

Data Sheet

TAF1 (BD2) Inhibitor Screening Assay Kit

Catalog # 32624
Size: 384 reactions

DESCRIPTION: The *TAF1 (BD2) Inhibitor Screening Assay Kit* is designed to measure the inhibition of TAF1 (BD2) binding to its substrate. The *TAF1 (BD2) Inhibitor Screening Assay Kit* comes in a convenient AlphaLISA[®] format, with biotinylated histone peptide substrate, assay buffer, detection buffer and purified GST-tagged TAF1 bromodomain 2 to perform a total of 384 enzyme reactions. The key to the *TAF1 (BD2) Inhibitor Screening Assay Kit* is the specific binding of the TAF1 bromodomain 2 (a.a. 1519-1651) to the acetylated histone substrate. With this kit, only three simple steps on a microtiter plate are required. First, a sample containing *TAF1 (BD2)* and an inhibitor of choice is incubated with the biotinylated substrate for thirty minutes. Next, acceptor beads are added, then donor beads, followed by reading the Alpha-counts.

COMPONENTS:

Catalog #	Component	Amount	Storage	
31126	TAF1-BD2 (1519-1651), GST-tag	24 µg	-80 °C	(Avoid freeze/thaw cycles!)
	BET Ligand 1	400 µl	-80 °C	
	Non-acetylated BET Ligand 1	200 µl	-80 °C	
33007	3x BRD Homogeneous Assay Buffer 2	4 ml	-20 °C	
33006	3x BRD Homogeneous Detection Buffer 2	3 ml	-20 °C	

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Glutathione AlphaLISA[®] Acceptor Beads, 5 mg/ml (PerkinElmer #AL109C)
 AlphaScreen[®] Streptavidin-conjugated donor beads, 5 mg/ml (PerkinElmer #6760002)
 Optiplate-384 (PerkinElmer #6007290)
 AlphaScreen[®] microplate reader
 Adjustable micropipettor and sterile tips

APPLICATIONS: Useful for the study of bromodomain binding assays, screening inhibitors and selectivity profiling.

CONTRAINDICATIONS: Keep DMSO levels below 0.5%. Avoid 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

REFERENCE(S): Filippakopoulos, P., *et al.*, *Cell* 2012; **149**:214.

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate. Use slow shaking for all incubations.

Step 1:

- 1) Prepare the master mixture: N wells × (2.5 µl **3x BRD Homogeneous Assay Buffer 2** + 1 µl **BET Ligand 1** + 1.5 µl **H₂O**).
- 2) Add 5 µl of master mixture to each well designated for the “Positive Control”, “Test Inhibitor”, and “Blank”. For the “Substrate Control”, add 2.5 µl **3x BRD Homogeneous Assay Buffer 2** + 1.5 µl **H₂O** + 1 µl of **Non-acetylated Ligand 1**.
- 3) Thaw **TAF1-BD2** on ice. Upon first thaw, briefly spin tube containing protein to recover full content of the tube. Aliquot protein into single use aliquots. Store remaining undiluted protein in aliquots at -80°C immediately. *Note: TAF1-BD2 is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 4) Dilute **TAF1-BD2** in 1x **BRD Homogeneous Assay Buffer 2** at 24 ng/µl. Keep diluted protein on ice until use. Discard any unused diluted protein after use.

	Blank	Substrate Control	Positive Control	Test Inhibitor
3x BRD Homogeneous Assay Buffer 2	2.5 µl	2.5 µl	2.5 µl	2.5 µl
BET Ligand 1	1 µl	-	1 µl	1 µl
Non-acetylated BET Ligand 1	-	1 µl	-	-
H ₂ O	1.5 µl	1.5 µl	1.5 µl	1.5 µl
Test Inhibitor/Activator	-	-	-	2.5 µl
Inhibitor buffer (no inhibitor)	2.5 µl	2.5 µl	2.5 µl	-
1x BRD Homogeneous Assay Buffer 2	2.5 µl			
TAF1-BD2 (24 ng/µl)	-	2.5 µl	2.5 µl	2.5 µl
Total	10 µl	10 µl	10 µl	10 µl

- 5) Add 2.5 µl of **test inhibitor solution** to each well designated “Test Inhibitor”. For the “Positive Control”, “Substrate Control” and “Blank”, add 2.5 µl of the same solution without inhibitor (**inhibitor buffer**). *Note: Keep final DMSO concentration below 0.5%.*
- 6) Add 2.5 µl of **1x BRD Homogeneous Assay Buffer 2** to the well designated “Blank”.

