

TRAIL-R2 (human) monoclonal antibody (DJR2-2)

DR5 is a 55kDa member 10B of the TNF receptor superfamily (TNFRSF10B), also known as TRAIL-R2, TRICK2, KILLER, and CD262. It binds the cytotoxic ligand TRAIL and induces apoptosis. The DR5 receptor is broadly expressed on a variety of normal tissues and many tumors. DR5 expression has been reported to be upregulated in human cells by interferon- α , 2-methoxyestradiol, and paclitaxel, and downregulated by adenoviral E3 proteins.

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

Ordering Information

[Order Online »](#)

ALX-804-913-0100	100 μ g
------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TRAIL receptor 2, DR5, Death receptor 5, KILLER, CD262, Trick2, Ly98, TNFRSF10B, TNF-related apoptosis-inducing ligand receptor 2, Tumor necrosis factor receptor superfamily member 10B

Application Flow Cytometry, IHC (PS)

Clone DJR2-2

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

Gene/Protein Identifier 8795 (Entrez GeneID)

Host Mouse

Immunogen Recombinant human TRAIL-R2 extracellular domain IgG1 Fc fusion protein.

Isotype IgG1

Purity Detail Affinity purified.

Species Reactivity Human

UniProt ID O14763

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-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

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