

PRODUCT INFORMATION

IgG Isotype Control Rabbit Monoclonal Antibody (Clone MOPC-21)
Item No. 42414

Overview and Properties

Contents:	This vial contains 200 µg of protein A-affinity purified monoclonal antibody.
Synonym:	Immunoglobulin G
Cross Reactivity:	Non-reactive
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 0.02% Proclin 300
Clone:	MOPC-21
Host:	Rabbit
Isotype:	IgG
Application:	Negative Control

Description

The IgG isotype control is a non-specific rabbit IgG produced from hybridoma clone MOPC-21 with rabbit Fc. Cayman's IgG Isotype Control Rabbit Monoclonal Antibody (Clone MOPC-21) can be used as a negative control IgG.

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, 12/11/2024

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