



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

**Data Sheet**  
**TAF1 (BD2) TR-FRET Assay Kit**  
**Catalog # 32627**

**DESCRIPTION:**

The TAF1 (BD2) TR-FRET Assay Kit is designed to measure the inhibition of TAF1 (BD2) binding to its substrate in a homogeneous 384 reaction format. This FRET-based assay requires no time-consuming washing steps, making it especially suitable for high throughput screening applications. The assay procedure is straightforward and simple; a sample containing terbium-labeled donor, dye-labeled acceptor, TAF1 (BD2), substrate, and an inhibitor is incubated for 120 minutes. Then, the fluorescence intensity is measured using a fluorescence reader.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
31126	TAF1 (BD2), GST-tag	5 µg	-80°C	<b>(Avoid freeze/ thaw cycles!)</b>
	BET Bromodomain Ligand	50 µl	-80°C	
	Non-acetylated Ligand 1	15 µl	-80°C	
	Tb donor	20 µl	-20°C	
	Dye-labeled acceptor	20 µl	-20°C	
33012	3x BRD TR-FRET Assay Buffer 1	4 ml	-20°C	
Fisher 07-200-330	White, Nonbinding Corning, low volume, microtiter plate	1	Room temp.	

**MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

Fluorescent microplate reader capable of measuring Time Resolved Fluorescence Resonance Energy Transfer (TR-FRET)  
Adjustable micropipettor and sterile tips

**APPLICATIONS:** Great for screening small molecular inhibitors for drug discovery and HTS applications.

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

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

**ASSAY PROTOCOL:**

All samples and controls should be tested in duplicate.

- 1) Dilute one part **3x BRD TR-FRET Assay Buffer 1** with 2 parts distilled water (3-fold dilution) to make **1x BRD TR-FRET Assay Buffer 1**. Make only a sufficient quantity needed for the assay; store remaining stock solution in aliquots at -20°C.

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](mailto:info@bpsbioscience.com)

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



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

- Dilute **Tb-labeled donor** and **Dye-labeled acceptor** 100-fold in **1x BRD TR-FRET Assay Buffer 1**. Make only sufficient quantities needed for the assay; store remaining stock solution in aliquots at -20°C.
- Add 5 µl of diluted **Tb-labeled donor**, and 5 µl of diluted **Dye-labeled acceptor** to every well.
- Add 2 µl of inhibitor solution to each well designated "Test Inhibitor". Add 2 µl of the same solution without inhibitor (inhibitor buffer) to the wells labeled "Substrate Control", and "Positive Control".

	Positive Control	Negative* Control	Test Inhibitor
Tb-labeled donor	5 µl	5 µl	5 µl
Dye-labeled acceptor	5 µl	5 µl	5 µl
Test Inhibitor	-	-	2 µl
Inhibitor Buffer (no inhibitor)	2 µl	2 µl	-
BET Bromodomain Ligand	5 µl	-	5 µl
Non-acetylated Ligand 1	-	-	-
1x TAF1 Buffer	-	5 µl*	-
TAF1 (BD2) (0.5 ng/µl)	3 µl	3 µl	3 µl
<b>Total</b>	<b>20 µl</b>	<b>20 µl</b>	<b>20 µl</b>

\***Non-acetylated Ligand 1** may be used as a substrate control in place of the negative control.

- Thaw **BET Bromodomain Ligand** on ice. Upon first thaw, briefly spin tube containing ligand to recover the full contents of the tube. Aliquot each ligand into single-use aliquots. Store remaining undiluted ligand at -80°C immediately. *Note: each ligand is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots.*
- Individually dilute **BET Bromodomain Ligand** 50-fold in **1x BRD TR-FRET Assay Buffer 1**. Add 5 µl of diluted **BET Bromodomain Ligand** to each well designated as "Positive Control" and "Test Inhibitor". Add 5 µl of **1x BRD TR-FRET Assay Buffer 1** to the wells labeled "Negative Control". *Note: if using the **Non-acetylated Ligand 1**, dilute **Non-acetylated Ligand 1** 50-fold in **1x BRD TR-FRET Assay Buffer 1** and add 5 µl of diluted **Non-acetylated Ligand 1** to the "Negative Control" well in place of the 5 µl of **1x BRD TR-FRET Assay Buffer 1**.*
- Thaw **TAF1 (BD2)** bromodomain protein on ice. Upon first thaw, briefly spin tube containing protein to recover the full contents of the tube. Aliquot **TAF1 (BD2)** protein into single-use aliquots. Store remaining undiluted **TAF1 (BD2)** 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.*

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](mailto:info@bpsbioscience.com)

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



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

- 8) Dilute **TAF1 (BD2)** in **1x BRD TR-FRET Assay Buffer 1** to 0.5 ng/μl (1.5 ng/reaction). Initiate reaction by adding 3 μl of diluted **TAF1 (BD2)** to every well. Discard any remaining diluted TAF1 (BD2) protein after use.
- 9) Incubate at room temperature for 2 hours.
- 10) Read the fluorescent intensity in a microtiter-plate reader capable of TR-FRET.

