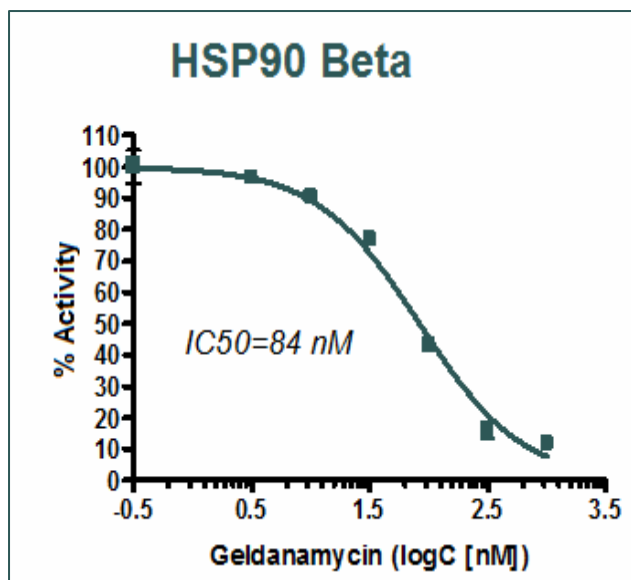
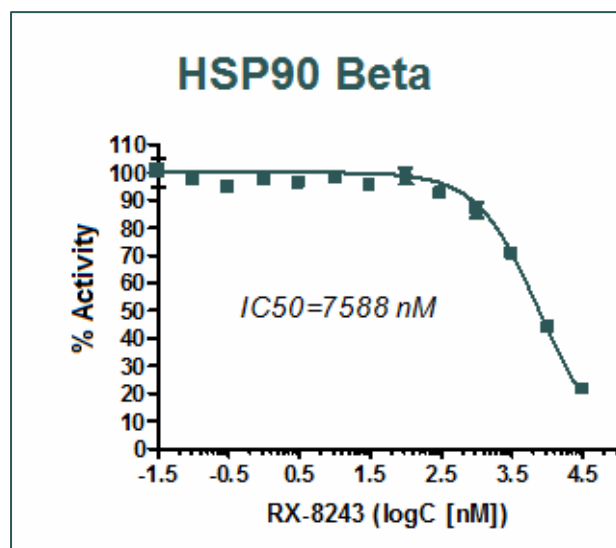
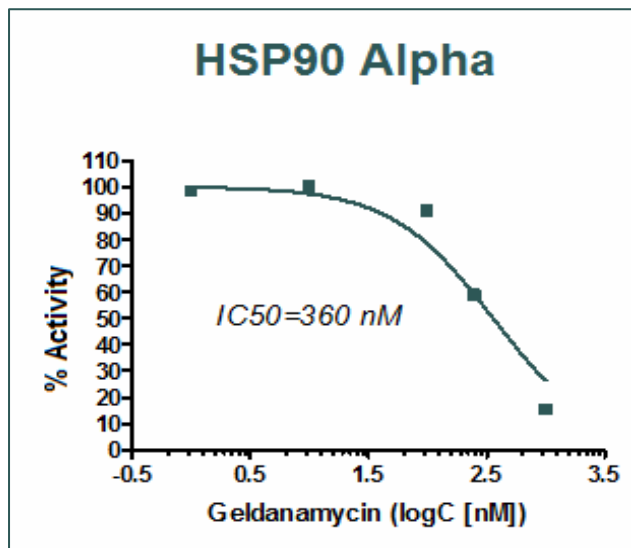


Hsp90 α and Hsp90 β Screening Services



A list of other Screening/Profiling Services offered by BPS:

Histone Methyltransferases
Histone Demethylases
DNA Methyltransferases
HDACs
DPPs
PDEs
PTPs
Pyruvate Kinases
Tumor Cell Proliferation Assay
MMP2 Screening
PARPs
EGFR
Cathepsins

Please contact us for more information on our screening and profiling services.

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