

# VEGFA recombinant monoclonal antibody (RM391)

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

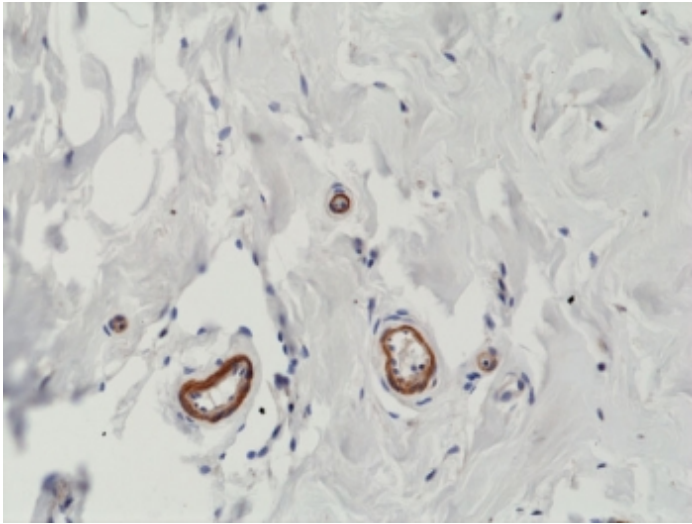
## Ordering Information

[Order Online »](#)

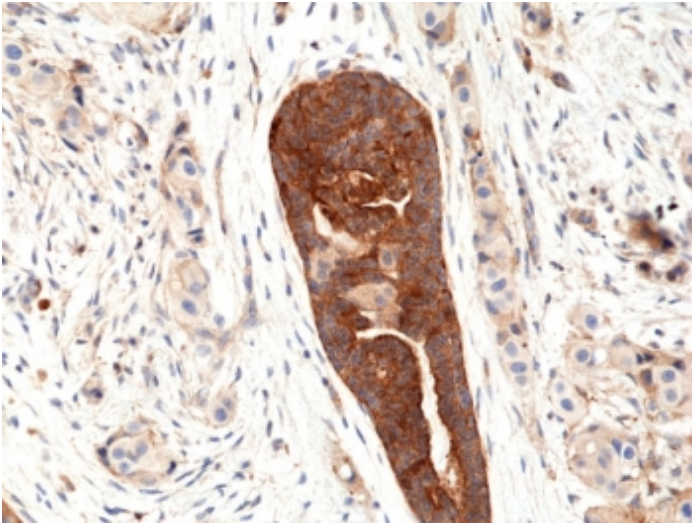
ENZ-ABS916-0100	100µl
-----------------	-------

## Manuals, SDS & CofA

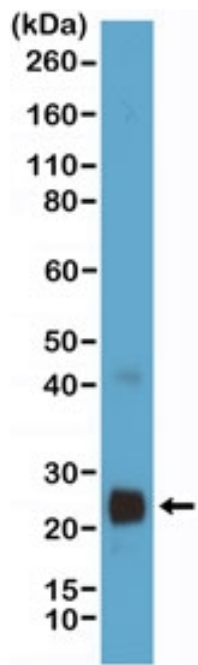
[View Online »](#)



IHC staining of FFPE human breast tissue with recombinant VEGFA antibody at 1:200.



IHC staining of FFPE human breast cancer tissue with recombinant VEGFA antibody at 1:100.



Western blot testing of VEGF protein-expressing human HEK293 cell lysate with recombinant VEGFA antibody.

## Handling & Storage

Long Term Storage                    -20°C

Shipping                                Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Vascular endothelial growth factor A, Vascular permeability factor, VPF
Application	IHC (PS), WB
Clone	RM391
Formulation	Liquid. In PBS containing 50% glycerol, 1% BSA and 0.09% sodium azide.
Host	Rabbit
Immunogen	Peptide corresponding to human VEGFA .
Isotype	IgG
Purity Detail	Protein A purified from animal origin-free supernatant.
Recommendation Dilutions/Conditions	Immunohistochemistry (FFPE) (1:100-1:200)Western blot (1:100-1:200)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P15692
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>



ENZO LIFE SCIENCES,  
INC.  
Phone: +1 949 949 1129  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

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

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

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)