

PRODUCT INFORMATION



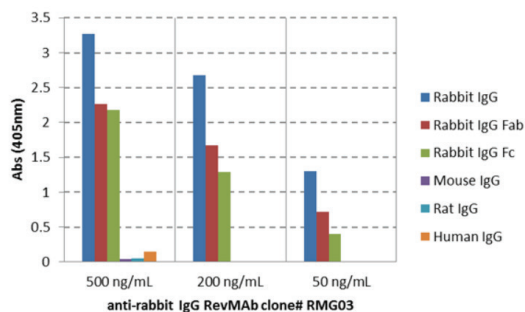
IgG (rabbit) Monoclonal Antibody - Biotinylated (RMG03)

Item No. 32360

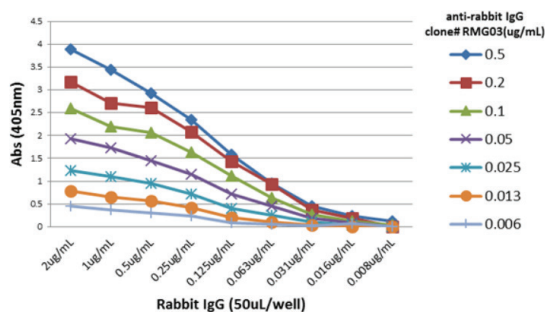
Overview and Properties

Contents:	This vial contains 50 µg of protein G-affinity purified monoclonal antibody.
Synonym:	Immunoglobulin G
Immunogen:	Rabbit IgG
Cross Reactivity:	(+) IgG; (-) Human, mouse, rat IgG
Species Reactivity:	(+) Rabbit
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:	RMG03
Host:	Goat
Isotype:	IgG
Applications:	ELISA and Western blot (WB); the recommended starting concentration is 0.01-0.5 µg/ml for ELISA and 0.1-0.5 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



ELISA of IgGs from different species. IgG (rabbit) Monoclonal Antibody - Biotinylated (RMG03) reacts only to rabbit IgG and not to human IgG, rat IgG, or mouse IgG. The plate was coated with 50 ng/well of different IgGs. IgG (rabbit) Monoclonal Antibody - Biotinylated (RMG03) was used as the primary antibody and an alkaline phosphatase-conjugated anti-goat IgG was used as the secondary antibody.



A Titer ELISA using IgG (rabbit) Monoclonal Antibody - Biotinylated (RMG03). The plate was coated with different amounts of rabbit IgG. A serial dilution of IgG (rabbit) Monoclonal Antibody - Biotinylated (RMG03) was used as the primary antibody and an alkaline phosphatase-conjugated anti-goat 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, 02/17/2024

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.^{1,2} It is produced by B cells and later secreted by plasma cells and is the most abundant circulating antibody in rabbit serum. IgG consists of two identical heavy chains of approximately 50 kDa each and two identical light chains of approximately 25 kDa each.¹ The heavy chains are linked together by a single disulfide bond to form an Fc region and also combine with the light chains through additional disulfide bonds to form the Fab region, which mediate receptor and antigen binding, respectively. 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.^{1,3,5} Rabbits express three light chains, IgG_K (K1), IgG_κ (K2), and IgG_γ, with IgG_K (K1)-containing IgGs comprising approximately 90% of the total IgG population. Unlike human and mouse IgG, rabbit IgG has a single subclass and a short upper and middle hinge length of only 11 residues that facilitates the hinge extension necessary for binding of the Fc receptor (FcR) and activating complement component 1q (C1q) to facilitate complement activation.^{2,4} Cayman's IgG (rabbit) Monoclonal Antibody - Biotinylated (RMG03) can be used for ELISA and Western blot (WB) applications.

References

1. Weber, J., Peng, H., and Rader, C. From rabbit antibody repertoires to rabbit monoclonal antibodies. *Exp. Mol. Med.* **49(3)**, e305 (2017).
2. Rayner, L.E., Kadkhodayi-Kholghi, N., Heenan, R.K., *et al.* The solution structure of rabbit IgG accounts for its interactions with the Fc receptor and complement C1q and its conformational stability. *J. Mol. Biol.* **425(3)**, 506-523 (2013).
3. Vidarsson, G., Dekkers, G., and Rispens, T. IgG subclasses and allotypes: From structure to effector functions. *Front. Immunol.* **5**, 520 (2014).
4. Allan, R. and Isliker, H. Studies on the complement-binding site of rabbit immunoglobulin G-11. The reaction of rabbit IgG and its fragments with Clq. *Immunochemistry* **11(5)**, 243-248 (1974).
5. Williams, R.C., Jr., Osterland, C.K., Margherita, S., *et al.* Studies of biologic and serologic activities of rabbit-IgG antibody depleted of carbohydrate residues. *J. Immunol.* **111(6)**, 1690-1698 (1973).

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