

Product Information

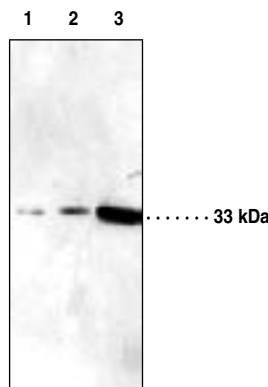


IGFBP5 Polyclonal Antibody

Item No. 10008207

- Synonyms:** Insulin-like Growth Factor Binding Protein 5, IBP-5, IGF-Binding Protein 5
- Contents:** This vial contains 200 µg of peptide affinity-purified IgG in 500 µl of TBS, pH 7.4, containing 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide.
- Host:** Rabbit
- Antigen:** Human IGFBP5 amino acids 192-206 (CRRHMEASLQELKAS)
- Cross-reactivity:** (+) Mouse and rat IGFBP5; other species not tested
Homology of peptide antigen with IGFBP5 from other species:
Human C R R H M E A S L Q E L K A S
Mouse C R R H M E A S L Q E f K A S
Rat C R R H M E A S L Q E f K A S
- Stability:** ≥1 year at -20°C
- Applications:** Recommended starting dilution for western blot is 1:200 (2 µg/ml). Other applications were not optimized so working dilutions should be determined empirically.

Insulin-like growth factor binding protein 5 (IGFBP5) is a secreted protein that binds IGF-1 and restricts it from accessing its cell-surface receptor (IGF-1R). This aids in regulation of cell growth, differentiation, and apoptosis.¹ IGFBP5 also displays IGF-1 independent transactivational activity in the nucleus.² Some tissues and cell types with higher levels of this protein include kidney, heart, placenta, and skeletal cells.³ Expression has also been reported in a variety of cancers, including mammary gland and neuroblastoma.^{1,4} IGFBP5 (272 amino acids) has a calculated molecular weight of 30.6 kDa. Reported glycosylation and phosphorylation sites may explain the band migration to 33 kDa.^{1,5} Cayman's IGFBP5 polyclonal antibody is directed against a synthetic peptide sequence that partially overlaps the IGFBP5 region identified as a nuclear localization signal.²



Lane 1: Murine skeletal muscle supernatant (25 µg)
Lane 2: Murine kidney supernatant (25 µg)
Lane 3: Rat heart supernatant (25 µg)

References

1. Beattie, J., Allan, G.J., Lochrie, J.D., *et al.* Insulin-like growth factor-binding protein-5 (IGFBP-5): A critical member of the IGF axis. *Biochem J.* **395**, 1-19 (2006).
2. Zhao, Y., Yin, P., Bach, L.A., *et al.* Several acidic amino acids in the N-domain of insulin-like growth factor-binding protein-5 are important for its transactivation activity. *J. Biol. Chem.* **281**(20), 14184-14191 (2006).
3. Shimasaki, S., Shimonaka, M., Zhang, H.-P., *et al.* Identification of five different insulin-like growth factor binding proteins (IGFBPs) from adult rat serum and molecular cloning of a novel IGFBP-5 in rat and human. *J. Biol. Chem.* **266**(16), 10646-10653 (1991).
4. Tanno, B., Negroni, A., Vitali, R., *et al.* Expression of insulin-like growth factor-binding protein 5 in neuroblastoma cells is regulated at the transcriptional level by c-Myb and B-Myb *via* direct and indirect mechanisms. *J. Biol. Chem.* **277**(26), 23172-23180 (2002).
5. Pampusch, M.S., Xi, G., Kamanga-Sollo, E., *et al.* Production of recombinant porcine IGF-binding protein-5 and its effect on proliferation of porcine embryonic myoblast cultures in the presence and absence of IGF-I and Long-R3-IGF-I. *J. Endocrinol.* **185**, 197-206 (2005).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10008207

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, 12/12/2011

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