

# HSP Screening Services

Heat shock or stress proteins (HSPs) are functionally related proteins found in unstressed cells. These molecular chaperones play a key role in the folding, unfolding and translocation of other proteins. Expression levels of HSPs increase when cells are exposed to heat or other stressors such as pH shift or hypoxia. Identification and optimization of small molecule inhibitors are ideal for studying cancer vaccines, the discovery of anticancer therapeutics, and the development of infectious disease treatments.

BPS Bioscience offers HSP90 screening services for the evaluation of lead compounds. Our unique assays can differentiate between inhibitors of the ATP-binding region (C-terminus) or the ATPase site (N-terminus) of HSP90. Inhibitors of the C-terminus of HSP90 prevent HSP90 autophosphorylation, while inhibitors of the N-terminus affect chaperoning activity.

HSP Assays	
<ul style="list-style-type: none"> <li>• HSP90<math>\alpha</math> (C-terminal) Inhibitor Screening</li> <li>• HSP90<math>\alpha</math> (N-terminal) Inhibitor Screening</li> <li>• HSP90<math>\beta</math> (C-terminal) Inhibitor Screening</li> <li>• HSP90<math>\beta</math> (N-terminal) Inhibitor Screening</li> </ul>	
HSP Inhibitors	
<ul style="list-style-type: none"> <li>• 17-AAG</li> <li>• BIIB021</li> <li>• Geldanamycin</li> <li>• Geldanamycin, FITC-labeled</li> </ul>	<ul style="list-style-type: none"> <li>• Novobiocin</li> <li>• Novo-TRX</li> <li>• NVP-AUY922</li> </ul>

\*To review sample data for the HSP assays listed above, click on the following link or copy and paste it into your web browser: [http://www.bpsbioscience.com/images/pdf/HSP\\_profiling\\_data.pdf](http://www.bpsbioscience.com/images/pdf/HSP_profiling_data.pdf)

- Fast turn-around time, typically under 2 weeks
- Fully customizable: Determine IC<sub>50</sub> values (10 point, duplicate) or screen a single concentration
- Enzyme activity is validated on every assay to ensure reliable results
- Choice of dose range and concentrations

### Sample Results: HSP90 Activity and Inhibition

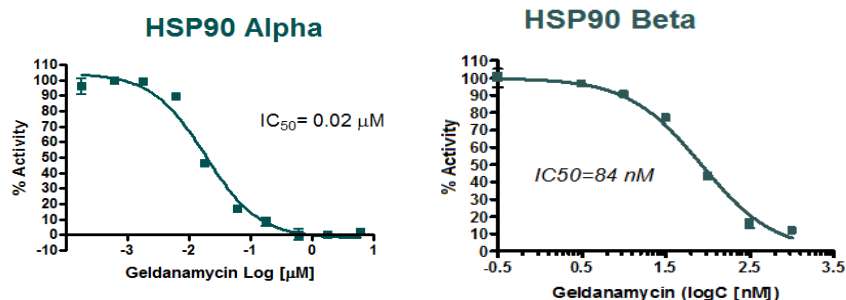


Figure 1. (Left) HSP90 $\alpha$  and (Right) HSP90 $\beta$  inhibition by Geldanamycin, measured using BPS Bioscience's HSP90 screening and profiling services, proteins and inhibitors.

Contact us for a free quotation or information on our HSP Screening Services by visiting:

[www.bpsbioscience.com/contact-us](http://www.bpsbioscience.com/contact-us)

BPS Bioscience, Inc. • 6044 Cornerstone Court West, Suite E, San Diego, CA 92121  
 Tel: 858.829.3082 • Fax: 858.481.8694 • Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com) • Website: [www.bpsbioscience.com](http://www.bpsbioscience.com)