

PRODUCT INFORMATION



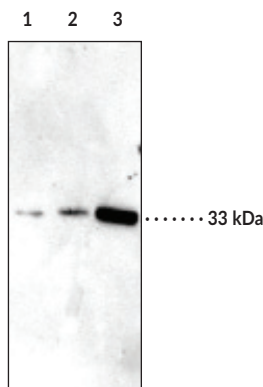
IGFBP5 Polyclonal Antibody

Item No. 10008207

Overview and Properties

Contents:	This vial contains peptide affinity-purified antibody lyophilized from 500 μ l.
Synonyms:	Insulin-like Growth Factor Binding Protein 5, IBP-5, IGF-Binding Protein 5
Immunogen:	Synthetic peptide from the internal region of human IGFBP5
Species Reactivity:	(+) Mouse and rat; other species not tested
Uniprot No.:	P24593
Form:	Solid
Storage:	-20°C (as supplied)
Stability:	\geq 3 years
Storage Buffer:	TBS, pH 7.4, when reconstituted in in 500 μ l double distilled water.
Host:	Rabbit
Application:	Western blot (WB); the recommended starting dilution 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Mouse skeletal muscle supernatant (25 μ g)

Lane 2: Mouse kidney supernatant (25 μ g)

Lane 3: Rat heart supernatant (25 μ g)

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, 10/17/2023

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

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} 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 IGFBP5 (272 amino acids) has a calculated molecular weight of 30.6 kDa. Cayman's IGFBP5 Polyclonal Antibody can be used for Western blot applications and recognizes IGFBP5 at 33 kDa from mouse and rat samples. Reported glycosylation and phosphorylation sites may explain the band migration to 33 kDa.^{1,5}

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).

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