



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

Data Sheet

3x BRD TR-FRET Assay Buffer 1

Catalog #: 33012

Lot #: 150409

DESCRIPTION: The *3x BRD TR-FRET Assay Buffer 1* is optimized for use with the following BPS Bioscience assay kits:

• ATAD2A TR-FRET Assay Kit	#32618	384 reactions
• ATAD2B TR-FRET Assay Kit	#32620	384 reactions
• BPTF/FALZ TR-FRET Assay Kit	#32632	384 reactions
• BRD1 TR-FRET Assay Kit	#32625	384 reactions
• BRD2 (BD1+BD2) TR-FRET Assay Kit	#32614	384 reactions
• BRD3 (BD1+BD2) TR-FRET Assay Kit	#32615	384 reactions
• BRD4 (BD1+BD2) TR-FRET Assay Kit	#32612	384 reactions
• BRD4 (BD1) TR-FRET Assay Kit	#32613	384 reactions
• BRD4 (BD2) TR-FRET Assay Kit	#32617	384 reactions
• BRD9 TR-FRET Assay Kit	#32621	384 reactions
• BRD9 TR-FRET Assay Kit	#32621	384 reactions
• BRPF3 TR-FRET Assay Kit	#32626	384 reactions
• CECR2 TR-FRET Assay Kit	#32622	384 reactions
• CREBBP TR-FRET Assay Kit	#32619	384 reactions
• TAF1 (BD2) TR-FRET Assay Kit	#32627	384 reactions
• TAF1 (BD1+BD2) TR-FRET Assay Kit	#32628	384 reactions
• TAF1L (BD1+BD2) TR-FRET Assay Kit	#32629	384 reactions
• WDR9/BRWD1 TR-FRET Assay Kit	#32631	384 reactions

APPLICATIONS: Studying enzyme kinetics and HTS applications in drug discovery.

SUPPLIED AS: HEPES-buffer solution.

QUANTITY: 10 ml

STORAGE: -20°C. Avoid freeze/thaw cycles.

STABILITY: Stable for 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



6044 Cornerstone Court West, Suite 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>
(+)-JQ1	27400	10 mg
BRD2 (339 – 459), His-tag	31020	100 µg
BRD2 (65 – 187), GST-tag	31021	100 µg
BRD3 (29 – 145), His-tag	31030	100 µg
BRD3 (29 – 145), GST-tag	31032	100 µg
BRD3 (306 – 417), His-tag	31031	100 µg
BRD3 (306 – 417), GST-tag	31033	100 µg
BRD4 (49 – 170), GST-tag	31040	100 µg
BRD4 (342 – 460), GST-tag	31041	100 µg
BRD4 (49 – 170), His-tag	31042	100 µg
BRD4 (342 – 460), His-tag	31043	100 µg
BRD9 (135 – 242), His-tag	31090	100 µg
BRDT (22 – 138), GST-tag	31108	100 µg
BRDT (22 – 138), His-tag	31101	100 µg
BRDT (257 – 382), His-tag	31100	100 µg
BET Bromodomain Ligand	33000	0.5 mL

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