#### Instrument Settings

Reading Mode	Time Resolved
Excitation Wavelength	340±20 nm
Emission Wavelength	620±10 nm
Lag Time	60 μs
Integration Time	500 μs
Excitation Wavelength	340±20 nm
Emission Wavelength	665±10 nm
Lag Time	60 μs
Integration Time	500 μs

#### CALCULATING RESULTS:

Two sequential measurements should be conducted. Tb-donor emission should be measured at 620 nm followed by dye-acceptor emission at 665 nm. Data analysis is performed using the TR-FRET ratio (665 nm emission/620 nm emission).

When percentage activity is calculated, the FRET value from the negative control can be set as zero percent activity and the FRET value from the positive control can be set as one hundred percent activity.

$$\% \text{ Activity} = \frac{FRET_s - FRET_{neg}}{FRET_p - FRET_{neg}} \times 100\%$$

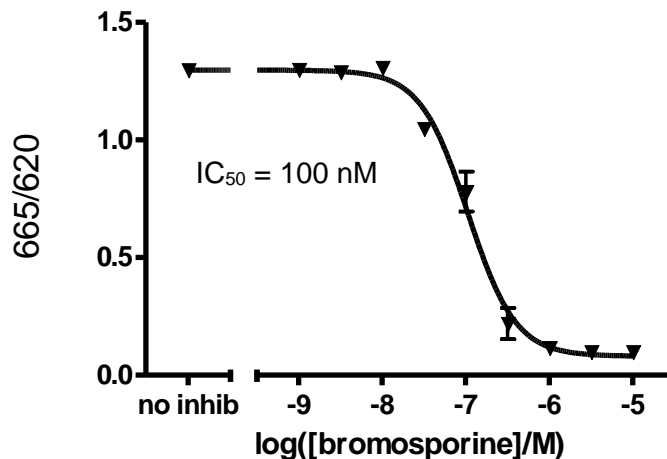
Where  $FRET_s$  = Sample FRET,  $FRET_{neg}$  = negative control FRET, and  $FRET_p$  = Positive control FRET.

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](mailto:info@bpsbioscience.com)

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

**EXAMPLE OF ASSAY RESULTS:**

Inhibition of TAF1 (BD2) (BPS Bioscience Cat. #31126) with Bromosporine (BPS Cat. #27612). Assay was done according to protocol for the TAF1 (BD2) TR-FRET Assay Kit (BPS Cat. #32627). Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto: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](mailto:info@bpsbioscience.com)

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



6044 Cornerstone Court W, Ste E  
San Diego, CA 92121  
**Tel:** 1.858.829.3082  
**Fax:** 1.858.481.8694  
**Email:** [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

**RELATED PRODUCTS:**

<b><u>Product Name</u></b>	<b><u>Catalog</u></b>	<b><u>Size</u></b>
BET Bromodomain Ligand	33000	0.5 ml
Bromodomain Non-acetylated Ligand 1	33005	0.5 ml
TAF1L (1398-1649), GST-tag	31107	100 µg
TAF1L (1398-1516), GST-tag*	31105	100 µg
TAF1L (1517-1649), GST-tag*	31106	100 µg
TAF1 (1519-1651), GST-tag*	31126	100 µg
TAF1, BD1 and BD2 (1400-1651), GST-tag	31124	100 µg
TAF1 (1400-1518), His-tag	31123	100 µg
TAF1 (BD1+BD2) Inhibitor Screening Kit	32604	384 rxns.
TAF1 (BD2) Inhibitor Screening Kit	32624	384 rxns.
TAF1L (BD1+BD2) Inhibitor Screening Kit	32603	384 rxns.
TAF1L (BD2) Inhibitor Screening Kit	32602	384 rxns.
Bromosporine	27612	1 mg
(+)-JQ1 Inhibitor	27401	1 mg

\*also available as His-tag

Note: Tb-labeled donor and dye-labeled acceptor are products of Cisbio Bioassays.

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](mailto:info@bpsbioscience.com)

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