

Product Information



TRAIL Polyclonal Antibody

Item No. 160750

Synonym:	Apo-2L
Contents:	This vial contains 200 µg IgG in 500 µl PBS, pH 7.4 containing 0.02% sodium azide.
Host:	Rabbit
Antigen:	Human TRAIL amino acids 261-277 (CTNEHLIDMDHEA-SFFGA) ^{1,2} Homology of the peptide antigen with TRAIL from other species: Human TNEHLIDMDHEASFFGAF Mouse TNEHLmD i DqEASFFGAF
Stability:	≥1 year at 4°C
Application:	This polyclonal antibody can be used for detection of TRAIL by Western blot at a dilution of 1:500. A whole cell lysate from HeLa cells can be used as a positive control and a 32 kDa band should be detected.

Apoptosis, or programmed cell death, occurs during normal cellular differentiation and development of multicellular organisms. Apoptosis is induced by certain cytokines including TNF and Fas ligand in the TNF family through their death domain containing receptors, TNFR-1 and Fas. A novel member in the TNF family was recently identified and designated TRAIL (TNF-related apoptosis-inducing ligand) and Apo-2L (Apo-2 ligand).^{1,2} TRAIL is a type II membrane protein and is expressed in a variety of human tissues.¹ Two novel death domain containing receptors DR4 and DR5 have been identified as the receptor for TRAIL.³⁻⁶ Like TNF and Fas ligand, TRAIL induces apoptosis and NF-κB activation in many tissues and cells.

References

1. Wiley, S.R., Schooley, K., Smolak, P.J., *et al.* Identification and characterization of a new member of the TNF family that induces apoptosis. *Immunity* **3**, 673-682 (1995).
2. Pitti, R.M., Marsters, S.A., Ruppert, S., *et al.* Induction of apoptosis by Apo-2 ligand, a new member of the tumor necrosis factor cytokine family. *J. Biol. Chem.* **271**, 12687-12690 (1996).
3. Pan, G., O'Rourke, K., Chinnaiyan, A.M., *et al.* The receptor for the cytotoxic ligand TRAIL. *Science* **276**, 111-113 (1997).
4. Schneider, P., Thome, M., Burns, K., *et al.* TRAIL receptors 1 (DR4) and 2 (DR5) signal FADD-dependent apoptosis and activate NF-κB. *Immunity* **7**, 831-836 (1997).
5. Pan, G., Ni, J., Wei, Y.-F., *et al.* An antagonist decoy receptor and a death domain-containing receptor for TRAIL. *Science* **277**, 815-817 (1997).
6. Sheridan, J.P., Marsters, S.A., Pitti, R.M., *et al.* Control of TRAIL-induced apoptosis by a family of signaling and decoy receptors. *Science* **277**, 818-821 (1997).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/160750

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 *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, 12/09/2011

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