

PRODUCT INFORMATION

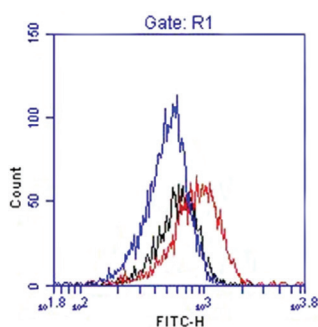


IgG1 Isotype Control FITC Monoclonal Antibody (Clone MOPC-21) Item No. 10343

Overview and Properties

Contents:	100 µg of protein G purified, fluorescein labeled monoclonal antibody
Synonyms:	IgG1κ Fluorescein Isotype Control, IgG1κ Isotype Control FITC
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	TBS, pH 7.4 with 50% glycerol
Clone:	MOPC-21
Host:	Mouse
Isotype:	IgG1
Applications:	Flow cytometry (FC) and immunofluorescence microscopy (IFM); the recommended working concentration for FC and IFM is 1 µg/test (100 tests provided). Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Image



Black: IgG1 FITC isotype control (1 µg)
Red: CD19 FITC (Clone CB19) Monoclonal Antibody (1 µg), stained 27% of lymphocytes; B-cells
Blue: Buffer control (no antibodies)

Fifty microliters of human whole blood stained with antibodies for 30 minutes on ice, lysed and analyzed on flow cytometer. 100,000 events collected and data gated on the lymphocyte cluster (R1).

Description

The IgG1 isotype control FITC is fluorescein conjugated non-specific mouse IgG1 produced from hybridoma clone MOPC-21. This conjugate should be used as a control antibody in direct fluorescence immunostaining experiments which utilize a mouse IgG1-fluorescein conjugate as the primary 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, 06/11/2020

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