



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 ***LSD1 Homogeneous Assay Kit*** Catalog #50108

DESCRIPTION: The *LSD1 Homogeneous Assay Kit* is designed to measure LSD1 activity for screening and profiling applications. LSD1, also known as AOF2 and KDM1A, is a chromatin-modifying enzyme that specifically removes methyl groups from H3-K₄Me² and H3-K₄Me. The *LSD1 Homogeneous Assay Kit* comes in a convenient AlphaLISA[®] format (Scheme 1), with biotinylated histone H3 peptide substrate, primary antibody, demethylase assay buffer, and purified LSD1 for 384 enzyme reactions. The key to the *LSD1 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 LSD1 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	
50100	LSD1 (KDM1A)	20 µg	-80°C	(Avoid freeze/thaw cycles!)
52140J	Primary antibody 10	5 µl	-80°C	
79855	Biotinylated histone H3 peptide substrate	500 rxns	-80°C	
79849	4x LSD1 Assay Buffer 2A (add DTT before use*)	3 ml	-80°C	
52031	4x Detection buffer	2 ml	-20°C	

* Adding DTT can affect potency of some inhibitors.

MATERIALS REQUIRED BUT NOT SUPPLIED:

DTT (Dithiothreitol), 0.5M (Sigma, Cat. # D0632)
AlphaLISA[®] anti-mIgG acceptor beads, 5 mg/ml (PerkinElmer #AL105C)
AlphaScreen[®] Streptavidin-conjugated donor beads, 5 mg/ml (PerkinElmer #6760002S)
Optiplate -384 (PerkinElmer #6007290)
AlphaScreen[®] microplate reader
Adjustable micropipettor and sterile tips

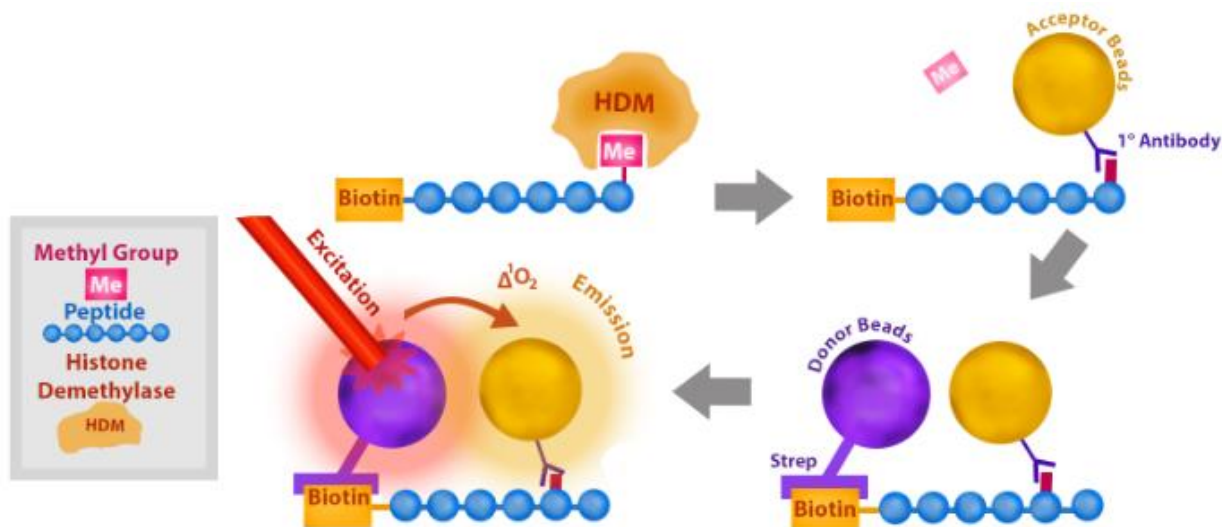
APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

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.

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 (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.

REFERENCE(S):

1. Forneris, F., Binda, C., Dall'Aglio, A., Fraaije, M.W., Battaglioli, E., and Mattevi, A. *J. Biol. Chem.* 2006; **281(46)**:35289-95.
2. Zhou, M., Diwu, Z., Panchuk-Voloshina, N., and Haugland, R.P. *Anal. Biochem.* 1997; **253(2)**:162-8.

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

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 LSD1 without the preincubation step.

