

# 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](#).

## Ordering Information

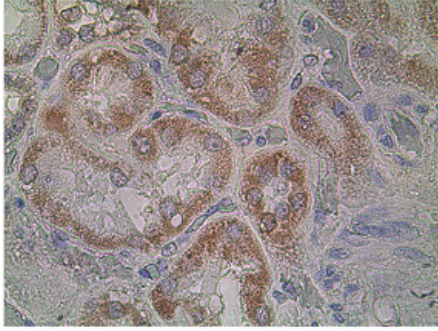
[Order Online »](#)

ALX-804-915-0100	100µg
------------------	-------

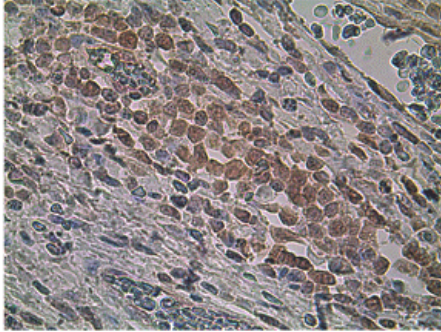
## Manuals, SDS & CofA

[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

<b>Alternative Name</b>	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
<b>Application</b>	Flow Cytometry, IHC (PS)
<b>Clone</b>	01
<b>Formulation</b>	Liquid. In PBS, pH 7.4. Contains no preservatives.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human TRAIL-R4 (aa 1-210) IgG1 Fc fusion protein.
<b>Isotype</b>	IgG1
<b>Recommendation Dilutions/Conditions</b>	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.
<b>Species Reactivity</b>	Human
<b>UniProt ID</b>	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-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
uk@enzolifesciences.com](mailto:info-uk@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**

For local distributors and detailed product information visit us online: [enzo.com](http://enzo.com) All product names, logos, and brands are the property of their respective owners.