

PRODUCT INFORMATION



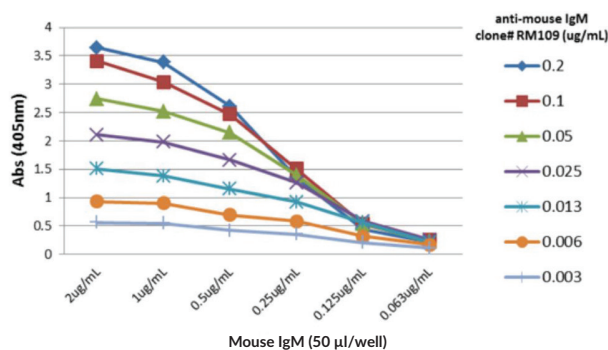
IgM (mouse) Rabbit Monoclonal Antibody (Clone RM109)

Item No. 32101

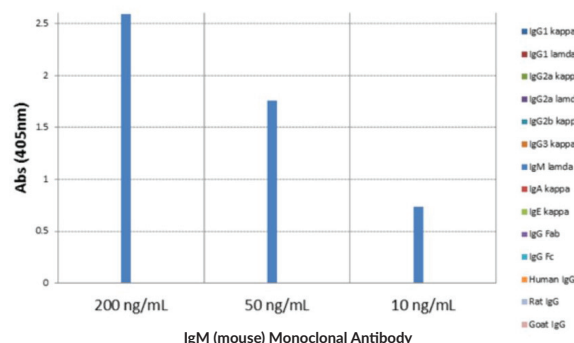
Overview and Properties

Contents: This vial contains 100 µg of protein A-purified monoclonal antibody.
Synonym: Immunoglobulin M
Immunogen: Mouse IgM
Cross Reactivity: (-) Mouse IgG1, IgG2a, IgG2b, IgG3, IgA, IgE; (-) Human, rat, goat IgG
Species Reactivity: (+) Mouse; other species not tested
Form: Liquid
Storage: -20°C (as supplied)
Stability: ≥1 year
Storage Buffer: PBS, with 50% glycerol, 1% BSA, and 0.09% sodium azide
Concentration: 1.0 mg/ml
Clone: RM109
Host: Rabbit
Isotype: IgG
Applications: ELISA, Flow cytometry (FC), and Immunoprecipitation (IP); the recommended starting concentration is 0.005-0.2 µg/ml for ELISA. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



A Titer ELISA of Mouse IgM. The plate was coated with different amounts of mouse IgM. A serial dilution of IgM (mouse) Rabbit Monoclonal Antibody was used as the primary antibody. An alkaline phosphatase-conjugated anti-rabbit IgG was used as the secondary antibody.



ELISA of Mouse Immunoglobulins (Igs). IgM (mouse) Rabbit Monoclonal Antibody reacts to mouse IgM. No cross reactivity with IgG1, IgG2a, IgG2b, IgG3, IgA, IgE, human IgG, rat IgG, or goat IgG. The plate was coated with 50 ng/well of different Igs. 200, 50, or 10 ng/ml of IgM (mouse) Rabbit Monoclonal Antibody was used as the primary antibody. An alkaline phosphatase-conjugated anti-rabbit IgG was used as the secondary antibody.

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, 11/08/2023

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

Immunoglobulin M (IgM) is a member of the immunoglobulin superfamily of glycoproteins that plays a central role in the adaptive immune response and in mucosal immunology.^{1,2} IgM consists of two light chains of approximately 25 kDa each, as well as two heavy chains of approximately 70 kDa each that contain C-terminal extensions, known as tailpieces, which allow for IgM oligomerization.^{2,3} The heavy chains are linked together by disulfide bonds to form an Fc region and also combine with the light chains to form the Fab region, which mediate receptor and antigen binding, respectively.⁴ Five IgM proteins oligomerize *via* disulfide bonds in the presence of a 15-kDa joining chain, a process that is required for transcytosis of IgM from plasma cells to mucosal epithelial cells *via* the polyimmunoglobulin receptor (pIgR).² Monomeric and oligomeric IgM are both ligands for the IgM and IgA-binding high affinity Ig α and the Ig μ Fc receptor (Fc α / μ -R) on dendritic cells, which mediates cellular uptake of IgM-conjugated antigens, and the IgM-binding Fc μ -R on B and T cells, which is important for B cell maturation among other functions.^{2,5-7} IgM is produced primarily in the plasma by naïve B cells and expressed in its monomeric low-affinity form on the cell surface.¹ Activated B cells secrete pentameric high-affinity IgM, which opsonizes antigens to target them for removal by phagocytes and to activate complement *via* the classical pathway.^{1,8} IgM antibodies are produced early in infection and have been used to determine exposure to a specific pathogen.¹ IgM levels are elevated in hyper-IgM syndromes, which are characterized by dysfunctions in Ig class switching recombination.⁹ Cayman's IgM (mouse) Monoclonal Antibody (Clone RM109) can be used for ELISA, flow cytometry (FC), and immunoprecipitation (IP) applications.

References

1. Schroeder, H.W., Jr. and Cavicini, L. Structure and function of immunoglobulins. *J. Allergy Clin. Immunol.* **125(2 Suppl. 2)**, S41-S52 (2010).
2. Li, Y., Wang, G., Li, N., *et al.* Structural insights into immunoglobulin M. *Science* **367(6481)**, 1014-1017 (2020).
3. Jacobin, M.-J., Santarelli, X., Laroche-Traineau, J., *et al.* Production of a human monoclonal IgM directed against human cardiac myosin in a hollow-fiber bioreactor for membrane anion exchange chromatography one-step purification. *Hum. Antibodies* **13(3)**, 69-79 (2004).
4. Vaillant A.A.J. and Ramphul K. Immunoglobulin. In: StatPearls [Internet]. Treasure Island (FL): StatPearls Publishing (2020). Available from: <https://www.ncbi.nlm.nih.gov/books/NBK513460/>
5. Shibuya, A., Sakamoto, N., Shimizu, Y., *et al.* Fc α / μ receptor mediates endocytosis of IgM-coated microbes. *Nat. Immunol.* **1(5)**, 441-446 (2000).
6. Kubagawa, H., Oka, S., Kubagawa, Y., *et al.* Identity of the elusive IgM Fc receptor (Fc μ R) in humans. *J. Exp. Med.* **206(12)**, 2779-2793 (2009).
7. Liu, J., Wang, Y., Min, Q., *et al.* Regulation of humoral immune responses and B cell tolerance by the IgM Fc receptor (Fc μ R). *B Cells in Immunity and Tolerance*. Wang, J.-Y., editor, 1st edition, Springer Nature (2020).
8. Janeway, C.A., Jr., Travers, P., Walport, M., *et al.* Antibodies of different isotype operate in distinct places and have distinct effector functions. *Immunobiology: The Immune System in Health and Disease*. Janeway, C.A., Jr., editor, 5th edition, Garland Science Publishing (2001).
9. Davies, E.G. and Thrasher, A.J. Update on the hyper immunoglobulin M syndromes. *Br. J. Haematol.* **149(2)**, 167-180 (2010).

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