Step 1:

- 1) Re-suspend lyophilized **Biotinylated histone H3 peptide substrate** in 500 μ l of distilled water.
- 2) Prepare serial dilutions of the test inhibitors in **1x LSD1 Assay Buffer 2A (Scheme 2)**. 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 LSD1 Assay Buffer 2A** with respective concentration of DMSO).
- 3) Thaw **LSD1** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **LSD1** enzyme into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C immediately. *Note: **LSD1** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 4) Dilute **LSD1** in **1x LSD1 Assay Buffer 2A** at 12.5 ng/ μ l (50 ng/4 μ l). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.
- 5) Preincubate 4 μ l of diluted **LSD1** 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 LSD1 Assay Buffer 2A**.
- 6) Prepare master mix: N wells \times (1.5 μ l **4x LSD1 Assay Buffer 2A** + 1 μ l **Biotinylated substrate** + 0.5 μ l **distilled water**).
- 7) 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.*

Scheme 2: The serial dilution of the compounds was first performed in 100% DMSO with the highest concentration at (X) mM. Each intermediate compound dilution (in 100% DMSO) will then get directly diluted 30x fold into **1x LSD1 Assay Buffer 2A** for 3.3x concentration (DMSO). From this intermediate step, 3 μ l of compound is added to 4 μ l of demethylase enzyme dilution is incubated for 30 minutes at room temperature. After this incubation, 3 μ l of peptide substrate is added. The final DMSO concentration is 1% for all wells.

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

Reagent	Blank	Positive Control	Test Inhibitor
1x LSD1 Assay Buffer 2A	4 μ l	–	–
4x LSD1 Assay Buffer 2A	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
Test Inhibitor/Activator	–	–	3 μ l
1x LSD1 Assay Buffer 2A (3.3% DMSO)	3 μ l	3 μ l	–
LSD1(12.5 ng/ μ l)	–	4 μ l	4 μ l
Total	10 μl	10 μl	10 μl

Step 2:

Note: Protect your samples from direct exposure to light!

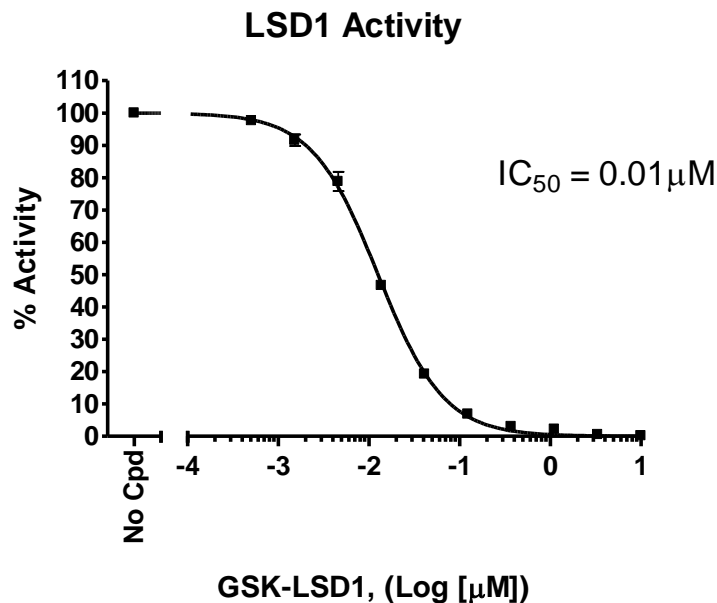
- 1) Dilute anti-Mouse Acceptor beads (PerkinElmer #AL105C) (1:500) and Primary antibody 10 (1:1600) with 1x Detection buffer in one step. Add 10 μ l of acceptor beads/antibody mixture per well. Incubate 30 min at room temperature.

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.

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

Example of Assay Results:


LSD1 enzyme activity, measured using the *LSD1 Homogeneous Assay Kit*, BPS Bioscience Cat. #50108. 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>
LSD2 Assay Kit, Homogeneous	50613	384 reactions
LSD2 Assay Kit, Homogeneous	50614	384 reactions
LSD1 Assay Kit, Fluorescent	50107	384 reactions
LSD1 Assay Kit, Fluorescent	50106	96 reactions
LSD1 Assay Kit, Chemiluminescence	50109	96 reactions
LSD1 recombinant protein	50100	50 μ g
LSD2 recombinant protein	50124	20 μ g
LSD1 recombinant protein	50097	50 μ g

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 **1.858.202.1401**, 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

TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Alpha-counts signal of positive control reaction is same as "blank" value.	LSD1 has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh JARID1A, BPS Bioscience #50100. Store enzyme in single-use aliquots. Increase time of enzyme incubation. Increase enzyme concentration.
	Streptavidin Donor beads or anti-mIgG acceptor beads fail to show significant signal.	Reorder Streptavidin Donor beads or anti-mIgG 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