

# Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric)

Base Catalog # P-3083

## PLEASE READ THIS ENTIRE USER GUIDE BEFORE USE

**Uses:** The Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric) is suitable for measuring activity or inhibition of total JARID using nuclear extracts or subtype JARID (JARID1A through JARID1D) purified enzymes from a broad range of species such as mammals, plant, fungal, and bacterial types, in a variety of forms including cultured cells and fresh tissues. Nuclear extracts can be prepared by using your own successful method. For your convenience and the best results, EpigenTek offers a nuclear extraction kit (Cat # OP-0002) optimized for use with this kit. Nuclear extracts can be used immediately or stored at  $-80^{\circ}\text{C}$  for future use. Purified enzymes can be active JARIDs from recombinant proteins or isolated from cell/tissues.

**Input Material:** Input materials can be nuclear extracts or purified JARID enzymes. The amount of nuclear extracts for each assay can be 0.5  $\mu\text{g}$  to 20  $\mu\text{g}$  with an optimal range of 2  $\mu\text{g}$  to 10  $\mu\text{g}$ . The amount of purified enzymes can be 10 ng to 500 ng, depending on the purity and subtypes of the enzymes.

**Internal Control:** A JARID assay standard (demethylated histone H3-K4) is provided in this kit for the quantification of JARID enzyme activity. Because JARID activity can vary from tissue to tissue, and from normal and diseased states, it is advised to run replicate samples to ensure that the signal generated is validated.

**Precautions:** To avoid cross-contamination, carefully pipette the sample or solution into the strip wells. Use aerosol-barrier pipette tips and always change pipette tips between liquid transfers. Wear gloves throughout the entire procedure. In case of contact between gloves and sample, change gloves immediately.

## KIT CONTENTS

Component	48 Assays Cat. #P-3083-48	96 Assays Cat. #P-3083-96	Storage Upon Receipt
<b>JD1</b> (10X Wash Buffer)	14 ml	28 ml	4°C
<b>JD2</b> (JARID Assay Buffer)	4 ml	8 ml	RT
<b>JD3</b> (JARID Substrate, 50 µg/ml)*	60 µl	120 µl	-20°C
<b>JD4</b> (JARID Assay Standard, 50 µg/ml)*	10 µl	20 µl	-20°C
<b>JD5</b> (Capture Antibody, 1000 µg/ml)*	5 µl	10 µl	4°C
<b>JD6</b> (Detection Antibody, 400 µg/ml)*	6 µl	12 µl	-20°C
<b>JD7</b> (Fluoro-Developer)*	10 µl	20 µl	-20°C
<b>JD8</b> (Fluoro-Enhancer)*	10 µl	20 µl	4°C
<b>JD9</b> ( Fluoro-Dilution)	4 ml	8 ml	RT
Co-factor 1*	30 µl	60 µl	4°C
Co-factor 2*	30 µl	60 µl	4°C
Co-factor 3*	30 µl	60 µl	4°C
8-Well Assay Strips (with frame)	6	12	4°C
Adhesive Covering Film	1	1	RT

\* Spin the solution down to the bottom prior to use.

## SHIPPING & STORAGE

The kit is shipped in two parts: the first part at ambient room temperature and the second and third parts on frozen ice packs at 4°C. Upon receipt: (1) Store **JD3**, **JD4**, **JD6**, and **JD7** at -20°C away from light; (2) Store **JD1**, **JD5**, **JD8**, **Co-factor 1**, **Co-factor 2**, **Co-factor 3**, and **8-Well Assay Strips** at 4°C away from light; (3) Store remaining components (**JD2**, **JD9**, and **Adhesive Covering Film**) at room temperature.

**Note:** (1) Check if **JD1** (10X Wash Buffer) contains salt precipitates before use. If so, warm (at room temperature or 37°C) and shake the buffer until the salts are re-dissolved.

All components of the kit are stable for 6 months from the date of shipment, when stored properly.

