

VEGFA recombinant monoclonal antibody (RM391)

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

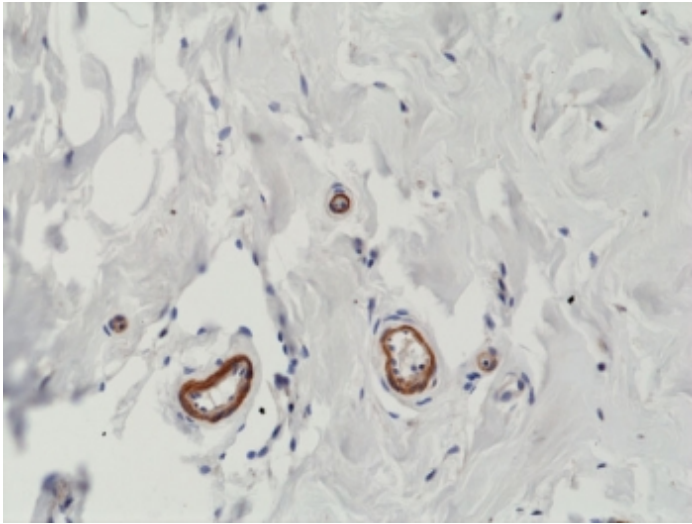
Ordering Information

[Order Online »](#)

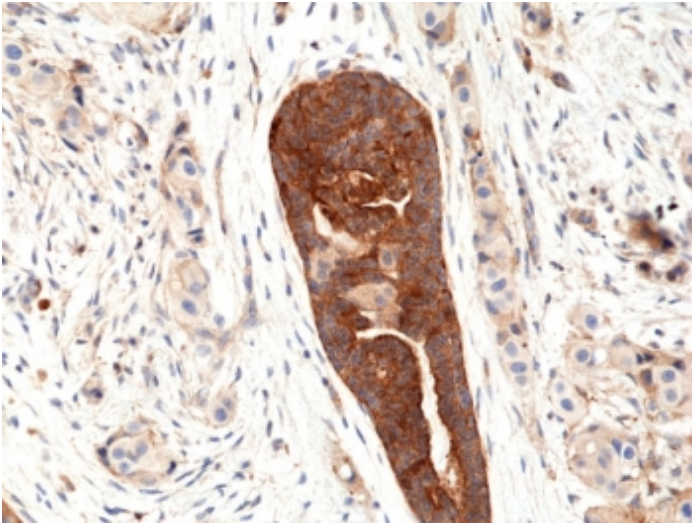
ENZ-ABS916-0100	100µl
-----------------	-------

Manuals, SDS & CofA

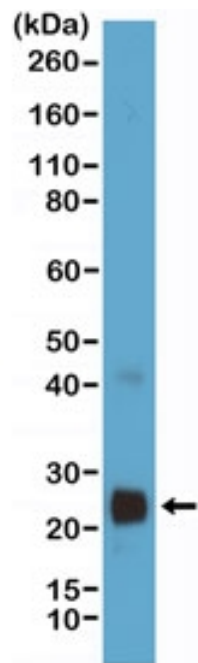
[View Online »](#)



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 [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)