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 GenelD)

Host Mouse

Immunogen Recombinant human TRAIL-R3 extracellular domain IgG1

Fc fusion protein.

lsotype lgG1

Purity Detail Affinity purified.

Species Reactivity Human

UniProt ID O14798

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

