

Product Information



CD5 Monoclonal FITC Antibody (Clone CRIS1)

Item No. 10232

- Contents:** This vial contains 300 µg of fluorescein-labeled, purified IgG_{2a} in 500 µl PBS, pH 7.2, containing 50% glycerol, 0.1% BSA, and 0.02% sodium azide.
- Synonyms:** Leu1, T1, Tp67
- Host:** Mouse, clone CRIS1 (clustered at the 1st International Workshop on Human Leukocyte Antigens, Paris, France (1982))
- Cross Reactivity:** (+) Human CD5; other species not tested
- Stability:** ≥1 year at -20°C
- Applications:** Flow cytometry (FC), immunofluorescence (IF), and immunohistochemistry (IHC) (frozen sections); the recommended starting dilution for FC is 1-3 µg/test and 1-10 µg/ml for IF and IHC (frozen sections). Not recommended for IHC (paraffin-embedded tissue) or western blot. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

This monoclonal antibody is labeled with fluorescein and intended for direct labeling of human CD5-expressing cells. CD5 is a type I transmembrane glycoprotein ($M_r = 67$ kDa), expressed on mature T-cells, the majority of thymocytes and B1a-subpopulation of B-cells.¹ CD5 is a marker of lymphocytes in autoimmune disorders (rheumatoid arthritis, systemic lupus erythematosus), B-cell chronic leukemia, T-cell acute lymphoblastic leukemia and small lymphocytic and mantle cell non-Hodgkin's lymphoma.

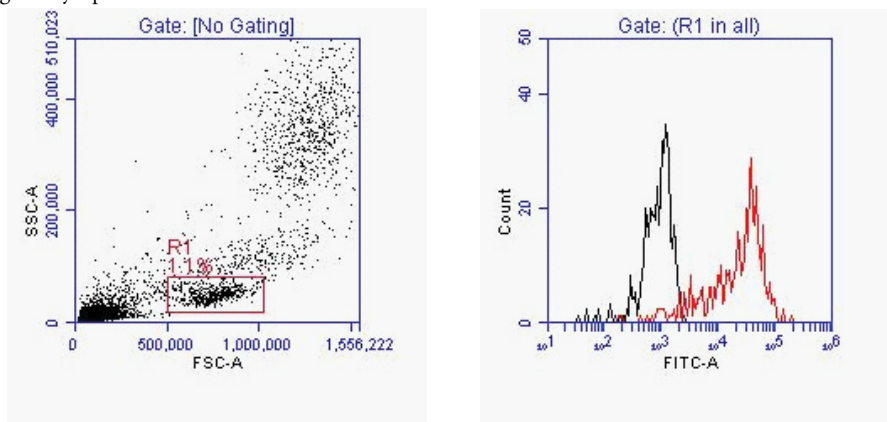


Figure 1: CD5 Monoclonal Antibody FITC (3 µg) was placed in a 1.5 ml tube followed by 50 µl human whole blood, mixed two seconds by vortex, and set at room temperature in the dark for 30 minutes. To this was added 1 ml RBC Lysis Buffer. The cell suspension was mixed again and incubated 10 minutes in the dark and read on an Accuri C6 cytometer. Fluorescein fluorescence (*Panel B*) from isotype control (left) and CD5 Monoclonal Antibody FITC (right) was generated from cell gate R1 (*Panel A*). CD5 Monoclonal Antibody FITC (3 µg) labeled 80% of the lymphocytes in the R1 gate. Enough antibody for 100 x 3 µg tests is provided per vial.

Reference

- Alberola-Ila, J., Places, L., Cantrell, D.A., *et al.* Intracellular events involved in CD5-induced human T cell activation and proliferation. *J. Immunol.* **148**, 1287-1293 (1992).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10232

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

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 Safety Data Sheet, which has been sent via email to 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, 01/13/2014

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