

PRODUCT INFORMATION



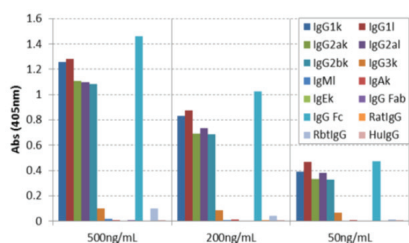
IgG Fc (mouse) Monoclonal Antibody - Biotinylated (RMG06)

Item No. 32362

Overview and Properties

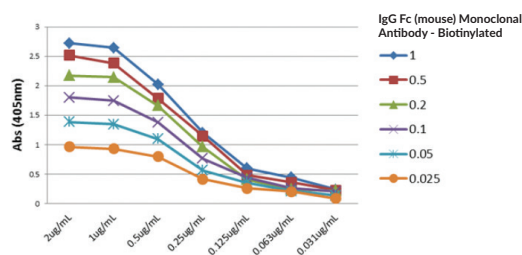
Contents:	This vial contains 50 µg of protein A-affinity purified monoclonal antibody.
Synonym:	Immunoglobulin G
Immunogen:	Mouse IgG
Cross Reactivity:	(+) IgG1, IgG2a, IgG2b, IgG3 (weakly); (-) IgM, IgA, IgE; (-) Human, rabbit, rat IgG
Species Reactivity:	(+) Mouse
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RMG06
Host:	Goat
Isotype:	IgG
Applications:	ELISA; the recommended starting concentration is 0.05-1 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



IgG Fc (mouse) Monoclonal Antibody - Biotinylated (RMG06)

ELISA of mouse immunoglobulins (Igs). IgG Fc (mouse) Monoclonal Antibody - Biotinylated (RMG06) reacts to the Fc region of mouse IgG1, IgG1k, IgG2ak, IgG2bk, and slightly to IgG3k, rat IgG, and rabbit IgG. No cross reactivity with IgM, IgE, IgG Fc, IgG1λ, IgG2aλ, IgAk, IgG Fab, and human IgG. The plate was coated with 50 ng/well of different immunoglobulins. 500 ng/ml, 200 ng/ml, or 50 ng/ml of IgG Fc (mouse) Monoclonal Antibody - Biotinylated (RMG06) was used as the primary antibody. An alkaline phosphatase-conjugated anti-rabbit IgG was used as the secondary antibody.



IgG Fc (mouse) Monoclonal Antibody - Biotinylated (RMG06)

A titer ELISA using IgG Fc (mouse) Monoclonal Antibody - Biotinylated (RMG06). The plate was coated with different amounts of mouse IgG1. A serial dilution of IgG Fc (mouse) Monoclonal Antibody - Biotinylated (RMG06) 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, 02/29/2024

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 G (IgG) is a member of the immunoglobulin superfamily of glycoproteins that plays a central role in the adaptive immune response.¹ It is produced by B cells and later secreted by plasma cells and is the most abundant circulating antibody in human and mouse serum.¹⁻³ IgG consists of two heavy chains of approximately 50 kDa each and two light chains of approximately 25 kDa each.¹ 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 IgG is produced following IgM class-switching in response to infection and is involved in numerous humoral host defense responses, including antibody-dependent cell-mediated cytotoxicity (ADCC), toxin neutralization, and pathogen opsonization.² IgG exists as four isotypes in mice: IgG1, IgG2b, IgG3, and, in a strain-specific manner, IgG2a or IgG2c.^{5,6} Formulations containing humanized, chimeric, or murine IgG monoclonal antibodies have been used in the treatment of inflammatory diseases, such as ulcerative colitis, rheumatoid arthritis, and asthma, as well as cancer.⁶ Cayman's IgG Fc (mouse) Monoclonal Antibody - Biotinylated (RMG06) can be used for ELISA. The antibody recognizes the Fc region of IgG from mouse samples.

References

1. Schroeder, H.W., Jr. and Cavicini, L. Structure and function of immunoglobulins. *J. Allergy. Clin. Immunol.* **125(2 Suppl. 2)**, S41-52 (2010).
2. Vidarsson, G., Dekkers, G., and Rispens, T. IgG subclasses and allotypes: From structure to effector functions. *Front. Immunol.* **5**, 520 (2014).
3. Mayumi, M., Kuritani, T., Kubagawa, H.M., *et al.* IgG subclass expression by human B lymphocytes and plasma cells: B lymphocytes precommitted to IgG subclass can be preferentially induced by polyclonal mitogens with T cell help. *J. Immunol.* **130(2)**, 671-677 (1983).
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. Collins, A.M. IgG subclass co-expression brings harmony to the quartet model of murine IgG function. *Immunol. Cell Biol.* **94(10)**, 949-954 (2016).
6. Martin, R.M., Brady, J.L., and Lew, A.M. The need for IgG2c specific antiserum when isotyping antibodies from C57BL/6 and NOD mice. *J. Immunol. Methods.* **212(2)**, 187-192 (1998).
7. Lu, R.-M., Hwang, Y.-C., Liu, I.-J., *et al.* Development of therapeutic antibodies for the treatment of diseases. *J. Biomed. Sci.* **27(1)**, 1 (2020).

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