

Product Information



CD27 Monoclonal Antibody (Clone 203.6)

Item No. 10004837

- Contents:** This vial contains 100 µg of affinity-purified IgG₃ in 1 ml PBS, pH 7.2, containing 0.02% sodium azide.
- Host:** Mouse, clone 203.6
- Cross Reactivity:** (+) Human CD27; other species not tested
- Stability:** ≥1 year at 4°C
- Applications:** Flow cytometry, immunocytochemistry, and immunohistochemistry (frozen sections). Not recommended for western blot. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

CD27 is a type II homodimeric transmembrane glycoprotein ($M_r = 55$ kDa) and a member of the nerve growth factor/tumor necrosis receptor superfamily.^{1,2} The antigen was identified on medullar thymocytes, mature T lymphocytes (CD45RA⁺CD45RO⁻ T-cell subpopulation), memory B cells, plasma, and NK cells. This antibody is well suited for immunohistochemical detection of CD27 in B cell non-Hodgkin's lymphoma samples.

References

1. Prasad, K.V.S., Ao, Z., Yoon, Y., *et al.* CD27, a member of the tumor necrosis factor receptor family, induces apoptosis and binds to Siva, a proapoptotic protein. *Proc. Natl. Acad. Sci. USA* **94**, 6346-6351 (1997).
2. Camerini, D., Walz, G., Loenen, W.A.M., *et al.* The T cell activation antigen CD27 is a member of the nerve growth factor/tumor necrosis factor receptor gene family. *J. Immunol.* **147**, 3165-3169 (1991).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10004837

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

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, 11/25/2013