

# Product Information



## DcR2 Polyclonal Antibody

Catalog No. 160755

<b>Synonyms:</b>	TRAIL-R4; TRUNDD
<b>Contents:</b>	This vial contains 100 µg of affinity-purified IgG in 200 µl PBS, containing 0.02% azide.
<b>Host:</b>	Rabbit
<b>Antigen:</b>	The anti-DcR2 polyclonal antibody was raised against a peptide corresponding to amino acids 249 to 263 of human DcR2 precursor.
<b>Cross Reactivity:</b>	(+) human DcR2; other species not tested.
<b>Stability:</b>	≥1 one year at 4°C
<b>Application:</b>	This polyclonal antibody can be used for detection of DcR2 by western blot at a dilution of 1:1,000. A whole cell lysate from HeLa cells can be used as a positive control.

Apoptosis is induced by certain cytokines including TNF and Fas ligand in the TNF family through their death domain containing receptors. TRAIL/Apo2L is a new member of the TNF family and induces apoptosis of a variety of tumor cell lines. DR4 and DR5 are the recently identified functional receptors for TRAIL, and DcR1/TRID is a decoy receptor.<sup>1-3</sup> Another member of the TRAIL receptor family was more recently identified and designated DcR2, TRAIL-R4, or TRUNDD.<sup>4-6</sup> DcR2 has an extracellular TRAIL-binding domain but lacks an intracellular death domain and does not induce apoptosis. Like DR4 and DR5, DcR2 transcript is widely expressed in normal human tissues. Overexpression of DcR2 attenuates TRAIL-induced apoptosis.

### References

1. Pan, G., O'Rourke, K., Chinnaiyan, A.M., *et al.* The receptor for the cytotoxic ligand TRAIL. *Science* **276**, 111-113 (1997).
2. 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).
3. 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).
4. Marsters, S.A., Sheridan, J.P., Pitti, R.M., *et al.* A novel receptor for Apo2L/TRAIL contains a truncated death domain. *Curr. Biol.* **7**, 1003-1006 (1997).
5. Degli-Esposti, M.A., Dougall, W.C., Smolak, P.J., *et al.* The novel receptor TRAIL-R4 induces NF- $\kappa$ B and protects against TRAIL-mediated apoptosis, yet retains an incomplete death domain. *Immunity* **7**, 813-820 (1997).
6. Pan, G., Ni, J., Yu, G.-L., *et al.* TRUNDD, a new member of the TRAIL receptor family that antagonizes TRAIL signalling. *FEBS Lett.* **424**, 41-45 (1998).

### Related Product

TRAIL Polyclonal Antibody - Cat. No. 160750

### 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 NOT INTENDED OR APPROVED FOR HUMAN OR VETERINARY USE. USE OF THIS PRODUCT FOR HUMAN OR ANIMAL TESTING IS EXTREMELY HAZARDOUS AND MAY RESULT IN DISEASE, SEVERE INJURY, OR DEATH.**

#### 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 under separate cover to the MSDS supervisor at 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, 09/25/2007