## MATERIALS REQUIRED BUT NOT SUPPLIED

- Adjustable pipette or multiple-channel pipette
- Multiple-channel pipette reservoirs
- Aerosol resistant pipette tips
- Fluorescence microplate reader capable of reading fluorescence at 530ex/590em nm
- 1.5 ml microcentrifuge tubes
- Incubator for 37°C incubation
- Plate seal

- Distilled water
- Nuclear extract or purified enzymes
- Parafilm M or aluminium foil

## GENERAL PRODUCT INFORMATION

**Quality Control:** Each lot of the Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric) is tested against predetermined specifications to ensure consistent product quality. EpigenTek guarantees the performance of all products in the manner described in our product instructions.

**Product Warranty:** If this product does not meet your expectations, simply contact our technical support unit or your regional distributor. We also encourage you to contact us if you have any suggestions about product performance or new applications and techniques.

**Safety:** Suitable lab coat, disposable gloves, and proper eye protection are required when working with this product.

**Product Updates:** EpigenTek reserves the right to change or modify any product to enhance its performance and design. The information in this User Guide is subject to change at any time without notice. Be sure to use the latest User Guide for this kit which can be accessed online at [www.epigentek.com/datasheet](http://www.epigentek.com/datasheet).

**Usage Limitation:** The Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric) is for research use only and is not intended for diagnostic or therapeutic application.

**Intellectual Property:** The Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric) and methods of use contain proprietary technologies by EpigenTek.

## A BRIEF OVERVIEW

Lysine histone methylation is one of the most robust epigenetic marks and is essential for the regulation of multiple cellular processes. The methylation of H3-K4 seems to be of particular significance, as it is associated with active regions of the genome. H3-K4 methylation was considered irreversible until the identification of a large number of histone demethylases indicated that demethylation events play an important role in histone modification dynamics. So far at least 2 classes of H3-K4 specific histone demethylase, LSD1(BHC110, KDM1) and JARIDs have been identified. The JARID family, except JARID2 (JARID1A, JARID1B, JARID1C and JARID1D), can remove tri-methylation from H3-K4. JARID demethylases are Jumonji-domain proteins and catalyze the removal of methylation by using a hydroxylation reaction with a requirement of iron and a-ketoglutarate as cofactors.

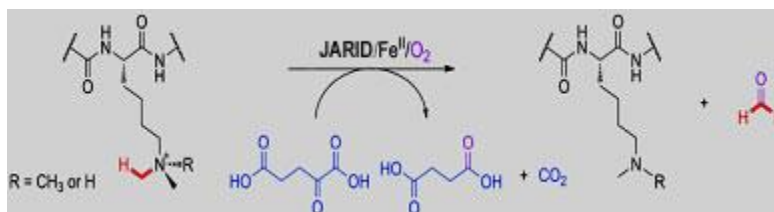


Fig 1. Histone H3-K4 demethylation reaction catalyzed by JARID demethylase.

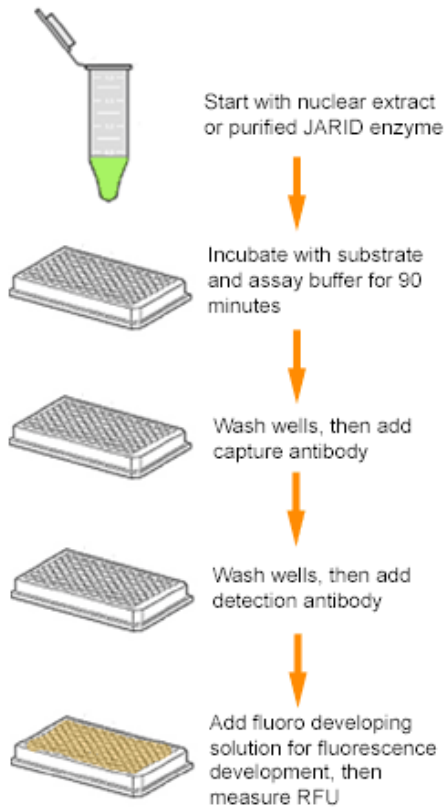
JARIDs function as transcription repressors and might participate in different biological processes through recruitment to different chromosomal regions and differing enzymatic activities. JARID demethylases are also found to have potential oncogenic functions. For example, all 5 members of JARID are overexpressed in bladder cancer, and may promote cancer progression by regulating E2F expression. Increased JARID1A expression is observed in gastric cancer and its inhibition triggers senescence of malignant cells. Detection of activity and inhibition of JARID would be important in elucidating mechanisms of epigenetic regulation of gene activation and silencing and benefiting cancer diagnostics and therapeutics.

