



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
BRD4 (BD1+BD2) TR-FRET Assay Kit
Catalog # 32612
Size: 384 reactions

DESCRIPTION: The BRD4 (BD1+BD2) TR-FRET Assay Kit is designed to measure the inhibition of BRD4 (BD1+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, BRD4, substrate, and an inhibitor is incubated for 120 minutes. Then, the fluorescence intensity is measured using a fluorescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
31044	BRD4 (BD1+BD2)	10 µg	-80°C	<i>Avoid freeze/ thaw cycles!</i>
33000	BET Bromodomain Ligand	50 µl	-80°C	
33005	Non-acetylated Ligand 1	15 µl	-80°C	
	Tb-labeled donor	2 x 10 µl	-20°C	
	Dye-labeled acceptor	2 x 10 µl	-20°C	
33012	3x BRD TR-FRET Assay Buffer 1	4 ml	-20°C	
	White, Nonbinding, low volume, microtiter plate	1	Room temp.	

MATERIALS 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):

1. Filippakopoulos, P., *et al.*, *Cell* 2012; **149**:214.

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**](mailto:info@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.

Protocol for BRD4 (BD1+BD2) Assay

- 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.
- 2) 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.
- 3) Add 5 µl of diluted **Tb-labeled donor**, and 5 µl of diluted **dye-labeled acceptor** to each well designated "Test Inhibitor," "Negative Control," and "Positive Control."
- 4) Add 2 µl of inhibitor solution to each well designated "Test Inhibitor." Add 2 µl of 10% DMSO in water (inhibitor buffer) to the wells labeled "Negative Control" and "Positive Control."

	Negative Control*	Positive 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
10% DMSO in water (Inhibitor Buffer)	2 µl	2 µl	–
BET Bromodomain Ligand		5 µl	5 µl
Non-acetylated Ligand 1	5 µl	–	–
BRD4 (BD1 + BD2) 6 ng/µl	3 µl	3 µl	3 µl
Total	20 µl	20 µl	20 µl

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

Note: The BRD4 (BD1+BD2) TR-FRET Assay Kit is compatible with up to 0.5% final DMSO concentration.

- 5) Thaw **BET Bromodomain Ligand** and **Non-acetylated Ligand 1** on ice. Upon first thaw, briefly spin tube containing enzyme 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.

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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

- 6) Individually dilute **BET Bromodomain Ligand** 40-fold in **1x BRD TR-FRET Assay Buffer 1**. Add 5 μ L of diluted **BET Bromodomain Ligand** to each well designated as "Positive Control" and "Test Inhibitor". Add 5 μ L of **1x BRD TR-FRET Assay Buffer 1** to the wells labeled as "Negative Control." *Note: if using the **Non-acetylated Ligand 1**, dilute **Non-acetylated Ligand 1** 40-fold in **1x BRD TR-FRET Assay Buffer 1** and add 5 μ L of diluted **Non-acetylated Ligand 1** to the "Negative Control" well in place of the 5 μ L of **1x BRD TR-FRET Assay Buffer 1**.*
- 7) Thaw **BRD4 (BD1 + BD2)** bromodomain protein on ice. Upon first thaw, briefly spin tube containing protein to recover the full contents of the tube. Aliquot **BRD4 (BD1 + BD2)** protein into single-use aliquots. Store remaining undiluted **BRD4** in aliquots at -80°C immediately. *Note: **BRD4** is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 8) Dilute **BRD4 (BD1 + BD2)** in **1x BRD TR-FRET Assay Buffer 1** to 6 ng/ μ L (18 ng/reaction). Initiate reaction by adding 3 μ L of diluted **BRD4 (BD1 + BD2)** to wells designated for the "Negative Control," "Positive Control," and "Test Inhibitor." Discard any remaining diluted BRD 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 \pm 20 nm
Emission Wavelength	620 \pm 10 nm
Lag Time	60 μ s
Integration Time	500 μ s
Excitation Wavelength	340 \pm 20 nm
Emission Wavelength	665 \pm 10 nm
Lag Time	60 μ s
Integration Time	500 μ s

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**](mailto:info@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

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.

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

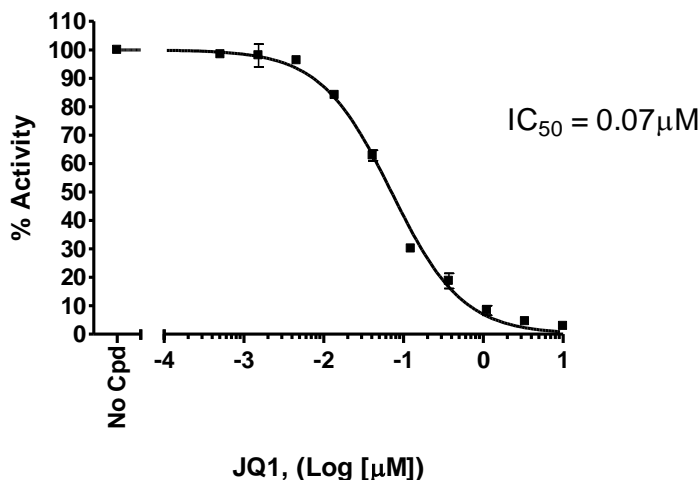
Where $FRET_s$ = Sample FRET, $FRET_{Neg}$ = Negative control FRET, and $FRET_p$ = Positive control FRET.

EXAMPLE OF ASSAY RESULTS:

TR-FRET Assay of BRD4 (BD1+BD2)

Inhibition by JQ1

BRD4 (BD1 + BD2) Activity



Inhibition of BRD4 (BD1 + BD2) by JQ1 (BPS Cat. #27401), measured using the *BRD4 TR-FRET Assay Kit*, BPS Bioscience # 32612. 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 **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@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

RELATED PRODUCTS:

<u>Product</u>	<u>Catalog #</u>	<u>Size</u>
BET Bromodomain Ligand	33000	0.5 mL
Bromodomain Non-acetylated Ligand 1	33005	0.5 mL
BRD1 (561 – 668), His-tag*	31010	100 µg
BRD2 (65 – 187), His-tag*	31022	100 µg
BRD2 (339 – 459), His-tag*	31020	100 µg
BRD2 (65 – 459), His-tag*	31025	100 µg
BRD3 (29 – 145), His-tag*	31030	100 µg
BRD3 (306 – 417), His-tag*	31031	100 µg
BRD4 (49 – 170), His-tag*	31042	100 µg
BRD4 (342 – 460), His-tag*	31043	100 µg
BRD4 (49 – 460), His-tag*	31045	100 µg
BRD1 Inhibitor Screening Kit	32521	384 rxns
BRD2 (BD2) Inhibitor Screening Kit	32522	384 rxns
BRD3 (BD1) Inhibitor Screening Kit	32513	384 rxns
BRD3 (BD2) Inhibitor Screening Kit	32523	384 rxns
BRD4 (BD1) Inhibitor Screening Kit	32514	384 rxns
BRD4 (BD2) Inhibitor Screening Kit	32524	384 rxns
(+)-JQ1 Inhibitor	27401	1 mg

*also available as GST-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 **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: [**info@bpsbioscience.com**](mailto:info@bpsbioscience.com)