



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet **JMJD1B Homogeneous Assay Kit** Catalog #79839

DESCRIPTION: The *JMJD1B Homogeneous Assay Kit* is designed to measure JMJD1B activity for screening and profiling applications. JMJD1B, also known as JHDM2B and KDM3B, is a JumonjiC (JmjC) domain containing histone lysine demethylase that exhibits demethylation activity toward H3-K₉Me² and H3-K₉Me¹. The *JMJD1B Homogeneous Assay Kit* comes in a convenient AlphaLISA[®] format (Scheme 1), with biotinylated histone H3 peptide substrate, primary antibody, demethylase assay buffer, and purified JMJD1B for 384 enzyme reactions. The key to the *JMJD1B Homogeneous Assay Kit* is a highly specific antibody that recognizes demethylated substrate. With this kit, only three simple steps on a microtiter plate are required for methyltransferase detection. First, a sample containing JMJD1B enzyme is incubated with the biotinylated substrate. Next, acceptor beads and primary antibody are added, then donor beads, followed by reading the Alpha-counts.

COMPONENTS:

Catalog #	Component	Amount	Storage	
50421	JMJD1B (KDM3B)	60 µg	-80°C	(Avoid freeze/thaw cycles!)
52140Q4	Primary antibody 17-4	20 µl	-80°C	
79845	Biotinylated histone H3 peptide substrate	500 rxns	-80°C	
52407	4x HDM Assay Buffer 2	3 ml	-80°C	
52301	4x Detection buffer	2 ml	-20°C	

MATERIALS REQUIRED BUT NOT SUPPLIED:

AlphaLISA[®] anti-RlgG acceptor beads, 5 mg/ml (PerkinElmer #AL104C)
AlphaScreen[®] Streptavidin-conjugated donor beads, 5 mg/ml (PerkinElmer #6760002S)
Optiplate -384 (PerkinElmer #6007290)
AlphaScreen[®] microplate reader
Adjustable micropipettor and sterile tips

APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

SAFETY: This product is for research purposes only and not for human or therapeutic use. This product should be considered hazardous. Do not ingest, inhale, get in eyes, on skin, or on clothing. If so, wash thoroughly.

STABILITY: At least one year from date of receipt when stored as directed.

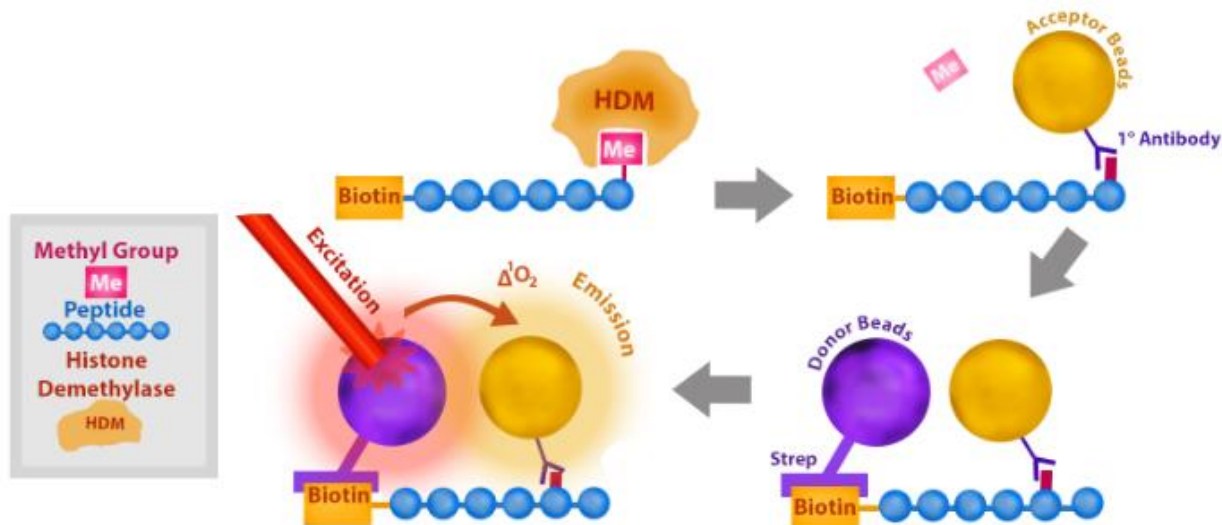
REFERENCE:

Horton, J.R., *et al. Nature Structural & Molecular Biology* 2010; **17(1)**: 38-43.

