

HSP90 α (C-terminal), Biotin-Labeled, His-Tag, Avi-Tag Recombinant

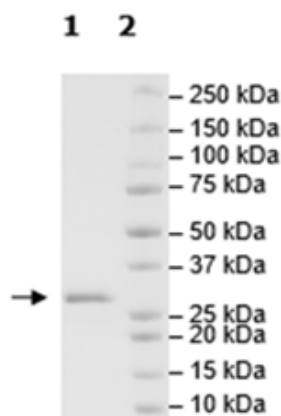
Catalog: 50316
Lot: 210628G

Product Information

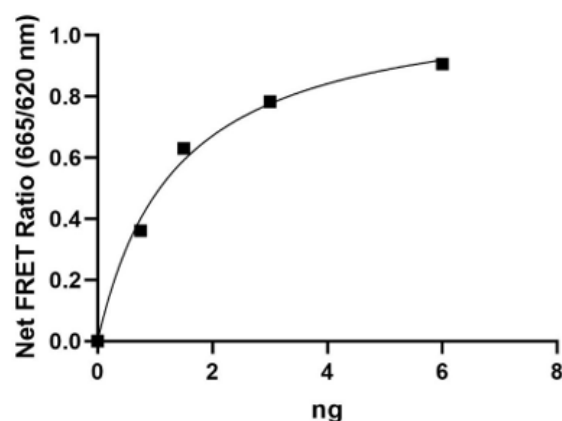
Description:	Recombinant human HSP90 α (Heat shock protein 90) encompassing amino acids 535-732(end). This construct contains an N-terminal His-tag (6xHis) followed by an Avi-Tag™. The recombinant protein was enzymatically biotinylated using the Avi-Tag™ and affinity purified.
Species	Human
Construct:	HSP90 α (His-Avi-535-732(end))-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be $\geq 90\%$.
Concentration:	0.63 mg/ml
Expression System:	<i>E. coli</i>
Purity:	$\geq 90\%$
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol
MW:	25 kDa
Genbank Accession:	NM_005348
Stability:	At least 6 months at -80°C .
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Assay was performed according to the HSP90 α C-Terminal Domain TR-FRET Assay Kit (BPS Bioscience #50289) with HSP90 α titrated from 48 ng/rxn – 0 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 α Activity



Quality Control Data

Biotin-Avidin Pulldown

