

PRODUCT INFORMATION



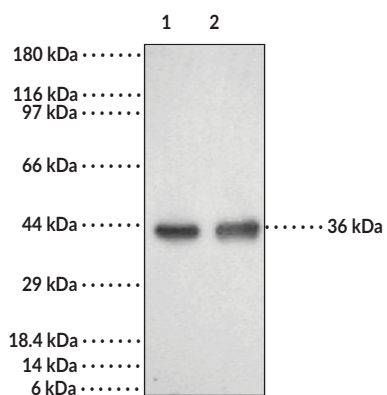
DcR2 Polyclonal Antibody

Item No. 160755

Overview and Properties

Contents:	This vial contains 200 µl of peptide affinity purified polyclonal antibody
Synonyms:	TRAIL-R4; TRUNDD
Immunogen:	Synthetic peptide from an internal region of human DcR2
Species Reactivity:	(+) Human, other species not tested
Uniprot No.:	Q9UBN6
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2 with 0.02% sodium azide
Host:	Rabbit
Applications:	Western Blot (WB); the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: HeLa Cell Lysate (25 µg)
Lane 2: HeLa Cell Lysate (50 µg)

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, 10/24/2023

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

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.¹⁻³ Another member of the TRAIL receptor family was more recently identified and designated DcR2, TRAIL-R4, or TRUNDD.⁴⁻⁶ 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(5309)**, 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(5327)**, 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(5327)**, 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(12)**, 1003-1006 (1997).
5. Degli-Esposti, M.A., Dougall, W.C., Smolak, P.J., *et al.* The novel receptor TRAIL-R4 induces NF- κ B and protects against TRAIL-mediated apoptosis, yet retains an incomplete death domain. *Immunity* **7(6)**, 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(1-2)**, 41-45 (1998).

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