

PRODUCT INFORMATION



Goat Anti-Rabbit IgG FITC

Item No. 10006588

Overview and Properties

Contents:	This vial contains Fluorescein-labeled, purified IgG.
Synonym:	Fluorescein-Conjugated Goat Anti-Rabbit IgG
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	1 year
Storage Buffer:	0.015 M sodium phosphate, 0.375 M sodium chloride, pH 7.6, with 50% glycerol, 22.5 mg/ml BSA, and 0.075% sodium azide
Host:	Goat
Isotype:	IgG
Applications:	Immunofluorescent staining (IF); the recommended starting dilution is 1:100. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

This goat anti-rabbit IgG FITC is used as the "secondary antibody" for immunofluorescent staining on tissues or culture cells where the primary antibody was generated in rabbits. 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, 07/18/2023

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