Prior to this kit, there was only one method used for detecting JARID activity/inhibition. This method is based on the measurement of formaldehyde release, a by-product of JARID enzymatic reaction and has significant weaknesses: (1) a large amount (at  $\mu\text{g}$  level) of substrate and enzyme are required; (2) nuclear extracts from cell/tissues cannot be used; (3) redox-sensitive JARID inhibitors are not suitable for testing with this method; (4) high interference by SDS, DMSO, thiol-containing chemicals, and ions, which are often contained in enzyme solution, tested compound solvents and assay buffers; and (5) Less accuracy than direct measurement of JARID-converted demethylated products. The Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric) addresses all of these issues. Compared to a formaldehyde release-based method, this kit has the following advantages:

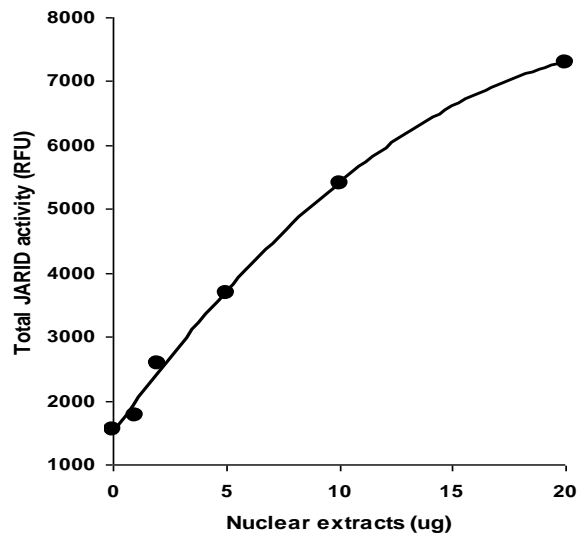
- 3-hour fluorometric procedure in a 96 stripwell microplate format allows for either manual or high throughput analysis.
- Directly measures JARID activity via a straightforward detection of JARID-converted demethylated products, rather than by-products, thus eliminating assay interference caused by thiol-containing chemicals such as DTT, GSH and 2-mercaptoethanol, or caused by detergents/ions such as tween-20, SDS, triton X-100, Fe, and Na.
- Both cell/tissue extracts and purified JARID proteins (including JARID1A, JARID1B, JARID1C, and JARID1D) can be used, which allows for the detection of inhibitory effects of JARID inhibitors in vivo and in vitro.
- Sensitivity is up to 2,000 times higher than formaldehyde release-based JARID assays, allowing activity to be fluorometrically detected from as low as 5 ng of purified JARID enzyme.
- Demethylated H3-K4 standard is included, allowing specific activity of JARID to be quantified.
- Accurate, reliable, and consistent with extremely low background signals.

## PRINCIPLE & PROCEDURE

In this assay, a tri-methylated histone H3-K4 substrate is stably coated onto microplate wells. Active JARIDs bind to the substrate and remove methyl groups from the substrate. The JARID-demethylated products can be recognized with a specific antibody. The ratio or amount of demethylated products, which is proportional to enzyme activity, can then be fluorometrically measured by reading the fluorescence in a fluorescent microplate reader at 530 excitation and 590 emission. The activity of the JARID enzyme is in turn proportional to the relative fluorescent units measured.



