

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](#).

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