

PRODUCT INFORMATION



Normal Rabbit IgG-FITC

Item No. 10010322

Overview and Properties

Contents:	This vial contains 100 µg of fluorescein-labeled, protein A-purified antibody.
Synonym:	Pre-immune IgG-FITC
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Applications:	Suitable for Flow cytometry (FC), immunofluorescence (IF), and Western blotting (WB), working concentration/dilution should be determined empirically.

Description

This vial contains purified rabbit pre-immune IgG conjugated to fluorescein (*via* isothiocyanate). The antibody exhibits low or no binding to most cells or protein lysates as demonstrated by FC, IF, and WB and thereby acts as an appropriate negative control for polyclonal IgG-FITC.

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, 04/15/2019

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