Court W, Ste E
 San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

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 **4x HMT assay buffer 2** + 17.5 μ l **H₂O**

| | Blank | Substrate Control | Positive Control | Test Inhibitor |
|----------------------------------|-----------------------------|-----------------------------|-----------------------------|-----------------------------|
| 4x HMT assay buffer 2 | 7.5 μ l | 7.5 μ l | 7.5 μ l | 7.5 μ l |
| 400 μ M S-adenosylmethionine | 1.25 μ l | – | 1.25 μ l | 1.25 μ l |
| H ₂ O | 16.25 μ l | 17.5 μ l | 16.25 μ l | 16.25 μ l |
| Test Inhibitor/Activator | – | – | – | 5 μ l |
| Inhibitor buffer (no inhibitor) | 5 μ l | 5 μ l | 5 μ l | – |
| 1x HMT assay buffer 2 | 20 μ l | – | – | – |
| EZH2 (25 ng/ μ l) | – | 20 μ l | 20 μ l | 20 μ l |
| Total | 50 μl | 50 μl | 50 μl | 50 μl |

- 6) 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).
- 7) Add 20 μ l of 1 x HMT assay buffer 2 to the well designated “Blank”.
- 8) Initiate reaction by adding 20 μ l of diluted EZH2 prepared as described above. Incubate at room temperature for one hour.
- 9) Wash the plate three times with 200 μ l TBST buffer. Blot dry onto clean paper towels.
- 10) Add 100 μ l of Blocking buffer 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Step 2:

- 1) Dilute "Primary antibody 3" 100-fold with Blocking buffer.
- 2) Add 100 μ l per well. Incubate 1 hour at room temperature with slow shaking.
- 3) Wash plate three times with TBST buffer and incubate in Blocking buffer as in steps 1-9 and 1-10.

Step 3:

- 1) Dilute "Secondary HRP-labeled antibody 1" 1,000-fold with Blocking buffer.
- 2) Add 100 μ l per well. Incubate for 30 min. at room temperature with slow shaking.
- 3) Wash plate with TBST buffer and incubate in Blocking buffer as in step 1-9 and 1-10.
- 4) 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.

Step 4:

Immediately read sample in a luminometer or microtiter-plate 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 **1.858.829.3082** Fax **1.858.481.8694**

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

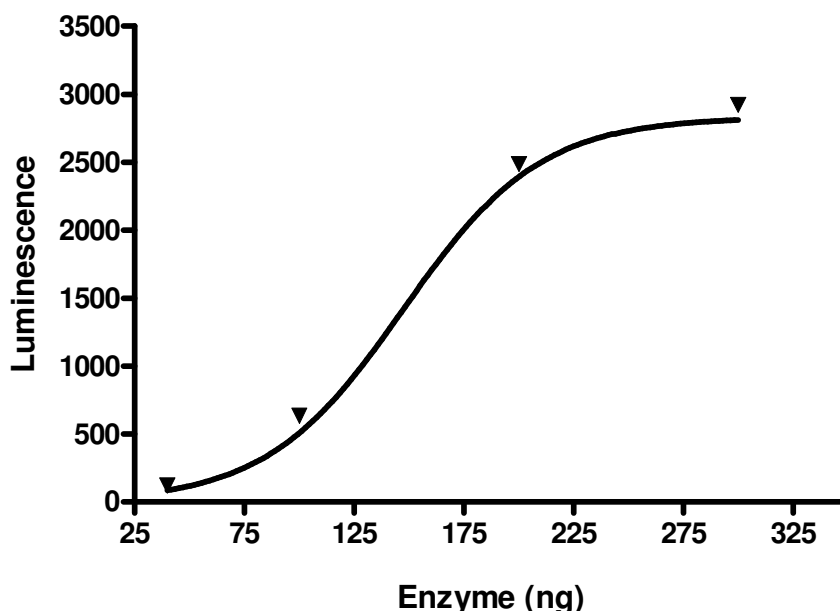
Please visit our website at: [**www.bpsbioscience.com**](http://www.bpsbioscience.com)



6044 Cornerstone Court W, Ste E
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/EED/SUZ12/RbAp48/AEBP2 enzyme activity, measured using the EZH2 Direct Assay Kit, BPS Bioscience # 52009L. 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 | CATALOG # | SIZE |
|-----------------------------|------------------|--------------|
| EZH1/EED/SUZ12 | 51006 | 50 µg |
| EZH1/EED/SUZ12/RbAp48/AEBP2 | 51007 | 50 µg |
| EZH2/EED inactive | 51002 | 20 µg |
| EZH2/EED/SUZ12 | 51003 | 50 µg |
| EZH2/EED/SUZ12/RbAp48/AEBP2 | 51004 | 50 µg |
| 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 **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
 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 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. |
| | 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 EZH2 Complex, BPS Bioscience #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