

# TRAIL-R4 (human) monoclonal antibody (01)

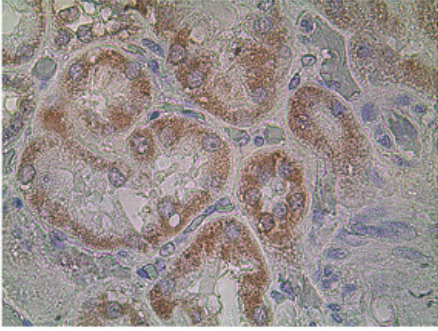
TRAIL-R4 is a type 1 transmembrane receptor that exhibits a truncated cytoplasmic domain.

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

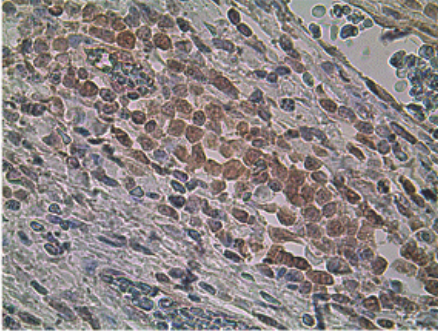
Manuals, SDS & CofA

[View Online »](#)

**A**



**B**



Immunohistochemistry staining of formalin fixed, paraffin-embedded human prostate tumor tissue using TRAIL-R2 (human), mAb (DJR2-2) at 1:50 dilution showing the positive staining in prostate carcinoma cells (63x).

## Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	TRAIL receptor 4, DcR2, Death Receptor 2, TRUNDD, CD264, TNFRSF 10D, TNF-related apoptosis-inducing ligand receptor 4, Tumor necrosis factor receptor superfamily member 10D
Application	Flow Cytometry, IHC (PS)
Clone	01
Formulation	Liquid. In PBS, pH 7.4. Contains no preservatives.
Host	Mouse
Immunogen	Recombinant human TRAIL-R4 (aa 1-210) IgG1 Fc fusion protein.
Isotype	IgG1
Recommendation Dilutions/Conditions	Flow Cytometry (1:10-1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q9UBN6

### 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