

PRODUCT INFORMATION



TRAF5 Monoclonal Antibody (Clone 55A219)

Item No. 10873

Overview and Properties

Contents: This vial contains 100 µg of protein G-purified IgG in 200 µl PBS, with 0.05% BSA and 0.05% sodium azide.

Immunogen: Fusion protein corresponding to amino acids 77-186 of human TRAF5

Species Reactivity: (+) Human and mouse TRAF5

Storage: -20°C (as supplied)

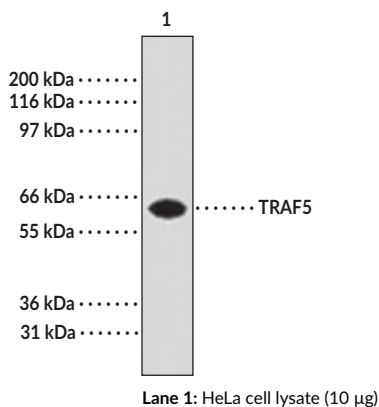
Stability: ≥6 months

Clone: 55A219

Host: Mouse

Applications: Western blot (WB); the recommended starting concentration is 2 µg/ml. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



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, 11/13/2024

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

PRODUCT INFORMATION



Description

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).¹⁻⁵ Human TRAF5 is a 557-amino acid protein. TRAF5 is implicated in NF- κ B 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.⁵

References

1. 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)**, 9803-9808 (1999).
2. Cao, Z., Xiong, J., Takeuchi, M., *et al.* TRAF6 is a signal transducer for interleukin-1. *Nature* **383(6599)**, 443-446 (1996).
3. 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).
4. 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).
5. Nakano, H., Oshima, H., Chung, W., *et al.* TRAF5, an activator of NF- κ B and putative signal transducer for the lymphotoxin- β receptor. *J. Biol. Chem.* **271(25)**, 14661-14664 (2011).

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