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
Please visit our website at: www.bpsbioscience.com

Scheme 1: Our histone demethylase assays utilize highly specific antibodies that recognize demethylated products. First, a sample containing the enzyme is incubated with a biotinylated substrate. Next, acceptor beads and primary antibody are added, then donor beads, followed by reading the Alpha-counts, as shown below.



CONTRAINDICATIONS: Green and blue dyes that absorb light in the AlphaScreen signal emission range (520-620 nm), such as Trypan Blue. Avoid the use of the potent singlet oxygen quenchers such as sodium azide (NaN_3) or metal ions (Fe^{2+} , Fe^{3+} , Cu^{2+} , Zn^{2+} and Ni^{2+}). The presence of >1% RPMI 1640 culture medium leads to a signal reduction due to the presence of excess biotin and iron in this medium. MEM, which lacks these components, does not affect AlphaScreen assays.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate. We recommend preincubating the enzyme with inhibitor, however, it is acceptable to add the substrate mixture and inhibitor followed by diluted JMJD1B without the preincubation step.

Step 1:

- 1) Re-suspend lyophilized **Biotinylated histone H3 peptide substrate** in 500 μl of distilled water.
- 2) Prepare serial dilutions of the test inhibitors in **1x HDM Assay Buffer 2**. Note: Final DMSO concentration in the assay should be $\leq 1\%$. If test inhibitor is dissolved in DMSO, make a 100x stock solution and dilute 1:30 in **1x HDM Assay Buffer 2** (DMSO)

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

Please visit our website at: 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

concentration will be 3.3%). If serial dilutions are needed, use **1x HDM Assay Buffer 2** with 3.3% DMSO to maintain the same DMSO concentration in all samples.

- 3) Add 3 μ l of inhibitor solution to each well designated "Test Sample". For the wells designated "Blank" and "Positive Control" add 3 μ l of the same solution without inhibitor (typically **1x HDM Assay Buffer 2** with respective concentration of DMSO).
- 4) Thaw **JMJD1B** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **JMJD1B** enzyme into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: **JMJD1B** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 5) Dilute **JMJD1B** in **1x HDM Assay Buffer 2** at 25 ng/ μ l (100 ng/4 μ l). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.
- 6) Preincubate 4 μ l of diluted **JMJD1B** with 3 μ l of diluted inhibitor(s) for up to 30 minutes at room temperature, with slow shaking. For the wells designated as "Blank", add 4 μ l **1x HDM Assay Buffer 2**.
- 7) Prepare master mix: N wells \times (1.5 μ l **4x HDM Assay Buffer 2** + 1 μ l **Biotinylated substrate** + 0.5 μ l **distilled water**).
- 8) Initiate reaction by adding 3 μ l of master mix prepared as described above. Incubate at room temperature for one hour. *Note: All incubations are done with slow shaking on a rotator platform.*

Reagent	Blank	Positive Control	Test Inhibitor
Test Inhibitor/Activator	–	–	3 μ l
1x HDM Assay Buffer 2 (3.3% DMSO)	3 μ l	3 μ l	–
JMJD1B (25 ng/ μ l)	–	4 μ l	4 μ l
1x HDM Assay Buffer 2	4 μ l	–	–
4x HDM Assay Buffer 2	1.5 μ l	1.5 μ l	1.5 μ l
Biotinylated Substrate	1 μ l	1 μ l	1 μ l
Distilled water	0.5 μ l	0.5 μ l	0.5 μ l
Total	10 μ l	10 μ l	10 μ l

