

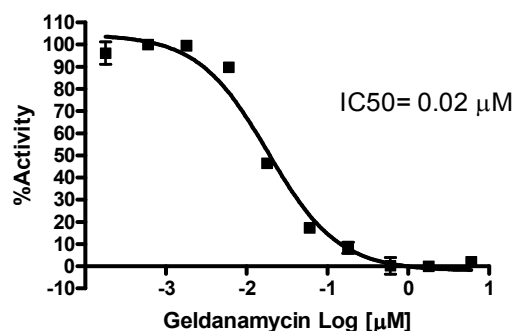
Heat Shock Proteins

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. HSPs are ideal for the study and effectiveness of cancer vaccines, the discovery of anticancer therapeutics, and the development of infectious disease treatments. BPS offers several HSP proteins, assay kits and inhibitors as well as HSP90 screening services.



HSP Assay Kits

The HSP Assay kits are designed for screening for various HSP inhibitors. The kits include all of the reagents necessary to measure HSP inhibition in a convenient 96-well or 384-well format.



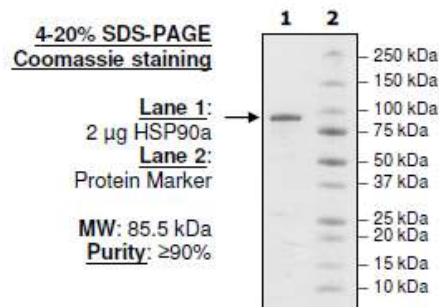
Determination of the IC₅₀ of FITC-labeled Geldanamycin (Cat. #50312) on HSP90α (Cat. #50290), measured with the HSP90α Assay Kit (Cat. #50293/50298).

Assay Kit	Catalog #	Amount
5x HSP90 Assay Buffer	50311	20 ml
HSP90α (N-terminal) Inhibitor Screening Assay Kit	50293	96 rxns
	50298	384 rxns
HSP90α (C-terminal) Inhibitor Screening Assay Kit	50317	384 rxns
HSP90β (N-terminal) Inhibitor Screening Assay Kit	50294	96 rxns
	50299	384 rxns
HSP90β (C-terminal) Inhibitor Screening Assay Kit	50314	384 rxns
HSP90α (C-terminal) TR-FRET Kit	50261	384 rxns
HSP90β (C-terminal) TR-FRET Kit	50262	384 rxns

Heat Shock Proteins

BPS Bioscience offers a wide range of high purity HSPs for drug discovery research.

Protein	Catalog #	Amount
HSP40	50285	100 µg
HSP70	50287	200 µg
HSP70 (Q435A)	50288	100 µg
HSP90α	50290	200 µg
HSP90β	50292	200 µg
HSP90α (C-terminal), Biotin-labeled	50316	100 µg
HSP90β (C-terminal), Biotin-labeled	50313	100 µg
Hop/STIP1	50322	100 µg
SNX-2112	27653-1	5 µg
	27653-2	10 µg
	27653-3	50 µg



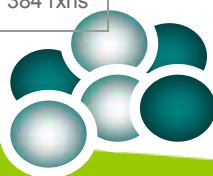
Purity and gel image of HSP90α (Cat. #50290).

HSP Inhibitors

Inhibitor	Catalog #	Amount
17-AAG	27026	5 mg
BIIB02	27046	5 mg
Geldanamycin	27008	5 mg
Geldanamycin, FITC-labeled	50312	30 µl
MS-275 (Entinostat)	27011	25 mg
Novobiocin	27501	250 µl
Novo-TRX	27500	20 µl
NVP-AUY922, HSP90 Inhibitor	27025	10 mg

6044 Cornerstone Court West, Suite E
San Diego, CA 92121
858.829.3082 ph
858.481.8694 fax
www.bpsbioscience.com

Rev. Jan 2015





The #1 Source for Active Enzymes

