



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 Homogeneous Assay Buffer 3

Catalog #: 33014

Lot #: 150414

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

- BAZ2B Inhibitor Screening Assay Kit #32600 384 reactions

NOTE: This product was formerly known as *BAZ2B Assay Buffer*.

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.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
(+)-JQ1	27400	10 mg
BAZ2B, His-tag	31113	100 µg
BAZ2B Inhibitor Screening Assay Kit	32600	384 reactions
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
Bromodomain Ligand 2	33003	0.5 ml
Bromodomain Non-Acetylated Ligand 2	33004	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