

# PRODUCT INFORMATION



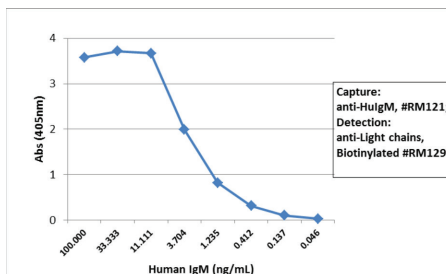
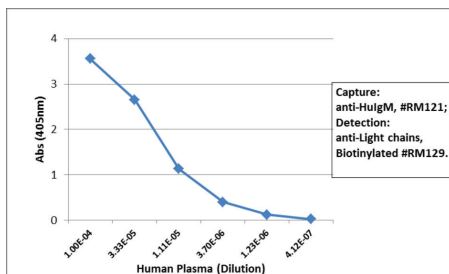
## IgM (human) Monoclonal Antibody (Clone RM121)

Item No. 32113

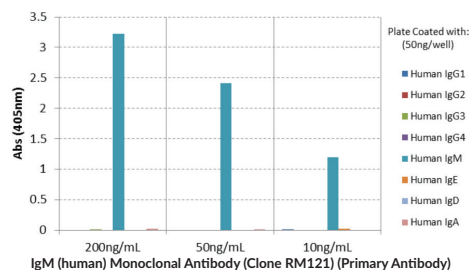
### Overview and Properties

**Contents:** This vial contains 100 µg of protein A-affinity purified monoclonal antibody.  
**Synonym:** Immunoglobulin M  
**Immunogen:** Human IgM  
**Cross Reactivity:** (+) Human IgM; (-) Human IgA, IgD, IgE, IgG  
**Species Reactivity:** (+) Human  
**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 mg/ml  
**Clone:** RM121  
**Host:** Rabbit  
**Isotype:** IgG  
**Applications:** ELISA, immunocytochemistry (ICC), and immunohistochemistry (IHC); the recommended starting concentration for ELISA is 50-200 ng/well (for capture) and 0.05-0.2 µg/ml (for detection), 0.5-2 µg/ml for ICC, and 0.5-2 µg/ml for IHC. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

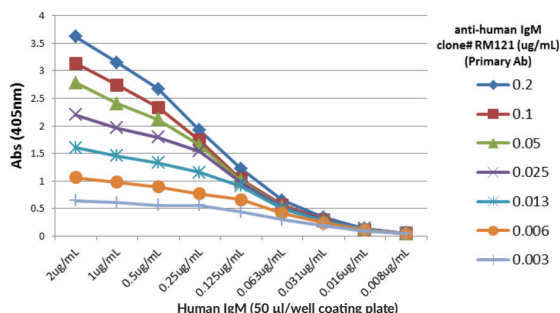
### Images



Sandwich ELISA using IgM (human) Monoclonal Antibody (Clone RM121) as the capture antibody (100 ng/well). Ig Light Chain (human) Monoclonal Antibody - Biotinylated (Item No. 32112) was used as the detection antibody, followed by alkaline phosphatase-conjugated streptavidin.



ELISA of human immunoglobulins (Igs). IgM (human) Monoclonal Antibody (Clone RM121) reacts only to human IgM and not to human IgG, IgE, IgD, or IgA. The plate was coated with 50 ng/well of different Igs. 200, 50 ng/ml of IgM (human) Monoclonal Antibody (Clone RM121) was used as the primary antibody. An alkaline phosphatase-conjugated anti-rabbit IgG was used as the secondary antibody.



A titer ELISA using IgM (human) Monoclonal Antibody (Clone RM121). The plate was coated with different amounts of human IgM. A serial dilution of IgM (human) Monoclonal Antibody (Clone RM121) 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, 11/09/2023

**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.<sup>1,2</sup> 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.<sup>2,3</sup> 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.<sup>4</sup> 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).<sup>2</sup> Monomeric and oligomeric IgM are both ligands for the IgM and IgA-binding high affinity Ig $\alpha$  and the Ig $\mu$  Fc receptor (Fc $\alpha$ / $\mu$ -R) on dendritic cells, which mediates cellular uptake of IgM-conjugated antigens, and the IgM-binding Fc $\mu$ -R on B and T cells, which is important for B cell maturation among other functions.<sup>2,5-7</sup> IgM is produced primarily in the plasma by naïve B cells and expressed in its monomeric low-affinity form on the cell surface.<sup>1</sup> 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.<sup>1,8</sup> IgM antibodies are produced early in infection and have been used to determine exposure to a specific pathogen.<sup>1</sup> IgM levels are elevated in hyper-IgM syndromes, which are characterized by dysfunctions in Ig class switching recombination.<sup>9</sup> Cayman's IgM (human) Monoclonal Antibody (Clone RM121) can be used for ELISA, immunocytochemistry (ICC), and immunohistochemistry (IHC) 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 $\alpha$ / $\mu$  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 $\mu$ 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 $\mu$ R). *B Cells in Immunity and Tolerance*. Wang, J.-Y., editor, 1<sup>st</sup> 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, 5<sup>th</sup> 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