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

- 7) Initiate reaction by adding 2.5 μ l of diluted **TAF1-BD2** prepared as described above. Incubate at room temperature for 30 minutes.

Step 2:

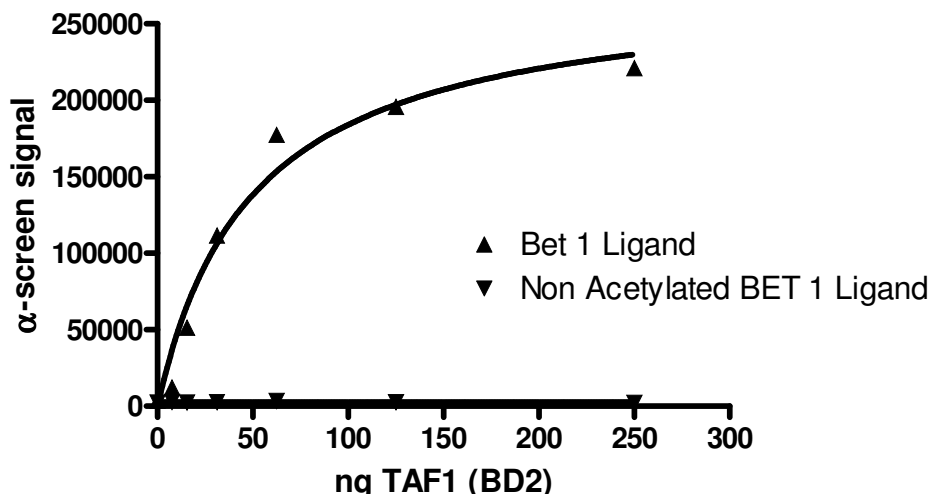
Note: Protect your samples from direct exposure to light!

- 1) Dilute Glutathione AlphaLISA[®] Acceptor Beads (PerkinElmer #AL109C) 250-fold with **1x BRD Homogeneous Detection Buffer 2**. Add 10 μ l per well. Shake plate briefly. Incubate at room temperature for 30 minutes.

Step 3:

- 1) Dilute Streptavidin-conjugated donor beads (PE #6760002S) 125-fold with **1x BRD Homogeneous Detection Buffer 2**. Add 10 μ l per well. Incubate at room temperature for 15 - 30 minutes.
- 2) Read Alpha-counts.

Example of Assay Results:



TAF1-BD2 binding activity, measured using the *TAF1 (BD2) Inhibitor Screening Assay Kit*, BPS Bioscience, Catalog # 32624. *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



6044 Cornerstone Ct. West, Ste. E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
TAF1 (1400-1518), His-tag	31123	100 µg
TAF1 (1519-1651), GST-tag	31126	100 µg
TAF1 (1519-1651), His-tag	31110	100 µg
TAF1, BD1 and BD2 (1400-1651), GST-tag	31124	100 µg
TAF1L (1398-1516), GST-tag	31105	100 µg
TAF1L (1398-1516), His-tag	31103	100 µg
TAF1L (1398-1649), GST-tag	31107	100 µg
TAF1L (1517-1649), GST-tag	31106	100 µg
TAF1L (1517-1649), His-tag	31104	100 µg
TAF1 (BD1 + BD2) Inhibitor Screening Kit	32604	384 rxns.
TAF1L (BD1 + BD2) Inhibitor Screening Kit	32603	384 rxns.
TAF1L (BD2) Inhibitor Screening Kit	32602	384 rxns.
(+)-JQ1 Bromodomain Inhibitor	27400	10 mg

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