Schematic procedure of Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric)



Demonstration of high sensitivity of JARID activity assay achieved by using A549 nuclear extracts with Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Fluorometric).

## PROTOCOL

For the best results, please read the protocol in its entirety prior to starting your experiment.

### Starting Materials

**Input Amount:** The amount of nuclear extracts for each assay can be 0.5  $\mu$ g to 20  $\mu$ g with an optimal range of 2-10  $\mu$ g. The amount of purified enzymes can be 10 ng to 500 ng, depending on the purity and catalytic activity of the enzymes.

**Nuclear Extraction:** You can use your method of choice for preparing nuclear extracts. EpigenTek offers a nuclear extraction kit (Cat # OP-0002) optimized for use with this kit.

**Nuclear Extract or Purified JARID Storage:** Nuclear extract or purified JARID enzyme should be stored in aliquots at  $-80^{\circ}\text{C}$  until use.

### 1. Working Buffer and Solution Preparation

- a. Prepare **Diluted JD1** 1X Wash Buffer:

**48-Assay Kit:** Add 13 ml of **JD1** (10X Wash Buffer) to 117 ml of distilled water and adjust pH to 7.2-7.5.



96-Assay Kit: Add 26 ml of **JD1** (10X Wash Buffer) to 234 ml of distilled water and adjust pH to 7.2-7.5.

This **Diluted JD1** 1X Wash Buffer can now be stored at 4°C for up to six months.

- b. Prepare **CJD2** Completed JARID Assay Buffer:

Add **Co-Factor 1**, **Co-Factor 2**, and **Co-Factor 3** to **JD2** JARID Assay Buffer at a ratio of 1:100 (ex: add 1 µl of each **Co-Factor** to every 100 µl of **JC2**).

- c. Prepare **Diluted JD5** Capture Antibody Solution:

Dilute **JD5** Capture Antibody with **Diluted JD1** 1X Wash Buffer at a ratio of 1:1000 (ex: add 1 µl of **JD5** to 1000 µl of **Diluted JD1** 1X Wash Buffer). 50 µl of **Diluted JD5** will be required for each assay well.

- d. Prepare **Diluted JD6** Detection Antibody Solution:

Dilute **JD6** Detection Antibody with **Diluted JD1** 1X Wash Buffer at a ratio of 1:2000 (ex: add 1 µl of **JD6** to 2000 µl of 1 X wash buffer). 50 µl of **Diluted JD6** will be required for each assay well.

- e. Prepare **Diluted JD4** Standard Solution:

Suggested Standard Curve Preparation: First, dilute **JD4** with **JD2** to 5 ng/µl by adding 1 µl of **JD4** to 9 µl of **JD2**. Then, further prepare five concentrations by combining the 5 ng/µl **Diluted JD4** with **JD2** into final concentrations of 0.2, 0.5, 1, 2, and 5 ng/µl according to the following dilution chart:

Tube	JD4 (5 ng/µl)	JD2	Resulting JD4 Concentration
1	1.0 µl	24.0 µl	0.2 ng/µl
2	1.0 µl	9.0 µl	0.5 ng/µl
3	1.0 µl	4.0 µl	1.0 ng/µl
4	2.0 µl	3.0 µl	2.0 ng/µl
5	4.0 µl	0.0 µl	5.0 ng/µl

- f. Prepare **Fluorescence Development Solution**:

Add 1 µl of **JD7** Fluoro Developer and 1 µl of **JD8** Fluoro Enhancer to each 500 µl of **JD9** Fluoro Dilutor.

**Note**: Keep each of the diluted solutions except **Diluted JD1** 1X Wash Buffer on ice until use. Any remaining diluted solutions other than **Diluted D1** should be discarded if not used within the same day.

