

PRODUCT INFORMATION



Goat Anti-Mouse (IgG+IgM) FITC

Item No. 10006617

Overview and Properties

Contents:	This vial contains 500 µl whole IgG affinity-purified antibody conjugated to FITC.
Synonym:	Fluorescein-Conjugated Goat Anti-Mouse (IgG+IgM)
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2 with 50% glycerol, 1.13% BSA, and 0.04% sodium azide
Host:	Goat
Applications:	Immunofluorescence (IF); the recommended starting concentration is 1:100. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

This goat anti-mouse (IgG+IgM) FITC is used as the 'secondary antibody' for immunofluorescent staining on tissues or culture cells where the primary antibody is generated in mice. It is highly recommended that proper negative controls, such as omitting the primary antibody, be included in experiments when conducting immunofluorescent staining.

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, 05/01/2019

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