

PRODUCT INFORMATION



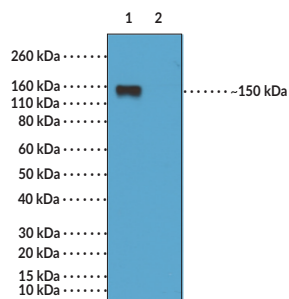
IgG2a (mouse) Monoclonal Antibody (Clone RM219)

Item No. 32005

Overview and Properties

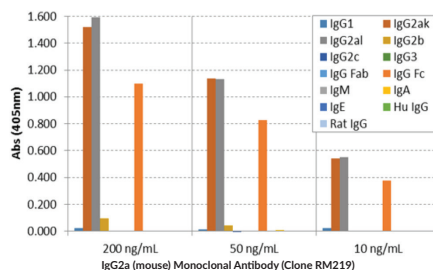
Contents:	This vial contains 100 µg of protein A-affinity purified monoclonal antibody.
Synonym:	Immunoglobulin G2a
Immunogen:	Mouse IgG2a
Cross Reactivity:	(-) Mouse IgG1, IgG2b, IgG2c, IgG3, IgM, IgA, IgE; (-) Human, rat IgG
Species Reactivity:	(+) Mouse IgG2a Fc region
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Concentration:	1.0 mg/ml
Clone:	RM219
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA and Western blot (WB; non-reduced); the recommended starting concentration for ELISA is 0.005-0.2 µg/ml and 0.5-2 µg/ml for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images

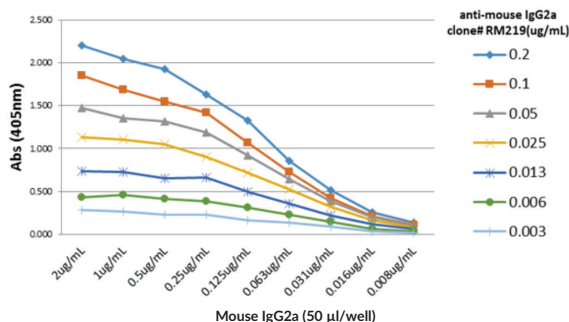


Lane 1: Non-reduced mouse IgG2a (0.5 µg/ml)
Lane 2: Reduced mouse IgG2a (0.5 µg/ml)

WB of non-reduced and reduced mouse IgG2a using 0.5 µg/ml of IgG2a (mouse) Monoclonal Antibody (Clone RM219). This antibody reacts to non-reduced IgG2a (~150 kDa).



ELISA of Mouse Immunoglobulins (Igs). IgG2a (mouse) Monoclonal Antibody (Clone RM219) reacts to the Fc region of mouse IgG2a. No cross reactivity with IgG1, IgG2b, IgG2c, IgG3, IgM, IgA, IgE, human IgG, or rat IgG. The plate was coated with 50 ng/well of different Igs. 200, 50, or 10 ng/ml of IgG2a (mouse) Monoclonal Antibody (Clone RM219) was used as the primary antibody. An alkaline phosphatase conjugated anti-rabbit IgG was used as the secondary antibody.



A Titer ELISA of Mouse IgG2a. The plate was coated with different amounts of mouse IgG2a. A serial dilution of IgG2a (mouse) Monoclonal Antibody (Clone RM219) 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 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} *In vivo*, class switching to the IgG2a isotype can happen via IFN- γ -dependent and -independent mechanisms, with the former resulting from the cognate interaction of B cells with T helper 1 (Th1) cells.⁷ IgG2a is the predominant isotype produced in response to infection with DNA or RNA viruses in mice.⁸ Cayman's IgG2a (mouse) Monoclonal Antibody (Clone RM219) can be used for ELISA and Western blot (WB; non-reducing conditions) applications. The antibody recognizes the Fc region of IgG2a from mouse samples at approximately 150 kDa.

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. Vidarsson, G., Dekkers, G., and Rispen, 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. Jegerlehner, A., Maurer, P., Bessa, J.M., *et al.* TLR9 Signaling in B Cells Determines Class Switch Recombination to IgG2a. *J. Immunol.* **178**(4), 2415-2420 (2007).
8. Coutlelier, J.P., Van Der Logt, J.T.M., Heessen, F.W.A., *et al.* IgG2a restriction of murine antibodies elicited by viral infections. *J. Exp. Med.* **165**(1), 64-69 (1987).

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