## 2. Enzymatic Reaction

- a. Predetermine the number of strip wells required for your experiment. It is advised to run replicate samples (include blank and positive control) to ensure that the signal generated is validated. Carefully remove un-needed strip wells from the plate frame and place them back in the bag (seal the bag tightly and store at 4°C).
- b. Blank Wells: Add 49 µl of **CJD2** and 1 µl of **JD3** to each blank well.

- c. **Standard Wells:** For a standard curve, add 49  $\mu\text{l}$  of **CJD2** and 1  $\mu\text{l}$  of **JD4** to each standard well with a minimum of five wells, each at different concentrations between 0.2 to 5 ng/ $\mu\text{l}$  (based on the dilution chart in Step 1e; see [Table 2](#) as an example).
- d. **Sample Wells Without Inhibitor:** Add 45 to 48  $\mu\text{l}$  of **CJD2**, 1  $\mu\text{l}$  of **JD3**, and 1 to 4  $\mu\text{l}$  of your nuclear extract or 1 to 4  $\mu\text{l}$  of purified JARID enzyme to each sample well without inhibitors. Total volume should be 50  $\mu\text{l}$  per well.
- e. **Sample Wells With Inhibitor:** Add 40 to 43  $\mu\text{l}$  of **CJD2**, 1  $\mu\text{l}$  of **JD3**, 1 to 4  $\mu\text{l}$  of nuclear extract or 1 to 4  $\mu\text{l}$  of purified JARID enzyme, and 5  $\mu\text{l}$  of inhibitor solution. Total volume should be 50  $\mu\text{l}$  per well.

**Note:** (1) Follow the suggested well setup diagrams; (2) It is recommended to use 2  $\mu\text{g}$  to 10  $\mu\text{g}$  of nuclear extract per well or 10 ng to 100 ng of purified enzyme per well; (3) The concentration of inhibitor to be added into the sample wells can be varied (e.g., 1  $\mu\text{M}$  to 1000  $\mu\text{M}$ ). However, the final concentration of the inhibitors before adding to the wells should be prepared with **JD2** at a 1:10 ratio (e.g., add 0.5  $\mu\text{l}$  of inhibitor to 4.5  $\mu\text{l}$  of **JD2**) so that the original solvent of the inhibitor can be reduced to 1% of the reaction solution or less. The Jumonji demethylase general inhibitor, N-Oxalylglycine can be used as the control inhibitor.

- f. Tightly cover the strip-well microplate with **Adhesive Covering Film** to avoid evaporation and incubate at 37°C for 60 to 120 min.

**Note:** (1) The incubation time may depend on intrinsic JARID activity. However, in general, 60-90 min incubation is suitable for active purified JARID enzyme and 90-120 min incubation is required for nuclear extract; (2) The **Adhesive Covering Film** can be cut to the required size to cover the strips based on the number of strips to be used.

- g. Remove the reaction solution from each well. Wash each well three times with 150  $\mu\text{l}$  of the **Diluted JD1** 1X Wash Buffer each time.

### **3. Antibody Binding & Signal Enhancing**

- a. Add 50  $\mu\text{l}$  of the **Diluted JD5** to each well, then cover with Parafilm M or aluminium foil and incubate at room temperature for 60 min.
- b. Remove the **Diluted JD5** solution from each well.
- c. Wash each well three times with 150  $\mu\text{l}$  of the **Diluted JD1** each time.
- d. Add 50  $\mu\text{l}$  of the **Diluted JD6** to each well, then cover with Parafilm M or aluminium foil and incubate at room temperature for 30 min.
- e. Remove the **Diluted JD6** solution from each well.
- f. Wash each well four times with 150  $\mu\text{l}$  of the **Diluted JD1** each time.

**Note:** Ensure any residual wash buffer in the wells is removed as much as possible at each wash step.

### **4. Signal Detection**

- a. Add 50  $\mu\text{l}$  of **Fluorescence Development solution** to each well and incubate at room temperature for 2 to 4 min away from light. **Fluorescence Development solution** will turn pink in the presence of sufficient demethylated products.

