

# TRAIL-R2 monoclonal antibody (DR5/3381)

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

## Ordering Information

[Order Online »](#)

|                 |       |
|-----------------|-------|
| ENZ-ABS924-0100 | 100µg |
|-----------------|-------|

## Manuals, SDS & CofA

[View Online »](#)

# Handling & Storage

|                   |                                                                                                       |
|-------------------|-------------------------------------------------------------------------------------------------------|
| Handling          | Store the DR5 antibody at 2-8°C (with azide) or aliquot and store at -20°C or colder (without azide). |
| Long Term Storage | -20°C                                                                                                 |
| Shipping          | Blue Ice                                                                                              |

## Regulatory Status

RUO - Research Use Only

# Product Details

|                                     |                                                                                                                                                                                         |
|-------------------------------------|-----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Alternative Name                    | TNF-related apoptosis-inducing ligand receptor 2, TRAIL-R2, DR5, Death receptor 5, Tumor necrosis factor receptor superfamily member 10B, TNFRSF10B, CD262                              |
| Application                         | IHC (PS)                                                                                                                                                                                |
| Clone                               | DR5/3381                                                                                                                                                                                |
| Formulation                         | Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced), 0.05% sodium azide.                                                                                                  |
| Host                                | Mouse                                                                                                                                                                                   |
| Immunogen                           | Recombinant TRAIL-R2 (aa 266-393).                                                                                                                                                      |
| Isotype                             | IgG1κ                                                                                                                                                                                   |
| Purity Detail                       | Protein G affinity chromatography purified.                                                                                                                                             |
| Recommendation Dilutions/Conditions | Immunohistochemistry (FFPE) (1-2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| Species Reactivity                  | Human                                                                                                                                                                                   |
| UniProt ID                          | O14753                                                                                                                                                                                  |
| Worry-free Guarantee                | This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .                                                                                                                  |



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)