



6405 Mira Mesa Blvd. Suite 100  
San Diego, CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: support@bpsbioscience.com

## Data Sheet

### ***FTO Chemiluminescent Assay Kit***

**Catalog # 79344**

**BACKGROUND:** The methylation of internal adenosines at the N<sup>6</sup> position (m<sup>6</sup>A) in RNA has been shown to play essential roles in various normal and pathologic processes, including various cancers, leukemia, Alzheimer's disease, and obesity. FTO (Fat mass and obesity-associated protein), also known as ALKBH9 and BMIQ14, is the first m<sup>6</sup>A demethylase identified.

**DESCRIPTION:** The *FTO Chemiluminescent Assay Kit* is designed to measure FTO activity for screening and profiling applications. The *FTO Chemiluminescent Assay Kit* comes in a convenient format, with a 96-well strip plate precoated with methylated FTO substrate, primary antibody, the secondary HRP-labeled antibody, demethylase assay buffer, and purified FTO for 96 enzyme reactions. The key to the *FTO Chemiluminescent Assay Kit* is a highly specific antibody that recognizes methylated substrate. Signal is inversely related to the activity. With this kit, only three simple steps on a microtiter plate are required for detection of demethylase activity. First, FTO enzyme is incubated with the substrate. Next, primary antibody is added. Finally, the plate is treated with an HRP-labeled secondary antibody followed by the addition of the HRP substrate to produce chemiluminescence that can be measured using a chemiluminescence reader.

#### **COMPONENTS:**

Catalog #	Component	Amount	Storage	
79306	FTO	100 µg	-80°C	<b>Avoid freeze/ thaw cycles!</b>
52140Z4	Primary antibody 29	100 µl	-80°C	
52131H	Secondary HRP-labeled antibody 2	10 µl	-80°C	
	4x FTO assay buffer	3 x 1 ml	-80°C	
79556	Blocking buffer 1	50 ml	+4°C	
79670	ELISA ECL substrate A (transparent bottle)	6 ml	RT	
	ELISA ECL substrate B (brown bottle)	6 ml	RT	
	White 8-well strip plate module precoated with RNA substrate*	1	+4°C	

**\* Make sure you are working in RNase-free conditions.**

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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



6405 Mira Mesa Blvd. Suite 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

**MATERIALS REQUIRED BUT NOT SUPPLIED:**

TBST buffer (1 x TBS, pH 8.0, containing 0.05% Tween-20)  
Luminometer or microplate reader capable of reading chemiluminescence  
Adjustable micropipettor and sterile tips  
Rotating or rocker platform

**APPLICATIONS:** Great for studying enzyme kinetics and HTS applications.

**CONTRAINDICATIONS:** DMSO >1%, strong acids or bases, ionic detergents, high salt

**STABILITY:** 6 months from date of receipt when stored as directed.

**REFERENCE:** Zhou Y., *et al. J Biomed Sci.* 2017; **24**(1):65.

**ASSAY PROTOCOL:**

All samples and controls should be tested in duplicate.

**Step 1:**

- 1) Rehydrate the microwells by adding 200 µl of TBST buffer (1 x TBS, pH 8.0, containing 0.05% Tween-20) to every well. Incubate 15 minutes at room temperature. Tap the strip plate onto clean paper towels to remove liquid.
- 2) Prepare master mix: N wells x (7.5 µl **4x FTO Assay Buffer** + 17.5 µl water). Add 25 µl of master mixture to each well.
- 3) Add 5 µl of inhibitor solution of each well designated "Test Inhibitor". For the "Positive Control" and "Blank" add 5 µl of the same solution without inhibitor (Inhibitor buffer). *Note: Keep final DMSO concentration ≤1%.*

	Blank	Positive Control	Test Inhibitor
4x FTO assay buffer	7.5 µl	7.5 µl	7.5 µl
Distilled water	17.5 µl	17.5 µl	17.5 µl
Test Inhibitor/Activator	-	-	5 µl
Inhibitor buffer (no inhibitor)	5 µl	5 µl	-
1x FTO buffer	20 µl	-	-
Diluted FTO (50 ng/µl)	-	20 µl	20 µl
Total	50 µl	50 µl	50 µl

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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)

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



6405 Mira Mesa Blvd. Suite 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

- 4) Prepare **1x FTO Buffer** by diluting 1 part **4x FTO Assay Buffer** with three parts distilled water. Prepare only enough **1x FTO Buffer** for the assay. Add 20  $\mu$ l of **1x FTO buffer** to wells designated as "Blank".
- 5) Thaw **FTO** on ice. Upon first thaw, briefly spin tube containing enzyme to recover full content of the tube. Aliquot **FTO** enzyme into single use aliquots. Store remaining undiluted enzyme in aliquots at  $-80^{\circ}\text{C}$ . *Note: FTO is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.*
- 6) Dilute **FTO** in **1x FTO Buffer** at 50 ng/ $\mu$ l (1000 ng/reaction). Keep diluted enzyme on ice until use. Discard any unused diluted enzyme after use.
- 7) Initiate reaction by adding 20  $\mu$ l of **diluted FTO** prepared as described above to wells designated "Positive Control" and "Test Inhibitor". Incubate overnight at room temperature with slow shaking. Glue the wells if necessary.
- 8) Wash the strip plate three times with TBST buffer. Blot dry onto clean paper towels.
- 9) Add 100  $\mu$ l of **Blocking buffer 1** to every well. Shake on a rotating platform for 10 minutes. Remove supernatant as described above.

