



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
ATAD2A TR-FRET Assay Kit
Catalog # 32618

DESCRIPTION: The ATAD2A TR-FRET Assay Kit is designed to measure the inhibition of ATAD2A 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, ATAD2A, substrate, and an inhibitor is incubated for 120 minutes. Then, the fluorescence intensity is measured using a fluorescence reader.

COMPONENTS:

Catalog #	Component	Amount	Storage	
31121	ATAD2A, GST	10 µg	-80 °C	(Avoid freeze/ thaw cycles!)
	BET Bromodomain Ligand	50 µl	-80 °C	
	Non-acetylated Ligand 1	15 µl	-80 °C	
	Tb 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 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.

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

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

Protocol for ATAD2A 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 the same solution without inhibitor (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
Inhibitor Buffer (no inhibitor)	2 µl	2µl	–
BET Bromodomain Ligand		5 µl	5 µl
Non-acetylated Ligand 1	5 µl	–	–
ATAD2A 0.9 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.

- 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.*
- 6) Individually dilute **BET Bromodomain** 40-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** 40-fold in **1x BRD TR-FRET Assay Buffer 1** and add 5 µl of diluted **Non-***
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 Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

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 **ATAD2A** bromodomain protein on ice. Upon first thaw, briefly spin tube containing protein to recover the full contents of the tube. Aliquot **ATAD2A** protein into single-use aliquots. Store remaining undiluted **ATAD2A** in aliquots at -80°C immediately. *Note: ATAD2A is very sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted protein.*
- 8) Dilute **ATAD2A** in **1x BRD TR-FRET Assay Buffer 1** to 0.9 ng/ μ l (2.7 ng/reaction). Initiate reaction by adding 3 μ l of diluted **ATAD2A** to wells designated for the “Negative Control” “Positive Control”, and “Test Inhibitor”. Discard any remaining diluted ATAD2A 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

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.

$$\% \text{ Activity} = \frac{\text{FRET}_s - \text{FRET}_{neg}}{\text{FRET}_p - \text{FRET}_{neg}} \times 100\%$$

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

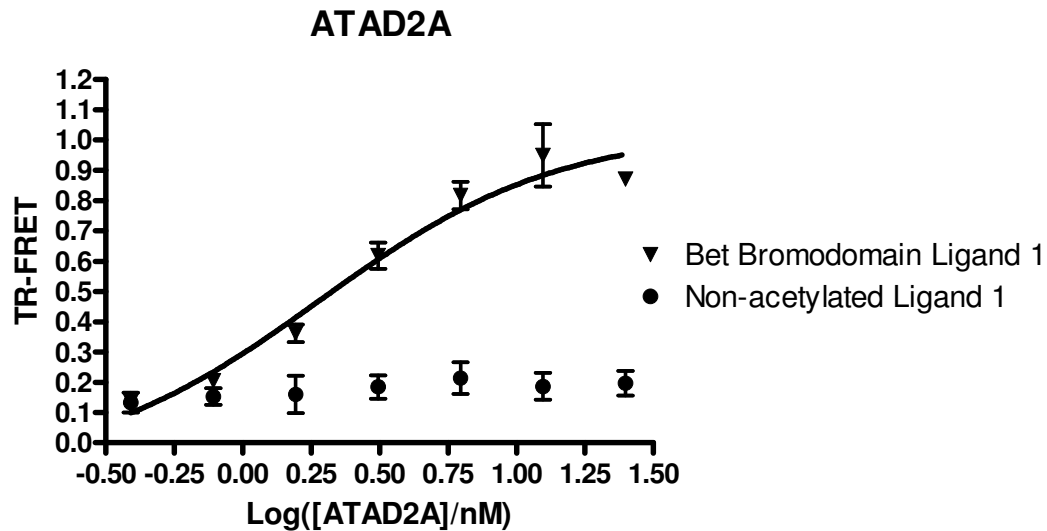
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

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

EXAMPLE OF ASSAY RESULTS:



Interaction of ATAD2A (BPS Bioscience # 31121) with peptide ligands. *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 Court W, 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>
BET Bromodomain Ligand	33000	0.5 mL
Bromodomain Non-acetylated Ligand 1	33005	0.5 mL
ATAD2A (981-1108), GST-tag*	31121	100 µg
ATAD2B (953 – 1080), GST-tag*	31125	100 µg
BRDT, BD1 and BD2 (22-382), GST-tag	31061	100 µg
BRD1 (561 – 668), GST-tag*	31011	100 µg
BRD2 (65 – 459), GST-tag*	31024	100 µg
BRD3 (29 – 417), GST-tag*	31034	100 µg
BRD4 (49 – 460), GST-tag*	31044	100 µg
ATAD2A Inhibitor Screening Assay Kit	32601	384 rxns.
ATAD2B Inhibitor Screening Assay Kit	32605	384 rxns.
BRD2 (BD1+BD2) TR-FRET Assay Kit	32614	384 rxns.
BRD3 (BD1+BD2) TR-FRET Assay Kit	32615	384 rxns.
BRD4 (BD1) TR-FRET Assay Kit	32613	384 rxns.
BRD4 (BD1+BD2) TR-FRET Assay Kit	32612	384 rxns.
(+)-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

Please visit our website at: www.bpsbioscience.com