- b. Read the fluorescence on a fluorescence microplate reader within 2 to 10 min at 530ex/590em nm.

**Note:** If the strip-well plate frame does not fit in the fluorescence microplate reader, transfer the solution to a standard 96-well microplate.

### 5. JARID Activity Calculation

- a. Calculate the average duplicate readings for the sample wells and blank wells.  
 b. Calculate JARID activity or inhibition using the following formulas:

For simple calculation:

$$\text{JARID Activity (OD/min/mg)} = \frac{(\text{Sample RFU} - \text{Blank RFU})}{(\text{Protein Amount } (\mu\text{g}) * \text{min}^{**})} \times 1000$$

\* Protein amount ( $\mu\text{g}$ ) added into the reaction at Step 2d.

\*\* Incubation time (minutes) at Step 2f.

Example calculation:

Average RFU of sample is 6800

Average RFU of blank is 800

Protein amount is 5  $\mu\text{g}$

Incubation time is 2 hours (120 min)

$$\text{JARID activity} = \frac{(6800 - 800)}{(5 \times 120)} \times 1000 = 10000 \text{ RFU/min/mg}$$

For accurate or specific activity calculation:

1. Generate a standard curve and plot RFU value versus amount of **JD4** at each concentration point.
2. Determine the slope as RFU/ng (you can use Microsoft Excel statistical functions for slope calculation), then calculate the amount of JARID-converted demethylated product using the following formulas:

$$\text{Demethylated product (ng)} = \frac{(\text{Sample RFU} - \text{Blank RFU})}{\text{Slope}}$$

$$\text{JARID Activity (ng/min/mg)} = \frac{\text{Demethylated Product (ng)}}{(\text{Protein Amount } (\mu\text{g}) \times \text{min}^*)} \times 1000$$

\* Incubation time (minutes) at Step 2f.

For inhibition calculation:

$$\text{Inhibition \%} = \left[ 1 - \frac{\text{Inhibitor Sample RFU} - \text{Blank RFU}}{\text{No Inhibitor Sample RFU} - \text{Blank RFU}} \right] \times 100\%$$



## SUGGESTED WORKING BUFFER AND SOLUTION SETUP

**Table 2.** Approximate amount of required buffers and solutions for defined assay wells.

Reagents	1 well	1 strip (8 wells)	2 strips (16 wells)	6 strips (48 wells)	12 strips (96 wells)
Diluted JD1	2.5 ml	20 ml	40 ml	120 ml	240 ml
CJD2	50 µl	400 µl	800 µl	2400 µl	4800 µl
JD3	1 µl	8 µl	16 µl	50 µl	120 µl
JD4	NA	NA	1 µl (optional)	2 µl	2 µl
Diluted JD5	50 µl	400 µl	800 µl	2400 µl	4800 µl
Diluted JD6	50 µl	400 µl	800 µl	2400 µl	4800 µl
Fluorescence Development Solution	0.05 ml	0.4ml	0.8 ml	2.4 ml	4.8 ml

## SUGGESTED STRIP WELL SETUP

**Table 3.** The suggested strip-well plate setup for JARID activity assay in a 48-assay format (in a 96-assay format, Strips 7 to 12 can be configured as Sample). The controls and samples can be measured in duplicate.

