



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

## Data Sheet

### **5x HSP90 Assay Buffer 1**

**Catalog #: 50311**  
**Lot #: 160210**

**SIZE:** 20 mL

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

- |                                       |        |               |
|---------------------------------------|--------|---------------|
| • HSP90 $\alpha$ Assay Kit (96 well)  | #50293 | 96 reactions  |
| • HSP90 $\alpha$ Assay Kit (384 well) | #50298 | 384 reactions |
| • HSP90 $\beta$ Assay Kit (96 well)   | #50294 | 96 reactions  |
| • HSP90 $\beta$ Assay Kit (384 well)  | #50299 | 384 reactions |

*Note: This buffer is not designed for use with the HSP90 $\beta$  (C-Terminal) Inhibitor Screening Assay Kit, Cat. # 50314.*

**APPLICATIONS:** Suitable for use with purified HSP90 enzymes.

**SUPPLIED AS:** aqueous buffer solution

**STORAGE/STABILITY:** >6 months at  $-20^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

#### **RELATED PRODUCTS:**

<u>Product Name</u>	<u>Catalog #</u>	<u>Size</u>
HSP90 $\alpha$ recombinant protein	50290	200 $\mu\text{g}$
HSP90 $\beta$ recombinant protein	50292	200 $\mu\text{g}$
HSP90 $\alpha$ Assay Kit (96 well)	50293	96 reactions
HSP90 $\alpha$ Assay Kit (384 well)	50298	384 reactions
HSP90 $\beta$ Assay Kit (96 well)	50294	96 reactions
HSP90 $\beta$ Assay Kit (384 well)	50299	384 reactions
Geldanamycin	27008	5 mg
Geldanamycin, FITC-labeled	50312	30 $\mu\text{l}$

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)