

Product Information



Hsp90 Complex Monoclonal Antibody (Clone 8D3)

Item No. 10011428

Contents:	This vial contains PEG purified IgG, at a concentration of 1 mg/ml in PBS, pH 7.2, containing 0.09% sodium azide and 50% glycerol.
Synonym:	Heat Shock Protein 90 Complex
Antigen:	Aryl hydrocarbon receptor Hsp90 complex ¹
Host:	Mouse, clone 8D3
Cross Reactivity:	(+) Human, mouse, rat, and rabbit Hsp90. Immunoprecipitates 90 kDa proteins corresponding to the molecular mass of Hsp90. Co-immunoprecipitates Hsp90 complexes, including Hsp70, Hop, aryl hydrocarbon receptors ¹ , glucocorticoid receptors ² , heme-regulated eukaryotic initiation factor 2 α (eIF-2 α) kinase (HRI). ³
Stability:	≥ 1 year at -20°C
Application:	Immunoprecipitation; best results for immunoprecipitation associated with use of goat anti-mouse IgM beads.

Hsp90 is a highly conserved and essential stress protein that is expressed in all eukaryotic cells. From a functional perspective, Hsp90 participates in the folding, assembly, maturation, and stabilization of specific proteins as an integral component of a chaperone complex.⁴⁻⁷ Despite its label of being a Hsp, Hsp90 is one of the most highly expressed proteins in unstressed cells (1-2% cytosolic protein). It carries out a number of housekeeping functions - including controlling the activity, turnover, and trafficking of a variety of proteins. Most of the Hsp90-regulated proteins that have been discovered to date are involved in cell signaling.^{8,9} The number of proteins now known to interact with Hsp90 is about 100. Target proteins include the kinases v-Src, Wee1, and c-Raf, transcriptional regulators such as p53 and steroid receptors, and the polymerases of the hepatitis B virus and telomerase.⁶ When bound to ATP, Hsp90 interacts with co-chaperones Cdc37, p23, and an assortment of immunophilin-like proteins, forming a complex that stabilizes and protects target proteins from proteasomal degradation.

In most cases, Hsp90-interacting proteins have been shown to co-precipitate with Hsp90 when carrying out immunoadsorption studies, and to exist in cytosolic heterocomplexes with it. In a number of cases, variations in Hsp90 expression or Hsp90 mutation has been shown to degrade signaling function *via* the protein or to impair a specific function of the protein (such as steroid binding and kinase activity) *in vivo*. Ansamycin antibiotics, such as geldanamycin and radicicol, inhibit Hsp90 function.¹⁰

References

1. Perdeew, G.H. *J. Biol. Chem.* **263**(27), 13802-13805 (1988).
2. Dalman, F.C., Bresnick, E.H., Patel, P.D., *et al. J. Biol. Chem.* **264**(33), 19815-19821 (1989).
3. Uma, S., Hartson, S.D., Chen, J.-J., *et al. J. Biol. Chem.* **272**(17), 11648-11656 (1997).
4. Arlander, S.J.H., Eapen, A.K., Vroman, B.T., *et al. J. Biol. Chem.* **278**(52), 52572-52577 (2003).
5. Pearl, L.H. and Prodromou, C. *Advances in Protein Chemistry* **59**, 157-172 (2001).
6. Neckers, L. *Trends in Molecular Medicine* **8**(4 Suppl.), S55-S61 (2002).
7. Pratt, W.B. and Toft, D.O. *Exp. Biol. Med.* **228**, 111-133 (2003).
8. Pratt, W.B. and Toft, D.O. *S Endocr. Rev.* **18**(3), 306-360 (1997).
9. Pratt, W.B. The Hsp90-based chaperone system: *Proc. Soc. Exp. Biol. Med.* **217**, 420-434 (1998).
10. Whitesell, L., Mimnaugh, E.G., De Costa, B., *et al. Proc. Natl. Acad. Sci. USA* **91**, 8324-8328 (1994).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10011428

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent *via* email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/13/2011

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com