

Product Information



Goat Anti-Rabbit IgG FITC

Catalog No. 10006588 • Lot No. XXXXXX

| | |
|----------------------|---|
| Synonym: | Fluorescein conjugated Goat Anti-Rabbit IgG |
| Contents: | This vial contains XXX µg purified IgG-FITC in XXX µl PBS, pH 7.2, containing 50% glycerol |
| Host: | Goat |
| Storage: | ≥1 year at -20°C |
| Applications: | The recommended starting dilution for immunofluorescent staining is 1:100. Other applications have not been tested so optimal concentration should be determined empirically. |

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.

Related Products

Goat Anti-Rabbit IgG HRP - Cat. No. 10004301 • Goat Anti-Mouse IgG HRP - Cat. No. 10004302

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY; NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSD supervisor at your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy** located on our website and in our catalog.

Copyright Cayman Chemical Company, 08/17/2009

Cayman Chemical

Mailing address

1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone

(800) 364-9897
(734) 971-3335

Fax

(734) 971-3640

E-Mail

custserv@caymanchem.com

Web

www.caymanchem.com