

# VEGF monoclonal antibody (6B7)

VEGF (Vascular endothelial growth factor) is a potent polypeptide regulator of blood vessel function. VEGF promotes an array of responses in the endothelium, including hyperpermeability, endothelial cell growth, angiogenesis, and enhanced glucose transport. Several alternate splice isoforms of VEGF are known. Two high-affinity receptors for VEGF, VEGFR1 and VEGFR2, have been identified and characterized.

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

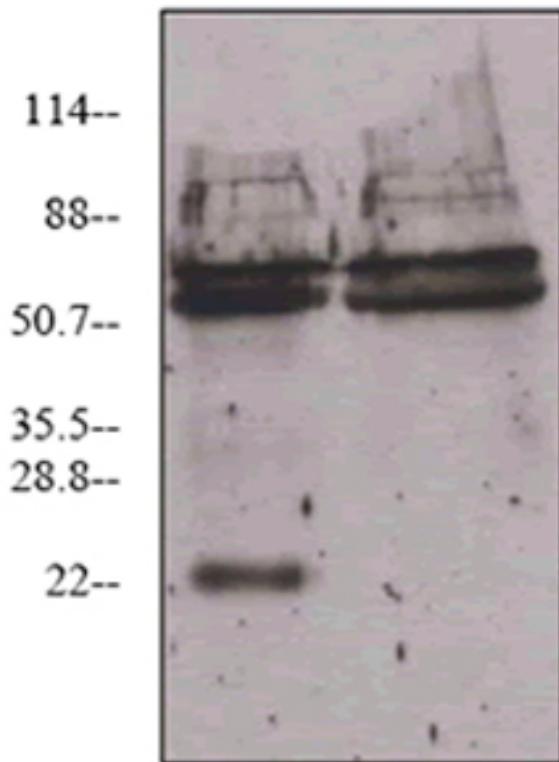
## Ordering Information

[Order Online »](#)

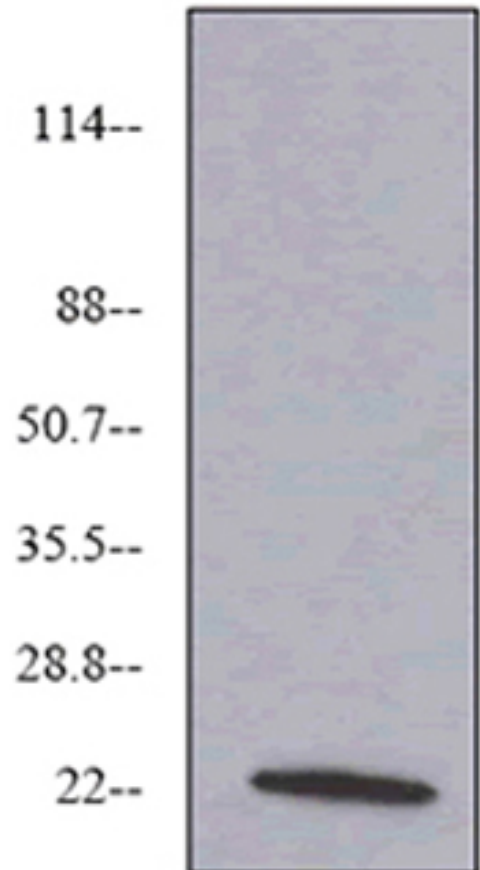
ADI-905-164-100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunoprecipitation analysis of VEGF. Lane 1- 1 $\mu$ g VEGF immunoprecipitated with 2 $\mu$ g antibody; Lane 2- HUVEC lysate immunoprecipitated with 2 $\mu$ g antibody.



Western blot analysis of VEGF probed with VEGF, mAb (6B7) (0.1 $\mu$ g).

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Vascular endothelial growth factor

**Application** IP, WB

**Application Notes** Detects a band of ~22kDa by Western blot.

**Clone** 6B7

**Formulation** Liquid. In PBS containing 0.09% sodium azide.

**Host** Mouse

**Immunogen** Human Vascular Endothelial Growth Factor 165.

**Isotype** IgG1

**Purity Detail** Protein G affinity purified.

**Source** Purified from hybridoma tissue culture supernatant.

**Species Reactivity** Human, Mouse

**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)