

PRODUCT INFORMATION



Rabbit Anti-Mouse IgA Polyclonal FITC

Item No. 10196

Overview

Contents: This vial contains 500 µl of purified fluorescein-labeled IgG antibody.
Immunogen: Mouse IgA
Cross Reactivity: (+) Mouse IgA, (-) Mouse IgG

Properties

Form: Liquid
Storage: -20°C (as supplied)
Storage Buffer: 0.02 M potassium phosphate, 0.15 M sodium chloride, pH 7.2, containing 50% glycerol, 2 mg/ml BSA, and 0.02% sodium azide
Stability: ≥1 year
Host: Rabbit

Application

Flow Cytometry: Suggested starting dilution is 1:50-100

Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

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, 10/23/2023

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