



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

LSD2(KDM1B) Homogeneous Assay Kit (Dimethyl)

Catalog #50614

DESCRIPTION: The *LSD2(KDM1B) Homogeneous Assay Kit (Dimethyl)* is designed to measure the activity of lysine-specific demethylase (LSD2) for screening and profiling applications. LSD2, also called KDM1B, is a chromatin-modifying enzyme that specifically removes methyl groups from mono- and dimethylated Lys of histone H3, thereby acting as a corepressor of transcription. LSD2 is a potential target for drug development. The *LSD2 Homogeneous Assay Kit (Dimethyl)* comes in a convenient AlphaLISA[®] format, with biotinylated, dimethylated histone H3 (2MeK4) peptide substrate, primary antibody to detect mono-methylated product, LSD2 assay buffer, and purified LSD2 for 384 enzyme reactions. The key to the *LSD2 Homogeneous Assay Kit (Dimethyl)* is a highly specific antibody that recognizes demethylated substrate. With this kit, only three simple steps on a microtiter plate are required for demethylase detection. First, a sample containing LSD2 enzyme is incubated with the biotinylated substrate. Next, acceptor beads are added, then donor beads, followed by reading the Alpha-counts.

COMPONENTS:

Cat. #		Amount	Storage	
50124	LSD2 (KDM1B)	40 µg	-80°C	(Avoid freeze/thaw cycles!)
52140V	Primary antibody 22	25 µl	-80°C	
	Biotinylated histone H3 peptide substrate (2MeK4)	500 rxns	-80°C	
	4x LSD2 assay buffer	3 ml	-20°C	
	4x Detection buffer	2 ml	-20°C	

MATERIALS REQUIRED BUT NOT SUPPLIED:

AlphaLISA anti-rIgG 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.

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 (NaN₃) or metal ions (Fe²⁺, Fe³⁺, Cu²⁺, Zn²⁺ and Ni²⁺). 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.

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

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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

REFERENCES: Binda, C., et al., J. Am. Chem. Soc. 2010; **132(19)**: 6827-6833.
Welte, M., et al., Cell Press. 2005; **15(14)**: 1266-1275.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Step 1:

- 1) Re-suspend lyophilized **Biotinylated histone H3 peptide substrate** in 500 μ l of distilled water.
- 2) Dilute 1 part **4x LSD2 Assay Buffer** with 3 parts distilled water (4-fold dilution) to make **1x LSD2 Assay Buffer**. Dilute **primary antibody 22** 10-fold using **1x LSD2 Assay Buffer**. Make only a sufficient quantity needed for the assay; store remaining undiluted buffer in aliquots at -20°C and undiluted antibody in aliquots at -80°C immediately.
- 3) Prepare master mix: N wells \times (1.5 μ l **4x LSD2 assay buffer** + 1 μ l **Biotinylated histone H3 peptide substrate** + 0.5 μ l **diluted primary antibody 22**). Add 3 μ l of master mix to each well.
- 4) Add 3 μ l of inhibitor solution to each well designated "Test Inhibitor". For the wells designated "Positive Control" and "Blank" add 3 μ l of the same solution without inhibitor (Inhibitor buffer).

	Positive Control	Test Sample	Blank
4x LSD2 assay buffer	1.5 μ l	1.5 μ l	1.5 μ l
Biotinylated substrate	1 μ l	1 μ l	1 μ l
Primary antibody 22 (10-fold diluted)	0.5 μ l	0.5 μ l	0.5 μ l
Test inhibitor	-	3	-
Inhibitor buffer (no inhibitor)	3	-	3
1x LSD2 assay buffer	-	-	4 μ l
LSD2 (20 ng/ μ l)	4 μ l	4 μ l	-
Total	10 μl	10 μl	10 μl

- 5) Add 4 μ l of **1x LSD2 assay buffer** to wells designated as "Blank".
- 6) Thaw LSD2 on ice. Upon first thaw, briefly spin tube containing enzyme to recover full contents of the tube. Aliquot LSD2 enzyme into single-use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. Note: LSD2 is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.

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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 7) Dilute **LSD2** in **1x LSD2 assay buffer** at -20 ng/ μ l. Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.
- 8) Initiate the reaction by adding 4 μ l of **diluted LSD2** prepared as described above to wells designated "Positive Control" and "Test Inhibitor". Incubate at room temperature for one hour. *Note: All incubations are done with slow shaking on a rotator platform.*

Step 2:

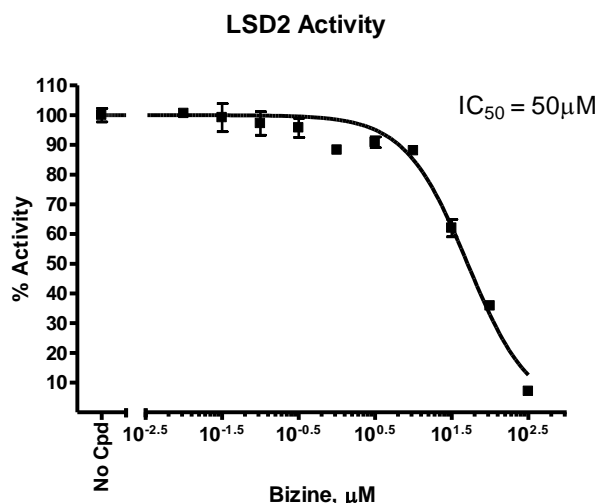
Note: Protect your samples from direct exposure to light!

- 1) Dilute anti-Rabbit Acceptor beads (PerkinElmer #AL104C) 1:500-fold with 1x Detection buffer. Add 10 μ l per well. Shake on a rotator platform for 30 minutes at room temperature.

Step 3:

- 1) Dilute Streptavidin-conjugated donor beads (PE #6760002S) 125-fold with 1x Detection buffer. Add 10 μ l per well. Shake on a rotator platform for 45 minutes at room temperature.
- 2) Read Alpha-counts.

Example of Assay Results:



LSD2 enzyme activity, measured using the LSD2 Homogeneous Assay Kit (Dimethyl), BPS Bioscience Cat. #50614. The compound was pre-incubated with the LSD2 enzyme for 30 min before the reaction was initiated with the addition of master mix. *Note: 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 **1.858.829.3082** 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: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS:

LSD2 (KDM1B) recombinant protein	#50124	20 µg
LSD1 recombinant protein	#50100	50 µg
LSD2 (KDM1B) Homogeneous Assay Kit (Monomethyl)	#50613	384 reactions
LSD1 Fluorescent Assay Kit	#50106	96 reactions
LSD1 Fluorescent Assay Kit	#50107	384 reactions
LSD1 substrate	#50101	500 µl
JMJD2A Homogeneous Assay Kit	#50413	384 reactions
JMJD2B Homogeneous Assay Kit	#50414	384 reactions
JMJD2C Homogeneous Assay Kit	#50415	384 reactions
JMJD2E Homogeneous Assay Kit	#50417	384 reactions
JMJD2C Assay Kit, Chemiluminescent	#50405	96 reactions
JMJD2A recombinant protein	#50103	20 µg
JMJD2B recombinant protein	#50104	20 µg
JMJD2C recombinant protein	#50105	20 µg
JMJD2E recombinant protein	#50118	20 µg

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

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