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
Please visit our website at: www.bpsbioscience.com

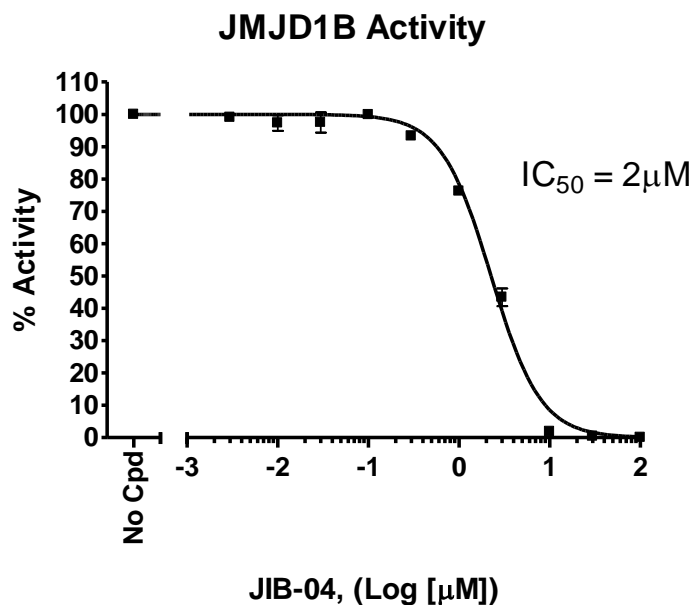
Step 2:

Note: Protect your samples from direct exposure to light!

- 1) Dilute anti-Rabbit Acceptor beads (PerkinElmer #AL104C) (1:500) and Primary antibody 17-4 (1:200) with 1x Detection buffer in one step. Add 10 μ l of acceptor beads/antibody mixture per well. Incubate 30 min at room temperature.

Step 3:

- 1) Dilute **Streptavidin-conjugated donor beads** (PE #6760002S) 125-fold with **1x Detection buffer**. Add 10 μ l of donor beads per well. Shake on a rotator platform for 30 minutes at room temperature.
- 2) Read Alpha-counts.

Example of Assay Results:

Inhibition of JMJD1B enzyme activity by JIB-04, measured using the *JMJD1B Homogeneous Assay Kit*, BPS Bioscience #79839. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at 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
Please visit our website at: 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

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
JMJD1A Assay Kit, Homogeneous	50412-2	384 reactions
JHDM1D Assay Kit, Homogeneous	50420-2	384 reactions
PHF8 Assay Kit, Homogeneous	50515-2	384 reactions
JMJD1A Assay Kit, Homogeneous	50420-1	96 reactions
PHF8 Assay Kit, Homogeneous	50515-1	96 reactions
JHDM1D Assay Kit, Homogeneous	50420-1	96 reactions
JHDM1D Assay Kit, Chemiluminescence	50612	96 reactions
JMJD1A recombinant protein	50130	20 µg
JMJD1B recombinant protein	50421	20 µg
JHDM1D recombinant protein	50419	20 µg
PHF8 recombinant protein	50131	20 µg

AlphaScreen® and AlphaLISA® are registered trademarks of PerkinElmer, Inc.

TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Alpha-counts signal of positive control reaction is same as “blank” value.	JMJD1B has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh JMJD1B, BPS Bioscience #50421. Store enzyme in single-use aliquots. Increase time of enzyme incubation. Increase enzyme concentration.
	Streptavidin Donor beads or anti-rIgG acceptor beads fail to show significant signal.	Reorder Streptavidin Donor beads or anti-rIgG acceptor beads from Perkin Elmer.
	Incorrect settings on instruments	Refer to instrument instructions for correct settings to increase sensitivity of light detection.
Alpha-counts 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.

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
 Please visit our website at: www.bpsbioscience.com