Well #	Strip 1	Strip 2	Strip 3	Strip 4	Strip 5	Strip 6
A	Blank	Blank	Sample	Sample	Sample	Sample
B	JD4 0.2 ng	JD4 0.2 ng	Sample	Sample	Sample	Sample
C	JD4 0.5 ng	JD4 0.5 ng	Sample	Sample	Sample	Sample
D	JD4 1.0 ng	JD4 1.0 ng	Sample	Sample	Sample	Sample
E	JD4 2.0 ng	JD4 2.0 ng	Sample	Sample	Sample	Sample
F	JD4 5.0 ng	JD4 5.0 ng	Sample	Sample	Sample	Sample
G	Sample	Sample	Sample	Sample	Sample	Sample
H	Sample	Sample	Sample	Sample	Sample	Sample

## TROUBLESHOOTING

Problem	Possible Cause	Suggestion
No signal or weak signal in both the positive control and sample wells	Reagents are added incorrectly.	Check if reagents are added in the proper order with the right amount, and if any steps in the protocol may have been omitted by mistake.
	The well is incorrectly washed before enzyme reaction.	Ensure the well is not washed prior to adding the positive control and sample.
	Incubation time and temperature are incorrect.	Ensure the incubation time and temperature described in the protocol are followed correctly.
	Incorrect fluorescence reading.	Check if appropriate fluorescence wavelength (530ex/590em) is used.
	Kit was not stored or handled properly.	Ensure all components of the kit were stored at the appropriate temperature and the cap is tightly closed after each opening or use.

No signal or weak signal in only the standard curve wells	Insufficient standard amount is added to the well in Step 2c.	Ensure a sufficient amount of standard is added.
	The standard is degraded due to improper storage.	Follow the Shipping & Storage guidance in this User Guide for storage of <b>JD4</b> (JARID Assay Standard).
High background present in the blank wells	Insufficient washing of wells.	Check if washing recommendations at each step is performed according to the protocol.
	Contaminated by sample or standard.	Ensure the well is not contaminated from adding sample or standard accidentally or from using contaminated tips.
	Incubation time with <b>Diluted JD6</b> is too long.	The incubation time at Step 3d should not exceed 45 min.
	Over-development of fluorescence	Decrease the development time in Step 4a.
No signal or weak signal only in sample wells	Protein sample is not properly extracted or purified.	Ensure your protocol is suitable for JARID protein extraction. For the best results, it is advised to use EpigenTek's Nuclear Extraction Kit (Cat. No. OP-0002). Also, use fresh cells or tissues for protein extraction, as frozen cells or tissues could lose enzyme activity.
	Sample amount added into the wells is insufficient.	Ensure a sufficient amount of purified enzymes or nuclear extracts is used as indicated in Step 2. The sample can be titrated to determine the optimal amount for the assay.
	Sample was not stored properly or has been stored for too long.	Ensure sample is stored in aliquots at $-80^{\circ}\text{C}$ , with no more than 6 weeks for nuclear extracts and 6 months for purified enzymes. Avoid repeated freezing/thawing.
	Little or no activity of JARID contained in the sample.	This problem may be a result of many factors. If the affecting factors cannot be determined, use new or re-prepared nuclear extracts or purified enzymes.
Uneven fluorescent development	Insufficient washing of the wells.	Ensure the wells are washed as described and residue washing buffer is removed as much as possible.
	Delayed fluorescence development in the wells.	Ensure fluorescence development is added sequentially and is consistent with the order you added the other reagents (e.g., from well A - G or from well 1 - 12).

## RELATED PRODUCTS

### Nuclear Extract Preparation

OP-0002-1 EpiQuik™ Nuclear Extraction Kit

### Histone Demethylase Activity/Inhibition Assay

P-3077 EpiQuik™ Histone Demethylase (H3-K9 Specific) Activity/Inhibition Fast Assay Kit  
 P-3078 Epigenase™ LSD1 Demethylase Activity/Inhibition Assay Kit (Colorimetric)  
 P-3079 Epigenase™ LSD1 Demethylase Activity/Inhibition Assay Kit (Fluorometric)  
 P-3080 Epigenase™ JMJD2 Demethylase Activity/Inhibition Assay Kit (Colorimetric)  
 P-3081 Epigenase™ JMJD2 Demethylase Activity/Inhibition Assay Kit (Fluorometric)  
 P-3082 Epigenase™ JARID Demethylase Activity/Inhibition Assay Kit (Colorimetric)