

HSP90b Heat Shock Protein Assay Service

HSP90b [Heat Shock Protein Assay](#) for compound screening and profiling via fluorescence polarization-based quantification of enzyme activity.

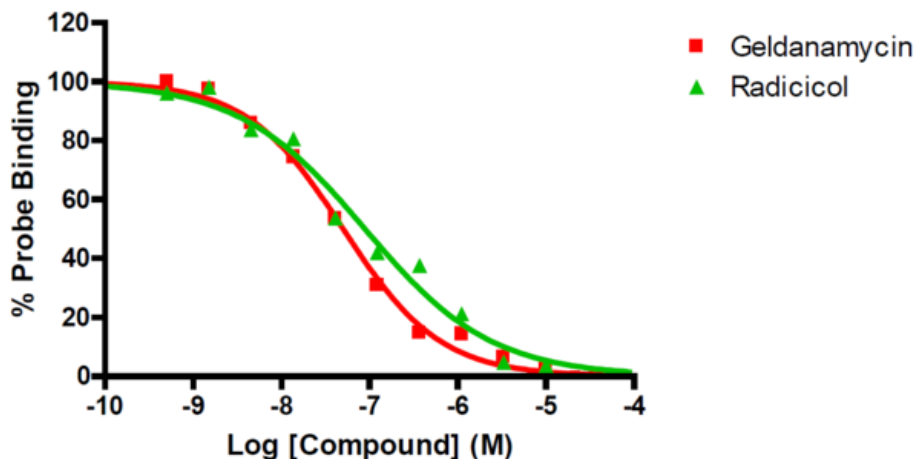
Target Overview

CAT#	HSP90b
Enzyme	Human HSP90b
Synonym	Heat shock protein HSP 90-beta, Heat shock 84 kDa, HSP90AB1, HSP90B, HSPC2, HSPCB, HS90B
Substrate	Geldanamycin, 5 nM

Assay Properties

Readout	Binding competition with fluorescent geldanamycin						
Reference Compound IC50s	<table border="1"> <thead> <tr> <th>Compounds</th> <th>IC50 (nM)</th> </tr> </thead> <tbody> <tr> <td>Geldanamycin</td> <td>50.3</td> </tr> <tr> <td>Radicicol</td> <td>89.0</td> </tr> </tbody> </table>	Compounds	IC50 (nM)	Geldanamycin	50.3	Radicicol	89.0
Compounds	IC50 (nM)						
Geldanamycin	50.3						
Radicicol	89.0						
Scientific Information	P08238 (http://www.uniprot.org/uniprot/P08238)						
Screening Location	Malvern, USA						
Further information	More information can be found on our website Heat Shock Protein Assays .						

Control Compound IC50 for HSP90b



Let's Discover Together.

© Oct 2022 **Reaction Biology**. All rights reserved.

https://www.reactionbiology.com/datasheet/hsp90b_hsp_malvern

