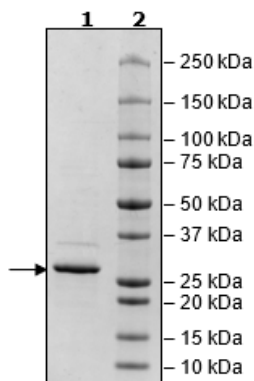


Product Information

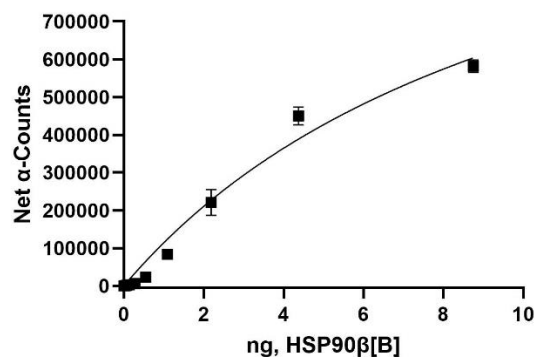
Construct:	HSP90 β (C-terminal) (His-Avi-527-724(end))-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be $\geq 90\%$.
Concentration:	2.57 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C . Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001271969
MW:	25 kDa
Purity:	$\geq 90\%$
Assay Conditions:	Assay was done according to HSP90 β (C-terminal) Inhibitor Screening Kit (BPS Bioscience #50314) with HSP90 β [B] at 35 ng/rxn and PPID-GST at 52.5 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

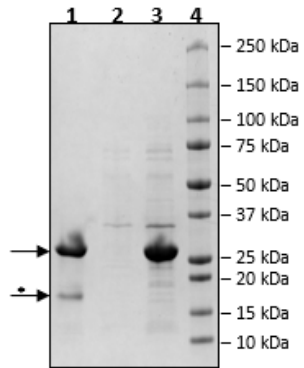
4-20% SDS-PAGE Coomassie Staining



HSP90 β [Biotinylated]: PPID-GST Activity



Biotin-Avidin Pulldown



1. Beads
2. Flow thru
3. Control
4. Standards

* Avidin from beads.