

PRODUCT INFORMATION



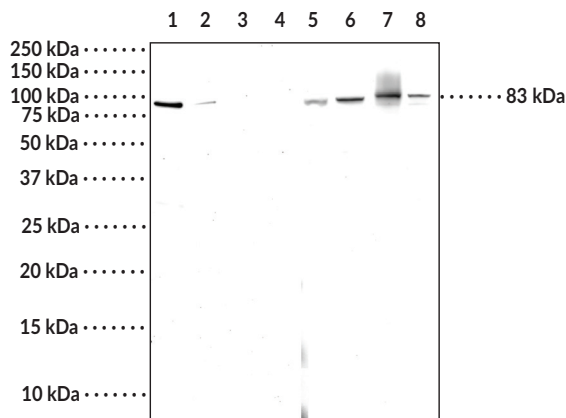
Hsp90 β Monoclonal Antibody (Clone 8D6)

Item No. 25695

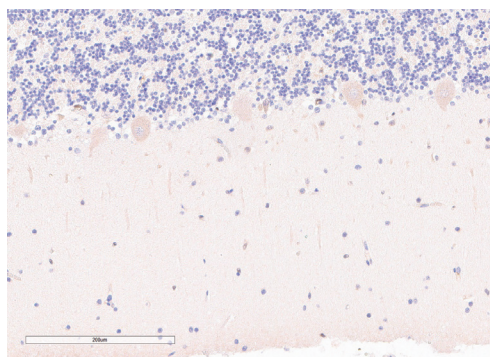
Overview and Properties

Contents:	This vial contains 100 μ g of protein G-purified antibody.
Synonyms:	Heat Shock 84 kDa, Heat Shock Protein Hsp 90-Beta, Hsp84, Hsp90 β
Immunogen:	Full length human recombinant Hsp90 β protein
Cross Reactivity:	(-) HSP90 α
Species Reactivity:	(+) Human, mouse, and rat; other species not tested
Uniprot No.:	P08238
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol, and 0.02% sodium azide
Clone:	8D6
Host:	Mouse
Isotype:	IgG2b
Applications:	ELISA, Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution for ELISA and WB is 1:1,000 and is 1:40 for IHC. Other applications were not tested, therefore optimal working dilution should be determined empirically.

Images



Lane 1: Hsp90 β Recombinant Protein (0.02 μ g)
Lane 2: Hsp90 β Recombinant Protein (0.005 μ g)
Lane 3: Hsp90 β Recombinant Protein (0.001 μ g)
Lane 4: Hsp90 α Recombinant Protein (0.1 μ g)
Lane 5: A549 Cell Lysate (50 μ g)
Lane 6: THP-1 Cell Lysate (50 μ g)
Lane 7: Mouse Kidney Lysate (40 μ g)
Lane 8: Rat Intestine Lysate (40 μ g)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human cerebellum tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with Hsp90 β Monoclonal Antibody (Clone 8D6) (Item No. 25695) at a 1:40 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

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, 12/04/2023

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 protein 90 β (Hsp90 β) is the constitutively active cytosolic isoform of Hsp90 that is encoded by *HSP90AB* in humans.¹ Hsp90 is a multidomain protein that functions as a molecular chaperone to assist in folding and activation of nascent peptides, refolding unfolded or misfolded proteins, and preventing protein aggregation.² C-terminal dimerization of Hsp90, coupled with ATPase molecular clamp activity induces a conformational change in the N-terminal nucleotide binding domain that facilitates substrate binding and initiates the chaperone cycle.³ Hsp90 interacts with many co-chaperones during its chaperone cycle including p23 and Sba1, which help recruit substrates to the Hsp90 complex, Hsp70 (Item Nos. 22739 | 23002), which loads nascent polypeptides onto the Hsp90 dimer, and the ATPase activator Aha1 that promotes ATP hydrolysis and substrate release.^{4,5} Hsp90 is overexpressed in cancer cells and stabilizes client proteins that promote oncogenesis, including transcription factors, signaling proteins, and kinases.^{1,5} Hsp90 also decreases α -synuclein fibril formation and toxicity as well as Q35 aggregation in *in vitro* models of Parkinson's and Huntington's disease, respectively, implying a role in neurodegenerative disease.⁶ Cayman's Hsp90 β Monoclonal Antibody (Clone8D6) can be used for Western blot and ELISA applications. This antibody recognizes Hsp90 β at 83 kDa from human, mouse, and rat samples.

References

1. 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(6)**, 627-637 (2005).
2. Fink, A.L. Chaperone-mediated protein folding. *Physiol. Rev.* **79(2)**, 425-449 (1999).
3. 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).
4. 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(7087)**, 1013-1017 (2006).
5. Li, J. and Buchner, J. Structure, function and regulation of the Hsp90 machinery. *Biomed. J.* **36(3)**, 106-117 (2013).
6. Lackie, R.E., Maciejewski, A., Ostapchenko, V.G., *et al.* The Hsp70/Hsp90 chaperone machinery in neurodegenerative diseases. *Front. Neurosci.* **11:254**, (2017).

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