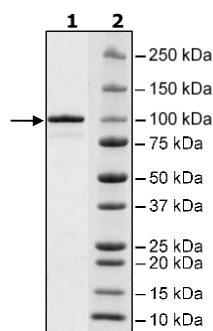


## Product Information

<b>Construct:</b>	HSP90α (1-732(end)-His)
<b>Concentration:</b>	0.87 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 10% glycerol, 3 mM DTT, 225 imidazole
<b>Expression System:</b>	<i>E. coli</i>
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_005348
<b>MW:</b>	86 kDa
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	Assay was performed according to the HSP90α N-Terminal Domain Assay Kit (BPS Bioscience #50293) with HSP90α titrated from 340 ng/rxn – 0 ng/rxn.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining



## HSP90α Activity

