

CERTIFICATE of ANALYSIS



IgG Fab (mouse) Monoclonal Antibody (Clone RMG05)

Item No. 21780 • Batch No. 0625122 • Expires 14AUG2022

Overview

Formulation:	100 µg of protein G affinity-purified monoclonal antibody in PBS with 50% Glycerol, 1% BSA and 0.09% sodium azide
Concentration:	1 mg/mL
Storage/Stability:	-20°C/1yr
Applications:	ELISA
Cross reactivity:	(+)Mouse IgG1, (+)Mouse IgG2a, (+)Mouse IgG2b, (+)Mouse IgG3, (-)Human IgG, (-)Mouse IgA, (-)Mouse IgE, (-)Mouse IgM, (-)Rabbit IgG, (-)Rat IgG
Species reactivity:	(+)Mouse
Association rate constant:	0.00
Dissociation rate constant (1/s):	0.00
Dissociation equilibrium constant:	0.00
Maximum analyte response (RU):	0.00

Reviewed and approved by: Kyle Bale

Approved date: 14AUG2021

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