

# Product Information

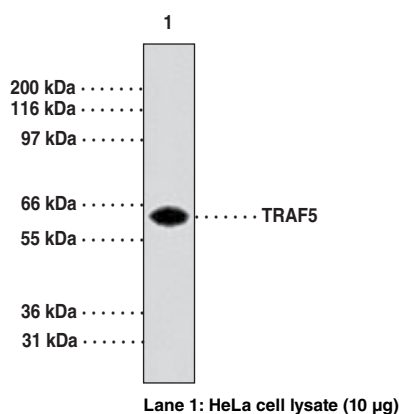


## TRAF5 Monoclonal Antibody (Clone 55A219)

Item No. 10873

- Contents:** This vial contains 100 µg of protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
- Antigen:** Fusion protein corresponding to amino acids 77-186 of human TRAF5
- Genbank Accession No.:** NP\_001029082
- Host:** Mouse
- Clone Designation:** 55A219
- Cross Reactivity:** (+) Human and mouse TRAF5
- Stability:** ≥6 months at -20°C
- Applications:** Western blot (WB); the recommended starting concentration for WB is 2 µg/ml. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Tumor necrosis factor (TNF)-induced signaling is mediated through association of TNF receptor (TNFR) with adaptor proteins, such as TNF receptor-associated factors (TRAFs). TRAFs form a family of cytoplasmic adapter proteins that mediate signal transduction from many members of the TNF-receptor superfamily (e.g., RANK, CD30, CD40, etc.) and the interleukin-1 receptor. The carboxy-terminal region of TRAFs is required for self-association and interaction with receptor cytoplasmic domains following ligand-induced oligomerization. Recent molecular cloning studies have led to identification of seven TRAFs (TRAF1-TRAF7).<sup>1-4</sup> Human TRAF5 is a 557-amino acid protein. TRAF5 is implicated in NF-κB and c-Jun NH(2)-terminal kinase/stress-activated protein kinase activation by members of the TNF receptor superfamily, including CD27, CD30, CD40, and lymphotoxin-β receptor. Targeted disruption of TRAF5 gene causes defects in CD40-CD27 mediated lymphocyte activation.<sup>5</sup>



### References

1. Cao, Z., Xiong, J., Takeuchi, M., *et al.* TRAF6 is a signal transducer for interleukin-1. *Nature* **383**, 443-446 (1996).
2. Rothe, M., Wong, S.C., Henzel, W.J., *et al.* A novel family of putative signal transducers associated with the cytoplasmic domain of the 75 kDa tumor necrosis factor. *Cell* **78**(4), 681-92 (1994).
3. Cheng, G., Cleary, A.M., Ye, Z.S., *et al.* Involvement of CRAF1, a relative of TRAF, in CD40 signaling. *Science* **267**(5203), 1494-8 (1995).
4. Nakano, H., Oshima, H., Chung, W., *et al.* TRAF5, an activator of NF-κB and putative signal transducer for the lymphotoxin-β receptor. *J. Biol. Chem.* **271**(25), 14661-4 (1996).
5. Nakano, H., Sakon, S., Koseki, H., *et al.* Targeted disruption of *traf5* gene causes defects in CD40- and CD27-mediated lymphocyte activation. *Proc. Natl. Acad. Sci. USA* **96**(17), 9803-8 (1999).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10873](http://www.caymanchem.com/catalog/10873)

**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 herof. 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, 05/02/2013

### 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](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)