

Product Information



GRP94 Monoclonal Antibody (Clone 9G10)

Item No. 10011424

Contents:	This vial contains protein G affinity-purified mouse IgG at a concentration of 1 mg/ml in PBS, pH 7.2, containing 0.09% sodium azide and 50% glycerol.
Synonym:	Glucose Regulated Protein, GP96
Antigen:	Purified recombinant GPR94 protein
Host:	Rat, clone 9G10
Isotype:	IgG _{2a}
Cross Reactivity:	(+) Human, mouse, rat, bovine, canine, chicken, guinea pig, hamster, horse, monkey, porcine, rabbit, ovine, and <i>Xenopus</i> GRP94. Detects an 98 kDa protein corresponding to the molecular mass of GRP94 on SDS-PAGE immunoblots. Does not detect human Hsp90, GRP74, or GRPE from <i>E. coli</i> .
Stability:	≥1 year at -20°C
Application:	Western blot, immunoprecipitation, and flow cytometry. ¹⁻⁷ The recommended starting concentration for western blot is 0.5 µg/ml, which was shown to be sufficient for the detection of GRP94 in 20 µg of HeLa lysate. Optimal working dilutions for other applications should be determined empirically.

Glucose regulated protein 94 (GRP94) is a constitutively expressed endoplasmic reticulum (ER) luminal protein that is up-regulated in response to cellular stress such as heat shock, oxidative stress or glucose depletion. GRP94 is thought to play a role in protein translocation to the ER, in their subsequent folding and assembly, and in regulating protein secretion.⁸ GRP94 also plays a role in antigen presentation by accessing the endogenous pathway and eliciting specific cytotoxic T lymphocyte (CTL) responses to chaperone bound peptides *via* the major histocompatibility complex (MHC) class I pathway.⁹

GRP94 is a member of the Hsp90 family of stress proteins and shares sequence homology with its cytosolic equivalent, Hsp90.¹⁰ Both Hsp90 and GRP94 are calcium binding proteins.¹¹ Despite sharing 50% sequence homology over its N domains and complete conservation in its ligand binding domains with Hsp90, GRP94, and Hsp90 differ in their interactions with regulatory ligands as GRP94 has weak ATP binding and hydrolysis activity.¹²

GRP94 exists as a homodimer and the two subunits interact at two distinct intermolecular sites the C-terminal dimerization domains and the N-terminal interacts with the middle domain of opposing subunits.¹³ GRP94 contains a carboxy terminal KDEL (Lys-Asp-Glu-Leu) sequence which is believed to aid in its retention in the ER.¹⁴

References

1. Allen, S., Abuzenadah, A.M., Hinks, J., *et al. Blood* **96**, 560-568 (2000).
2. Sato, K., Torimoto, Y., Tamura, Y., *et al. Blood* **98**, 1852-1857 (2001).
3. Yun, S.-W., Gärtner, U., Arendt, T., *et al. Brain Res. Bull.* **52(5)**, 371-378 (2000).
4. Choukhi, A., Ung, S., Wychowski, C., *et al. J. Virol.* **72(5)**, 3851-3858 (1998).
5. Hoshino, T., Wang, J., Devetten, M.P., *et al. Blood* **91(11)**, 4379-4386 (1998).
6. Riera, M., Roher, N., Miró, F., *et al. Mol. Cell. Biochem.* **191(1-2)**, 97-104 (1999).
7. Gusarova, V., Caplan, A.J., Brodsky, J.L., *et al. J. Biol. Chem.* **276(27)**, 24891-24900 (2001).
8. Ruddon, R.W. and Bedows, E. *J. Biol. Chem.* **272(6)**, 3125-3128 (1997).
9. Srivastava, P.K., Udono, H., Blachere, N.E., *et al. Immunogenetics* **39(2)**, 93-98 (1994).
10. Mazzarella, R.A. and Green, M. *J. Biol. Chem.* **262(18)**, 8875-8883 (1987).
11. Kang, H.S. and Welch, W.J. *J. Biol. Chem.* **266(9)**, 5643-5649 (1991).
12. Soldano, K.L., Jivan, A., Nicchitta, C.V., *et al. J. Biol. Chem.* **278(48)**, 48330-48338 (2003).
13. Chu, F., Maynard, J.C., Chiosis, G., *et al. Protein Science* **15**, 1260-1269 (2006).
14. Peter, F., Van, P.N., and Söling, H.-D. *J. Biol. Chem.* **267(15)**, 10631-10637 (1992).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10011424

WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG 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 under separate cover to the MSDS supervisor at 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, 12/14/2011

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