

CERTIFICATE of ANALYSIS



IgG Fc (mouse) Monoclonal Antibody (Clone RMG06)

Item No. 32106 • Batch No. 0825395 • Expires 25AUG2027

Overview

Formulation:	100 µg of protein G-affinity purified monoclonal antibody in 50% Glycerol/PBS with 1% BSA and 0.09% sodium azide
Concentration:	1 mg/mL *Buffer
Storage/stability:	-20°C/1yr
Applications:	ELISA
Cross reactivity:	(+)Mouse IgG1(+)Mouse IgG2a(+)Mouse IgG2b(+)Mouse IgG3 (weakly)(-)Human IgG(-)Mouse IgA(-)Mouse IgE(-)Mouse IgM(-)Rabbit IgG(-)Rat IgG
Species reactivity:	(+)Mouse

Reviewed and approved by: Justine Albertson

Approved date: 25AUG2026

A handwritten signature in black ink, appearing to read "Justine Albertson", written over a faint circular stamp.

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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