

# TRAIL-R1 (human) monoclonal antibody (DJR1)

DR4 is 56kDa member 10A of the TNFR superfamily (TNFRSF10A), also known as TRAIL-R1, Apo-2, and CD261. It is expressed at low levels by activated T cells and some tumors. After TRAIL engagement, DR4 (TRAIL-R1), through activation of NF- $\kappa$ B, induces apoptosis in the TRAIL ligated cell.

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

Citations: 1

[View Online »](#)

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 1, Apo-2, CD261, DR4, Death receptor 4, TNFRSF 10A, TNF-related apoptosis-inducing ligand receptor 1, Tumor necrosis factor receptor superfamily member 10A
Application	Flow Cytometry, WB
Clone	DJR1
Formulation	Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.
GenBank ID	U90875
Gene/Protein Identifier	8798 (Entrez GeneID)
Host	Mouse
Immunogen	Recombinant human TRAIL-R1 extracellular domain IgG1 Fc fusion protein.
Isotype	IgG1
Purity Detail	Affinity purified.
Species Reactivity	Human
UniProt ID	O00220

**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)

Last modified: May 29, 2024