

TRAIL-R4 polyclonal antibody

Decoy receptors (DcRs) are signaling-deficient receptors for certain members of the TNF family of cytokines, including TRAIL/Apo2L and FasL. Functional TRAIL receptors such as DR4 and DR5 transmit apoptotic signals via their intracellular death domains. In contrast, three identified TRAIL decoy receptors (DcR1, DcR2, and osteoprotegerin/OPG) completely or partially lack the signaling death domain, and/or are secreted from cells and unable to transmit intracellular signals. Thus, expression of DcRs promotes survival by sequestration of apoptotic ligands. A fourth identified decoy receptor, DcR3, binds both FasL and the vascular endothelial cell growth inhibitor (VEGI), inhibiting Fas and DR3 death signals.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[View Online »](#)

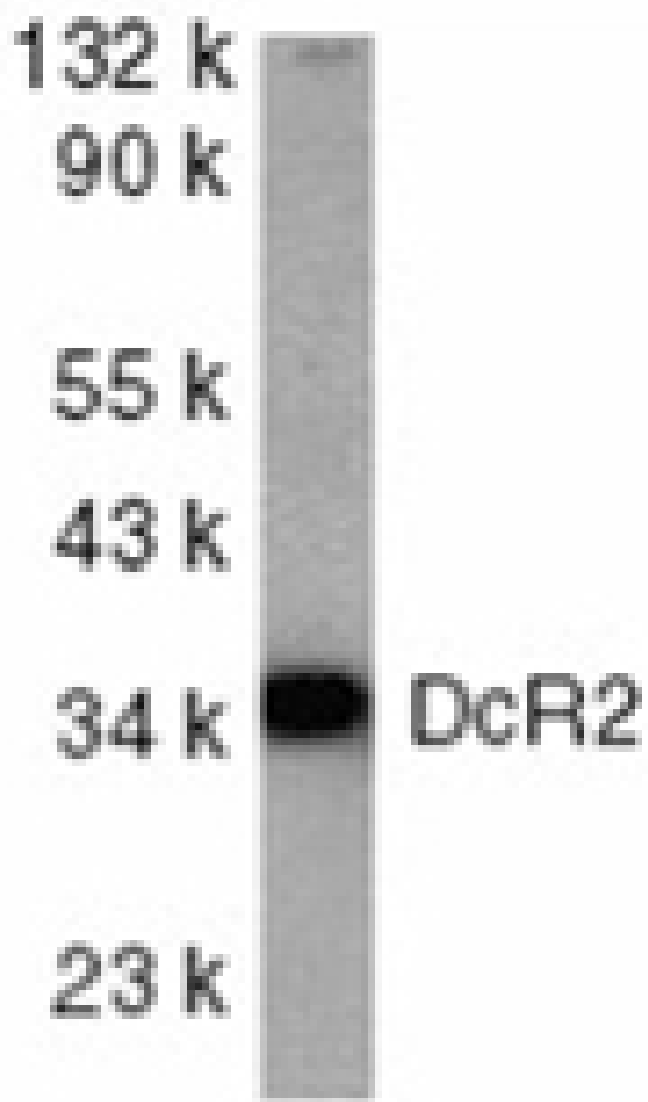
Ordering Information

[Order Online »](#)

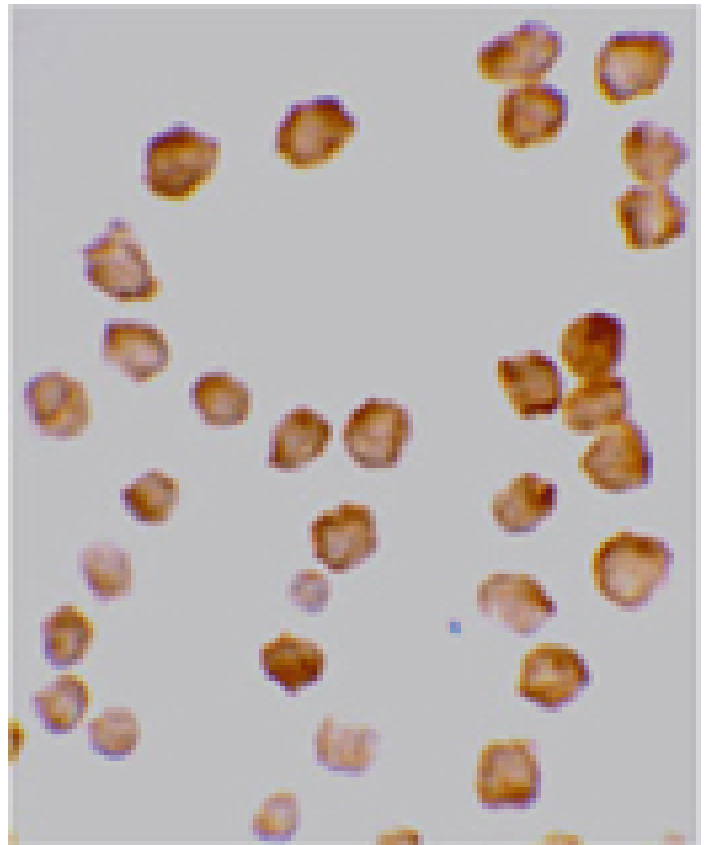
ADI-AAP-371-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of DcR2 in HeLa whole cell lysate with DcR2 antibody at 1:1000 dilution.



Immunocytochemistry analysis of HeLa cells using DcR2 antibody at 10 µg/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TRAIL receptor 4, DcR2, Death Receptor 2, TRUNDD, CD264, TNFRSF 10D, TNFRSF 10D, TNF-related apoptosis-inducing ligand receptor 4, Tumor necrosis factor receptor superfamily member 10D
Application	ICC, IHC, WB
Application Notes	Detects a band of ~36kDa by Western blot.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
GenBank ID	AF029761
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human TRAIL-R4.
Purity Detail	Peptide affinity purified.
Recommendation Dilutions/Conditions	Immunocytochemistry (10µg/ml)Western Blot (1:1,000-1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q9UBN6

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.