#### Step 2:

- 1) Dilute **Primary antibody 29** 100-fold with **Blocking Buffer 1**.
- 2) Add 100  $\mu$ l per well. Incubate 1 hour at room temperature with slow shaking.
- 3) Remove supernatant from the wells and wash the strip three times with 200  $\mu$ l of TBST buffer and incubate in **Blocking Buffer 1** as described in steps 1-8 and 1-9.

#### Step 3:

- 1) Dilute **Secondary HRP-labeled antibody 2** 1,000-fold with **Blocking Buffer 1**.
- 2) Add 100  $\mu$ l per well. Incubate for 30 min. at room temperature with slow shaking.
- 3) Remove supernatant from the wells and wash the strip three times with 200  $\mu$ l of TBST buffer and incubate in **Blocking Buffer 1** as described in steps 1-8 and 1-9
- 4) Just before use, mix on ice 50  $\mu$ l **ELISA ECL substrate A** and 50  $\mu$ l **ELISA ECL substrate B** and add 100  $\mu$ l per well. Discard any unused chemiluminescent reagent after use.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

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



6405 Mira Mesa Blvd. Suite 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

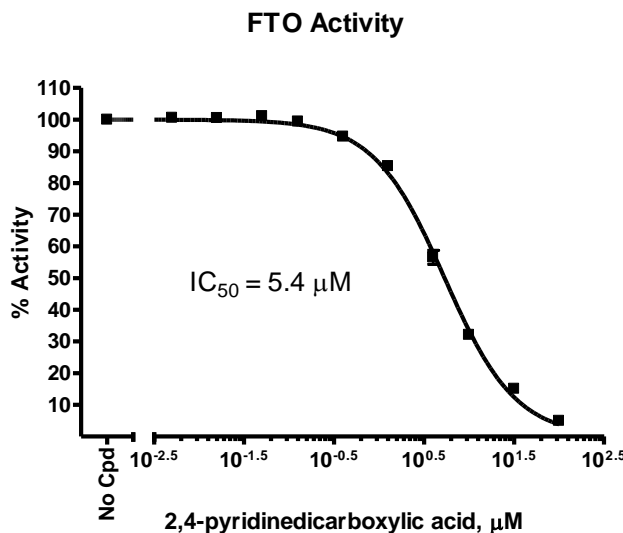
- 5) Immediately read sample in a luminometer or microtiter-plate capable of reading chemiluminescence.

### Reading Chemiluminescence:

Chemiluminescence is the emission of light (luminescence) which results from a chemical reaction. The detection of chemiluminescence requires no wavelength selection because the method used is emission photometry and is not emission spectrophotometry.

To properly read chemiluminescence, make sure the plate reader is set for LUMINESCENCE mode. Typical integration time is 1 second, delay after plate movement is 100 msec. Do not use a filter when measuring light emission. Typical settings for the Synergy 2 BioTek plate reader are: use the "hole" position on the filter wheel; Optics position: Top; Read type: endpoint. Sensitivity may be adjusted based on the luminescence of a control assay without enzyme (typically we set this value as 40000). Make sure signal decrease corresponds to increased activity.

### Examples of Assay Results:



FTO enzyme inhibition by 2-4-pyridinedicarboxylic acid (2,4-PC), measured using the *FTO Chemiluminescent Assay Kit*, BPS Bioscience #79344. Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [support@bpsbioscience.com](mailto:support@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: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6405 Mira Mesa Blvd. Suite 100  
San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

## RELATED PRODUCTS

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
JHDM1D (KDM7A) recombinant protein	50419	20 µg
JMJD1A (KDM3A) recombinant protein	50130	20 µg
JARID1A recombinant protein	50155	20 µg
JARID1B recombinant protein	50121	20 µg
JARID1C recombinant protein	50112	20 µg
JMJD2A recombinant protein	50123	100 µg
JMJD2B recombinant protein	50111	100 µg
JMJD2C recombinant protein	50105	100 µg
JMJD2E recombinant protein	50118	100 µg
LSD1 recombinant protein	50100	50 µg
JHDM1D (KDM7A) Homogeneous Assay Kit	50420	384 reactions
JARID1A Homogeneous Assay Kit	50510	384 reactions
JARID1B Homogeneous Assay Kit	50512	384 reactions
JARID1C Homogeneous Assay Kit	50511	384 reactions
JMJD2A Homogeneous Assay Kit	50413	384 reactions
JMJD2B Homogeneous Assay Kit	50414	384 reactions
JMJD2C Homogeneous Assay Kit	50415	384 reactions
JMJD2C Chemiluminescent Assay Kit	50405	96 reactions
JMJD2D Chemiluminescent Assay Kit	50418	96 reactions
JMJD3 Chemiluminescent Assay Kit	50406	96 reactions
LSD1 Chemiluminescent Assay Kit	50109	96 reactions

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

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

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



6405 Mira Mesa Blvd. Suite 100  
 San Diego, CA 92121  
**Tel:** 1.858.202.1401  
**Fax:** 1.858.481.8694  
**Email:** support@bpsbioscience.com

### TROUBLESHOOTING GUIDE

Problem	Possible Cause	Solution
Luminescence signal of positive control reaction is same as "blank" value.	FTO has lost activity	Enzyme loses activity upon repeated freeze/thaw cycles. Use fresh FTO, BPS Bioscience #79306. 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 FTO, BPS Bioscience #79306 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 **1.858.202.1401** Fax **1.858.481.8694**  
 Or you can Email us at: [support@bpsbioscience.com](mailto:support@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)