



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

3x HSP90 Assay Buffer 2

Catalog #: 50324
Lot #: 160810

SIZE: 10 mL

DESCRIPTION: Concentrated aqueous buffer optimized for use with the following BPS Bioscience HSP90 assay kits:

- HSP90 α (C-Terminal) Inhibitor Screening Assay Kit #50317 384 rxns
- HSP90 β (C-Terminal) Inhibitor Screening Kit #50314 384 rxns
- HSP90 α (C-Terminal Domain) TR-FRET Kit #50261 384 rxns
- HSP90 β (C-Terminal Domain) TR-FRET Kit #50262 384 rxns

Note: This buffer is not designed for use with the HSP90 N-Terminal Assay Kits.

APPLICATIONS: Suitable for use with purified HSP90 enzymes.

SUPPLIED AS: Aqueous buffer solution

STORAGE/STABILITY: At least 6 months at -20°C . Avoid freeze/thaw cycles.

RELATED PRODUCTS:

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
HSP90 α recombinant protein	50290	200 μg
HSP90 β recombinant protein	50292	200 μg
HSP90 α Assay Kit (96 well)	50293	96 reactions
HSP90 α Assay Kit (384 well)	50298	384 reactions
HSP90 β Assay Kit (96 well)	50294	96 reactions
HSP90 β Assay Kit (384 well)	50299	384 reactions
Geldanamycin	27008	5 mg
Geldanamycin, FITC-labeled	50312	30 μl

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