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-κB, induces apoptosis in the TRAIL ligated cell.

This antibody is covered by our Worry-Free Guarantee.

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-804-912-0100

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 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 GenelD)

Host Mouse

Immunogen Recombinant human TRAIL-R1 extracellular domain IgG1

Fc fusion protein.

lsotype lgG1

Purity Detail Affinity purified.

Species Reactivity Human

UniProt ID 000220

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

ENZO LIFE SCIENCES, INC.
Phone: 800.942.0430 infousa@enzolifesciences.cor

Last modified: May 29, 2024

European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
infode@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com