

# PRODUCT INFORMATION

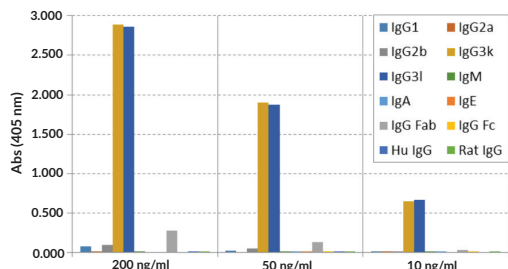


## IgG3 (mouse) Rabbit Monoclonal Antibody - Biotinylated (RM218) Item No. 32354

### Overview and Properties

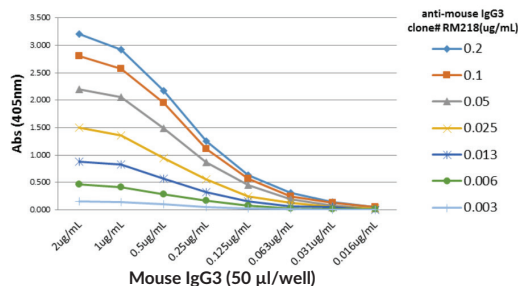
<b>Contents:</b>	This vial contains 50 µg of protein A-affinity purified monoclonal antibody.
<b>Synonym:</b>	Immunoglobulin G3
<b>Immunogen:</b>	Mouse IgG3
<b>Cross Reactivity:</b>	(+) IgG3; (-) Mouse IgG1, IgG2a, IgG2b, IgM, IgA, IgE; (-) Human, rat IgG
<b>Species Reactivity:</b>	(+) Mouse
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥1 year
<b>Storage Buffer:</b>	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
<b>Concentration:</b>	1 mg/ml
<b>Clone:</b>	RM218
<b>Host:</b>	Rabbit
<b>Isotype:</b>	IgG
<b>Application:</b>	ELISA; the recommended starting concentration is 0.005-0.2 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



#### IgG3 (mouse) Rabbit Monoclonal Antibody - Biotinylated (RM218)

ELISA of mouse immunoglobulins. IgG3 (mouse) Rabbit Monoclonal Antibody - Biotinylated (RM218) reacts to mouse IgG3k and IgG3λ and not to mouse IgG1, IgG2a, IgG2b, IgM, IgA, IgE, human IgG, or rat IgG. The plate was coated with 50 ng/well of different immunoglobulins. IgG3 (mouse) Rabbit Monoclonal Antibody - Biotinylated (RM218) was used as the primary antibody and an alkaline phosphatase-conjugated anti-rabbit IgG was used as the secondary antibody.



A Titer ELISA using IgG3 (mouse) Rabbit Monoclonal Antibody - Biotinylated (RM218). The plate was coated with different amounts of mouse IgG3. A serial dilution of IgG3 (mouse) Rabbit Monoclonal Antibody - Biotinylated (RM218) was used as the primary antibody and 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, 02/27/2024

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.<sup>1</sup> It is produced by B cells and later secreted by plasma cells and is the most abundant circulating antibody in human and mouse serum.<sup>1-3</sup> IgG consists of two heavy chains of approximately 50 kDa each and two light chains of approximately 25 kDa each.<sup>1</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> IgG exists as four isotypes in mice: IgG1, IgG2b, IgG3, and, in a strain-specific manner, IgG2a or IgG2c.<sup>5,6</sup> IgG3 production is driven by bacterial- or viral-associated antigens, including HIV-1 and *Staphylococcus* antigens, and occurs early in the immune response following IgM class-switching.<sup>2,7</sup> IgG3 binds to and neutralizes pathogens, as well as activates complement and opsonizes bacteria, leading to complement-dependent cytotoxicity (CDC) and antibody-dependent cell cytotoxicity (ADCC), respectively. Serum IgG3 levels are increased in patients with primary biliary cirrhosis, Sjögren's syndrome, systemic sclerosis, or systemic lupus erythematosus (SLE).<sup>8</sup> Cayman's IgG3 (mouse) Rabbit Monoclonal Antibody - Biotinylated (RM218) can be used for ELISA. This antibody recognizes the Fab region of IgG3 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-S52 (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. Damelang, T., Rogerson, S.J., Kent, S.J., et al. Role of IgG3 in infectious diseases. *Trends Immunol.* **40**(3), 197-211 (2019).
8. Zhang, H., Li, P., Wu, D., et al. Serum IgG subclasses in autoimmune diseases. *Medicine (Baltimore)* **94**(2), e387 (2015).

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