

PRODUCT INFORMATION



Hsp90 α (human recombinant)

Item No. 10202

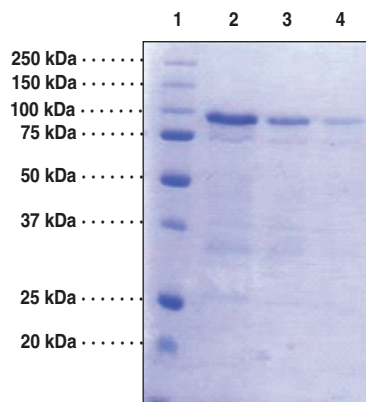
Overview and Properties

Synonyms: Heat Shock Protein 90 α , Hsp86
Source: Recombinant N-terminal hexahistidine-tagged protein, expressed in *E. coli*
Amino Acids: 2-732 (full-length)
Uniprot No.: P07900

Batch specific information can be found on the Certificate of Analysis or by contacting Technical Support

Molecular Weight: 87 kDa
Storage: -80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
Stability: As supplied, 6 months from the QC date provided on the Certificate of Analysis, when stored properly
Purity: *batch specific* ($\geq 80\%$ estimated by SDS-PAGE)
Supplied in: *batch specific*
Protein
Concentration: *batch specific* mg/ml
Additional Information: This protein has not been tested for enzyme activity.

Image



Lane 1: MW Markers
Lane 2: Hsp90 α (4 μ g)
Lane 3: Hsp90 α (2 μ g)
Lane 4: Hsp90 α (1 μ g)

Representative gel image shown; actual purity may vary between each batch but protein will be $\geq 80\%$ pure.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 07/18/2016

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Heat shock proteins (Hsps) are molecular chaperones that can account for upwards of 1-2% of all cellular protein.¹ The Hsp90 gene family consists of Hsp90A, Hsp90B, and TRAP subfamilies. The Hsp90A family is localized in the cytoplasm and can be further sub-divided into Hsp90AA (inducible) and Hsp90AB (constitutively expressed), yielding Hsp90 α and Hsp90 β proteins, respectively. Hsp90B codes for the ER-localized protein Grp94, while TRAP, producing TRAP1, is a mitochondrial chaperone.² Hsp90 acts as a dimer, which can be homo or heterodimers of the α and β isoforms, and functions to bind protein substrates that are unfolded and/or misfolded to assist in folding and prevent aggregation.³ C-terminus dimerization of Hsp90, coupled with the ATPase molecular clamp activity cause a conformational change in the N-terminal nucleotide binding domain thus facilitating substrate binding.⁴ Hsp90 associates with many co-chaperones including p23/Sba1, which help in recruiting substrates to the Hsp90 complex.⁵ Hsp90 works on many different substrates, however its activity seems to be more specific for protein kinases involved in signal transduction, steroid receptors, and cytoskeletal proteins.¹

References

1. Csermely, P., Schnaider, T., Soti, C., *et al.* The 90-kDa molecular chaperone family: Structure, function, and clinical applications. A comprehensive review. *Pharmacol. Ther.* **79**(2), 129-168 (1998).
2. Chen, B., Piel, W.H., Gui, L., *et al.* The HSP90 family of genes in the human genome: Insights into their divergence and evolution. *Genomics* **86**, 627-637 (2005).
3. Fink, A.L. Chaperone-mediated protein folding. *Physiol. Rev.* **79**(2), 425-449 (1999).
4. Prodromou, C., Panaretou, B., Chohan, S., *et al.* The ATPase cycle of Hsp90 drives a molecular 'clamp' via transient dimerization of the N-terminal domains. *EMBO J.* **19**(16), 4383-4392 (2000).
5. Ali, M.M.U., Roe, S.M., Vaughan, C.K., *et al.* Crystal structure of an Hsp90-nucleotide-p23/Sba1 closed chaperone complex. *Nature* **440**, 1013-1017 (2006).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM