

TRAIL-R3 (human) monoclonal antibody (DJR3)

DcR1 is a 65kDa GPI-anchored member of the TNFR superfamily (TNFRSF10C) also known as TRAIL-R3, TRID, CD263, and LIT. Although not expressed by most tumors, DcR1 is expressed at low levels on normal tissues, acting as a TRAIL decoy receptor, thus inhibiting apoptosis. DcR1 is termed a decoy receptor, as it does not have a functional death domain to induce apoptosis following TRAIL ligation.

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

Ordering Information

[Order Online »](#)

ALX-804-908-0100	100µg
------------------	-------

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 3, DcR1, Death Receptor 1, TRID, CD263, TNFRSF 10C, TNF-related apoptosis-inducing ligand receptor 3, Tumor necrosis factor receptor superfamily member 10C
Application	Flow Cytometry
Clone	DJR3
Formulation	Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.
Gene/Protein Identifier	8794 (Entrez GeneID)
Host	Mouse
Immunogen	Recombinant human TRAIL-R3 extracellular domain IgG1 Fc fusion protein.
Isotype	IgG1
Purity Detail	Affinity purified.
Species Reactivity	Human
UniProt ID	O14798
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee