

# TL1A recombinant monoclonal antibody (TLRM1-2R)

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

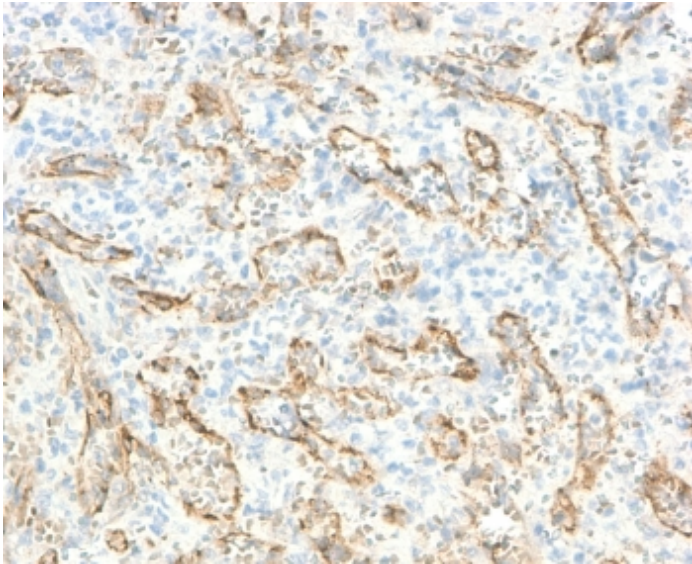
## Ordering Information

[Order Online »](#)

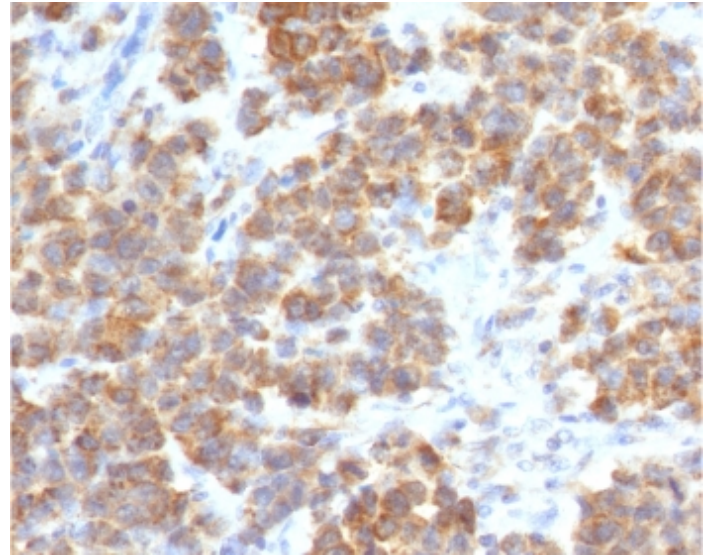
ENZ-ABS923-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



IHC staining of FFPE human spleen with recombinant TNFSF15 antibody (clone TLRM1-2R). Required HIER: boil sections in 10mM Tris with 1mM EDTA, pH9, for 10-20 min followed by cooling at RT for 20 min.



IHC staining of FFPE human parathyroid mass with recombinant VEGI antibody (clone TLRM1-2R). Required HIER: boil sections in 10mM Tris with 1mM EDTA, pH9, for 10-20 min followed by cooling at RT for 20 min.

# Handling & Storage

Handling	Store the recombinant VEGl 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	VEGl, Vascular endothelial cell growth inhibitor, Tumor necrosis factor ligand superfamily member 15, TNFSF15
Application	IHC (PS)
Clone	TLRM1-2R
Formulation	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.
Host	Rabbit
Immunogen	Full length human recombinant VEGl.
Isotype	IgGk
Purity Detail	Protein A 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	O95150
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)