

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](#).

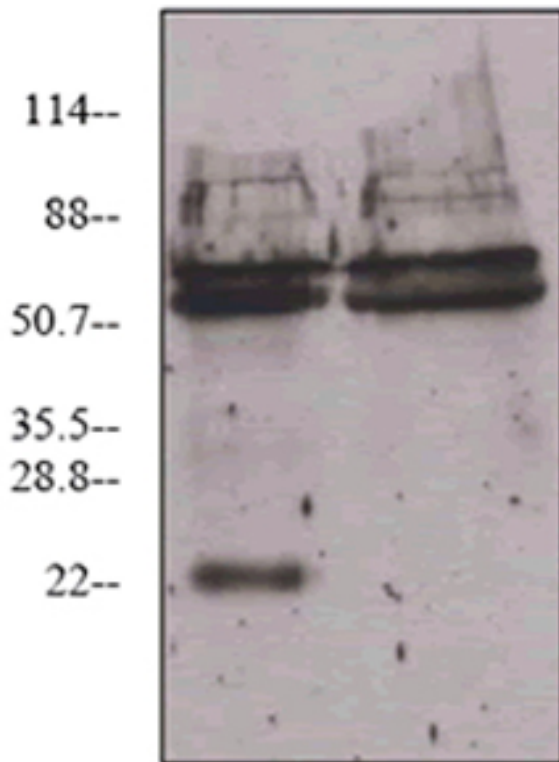
Ordering Information

[Order Online »](#)

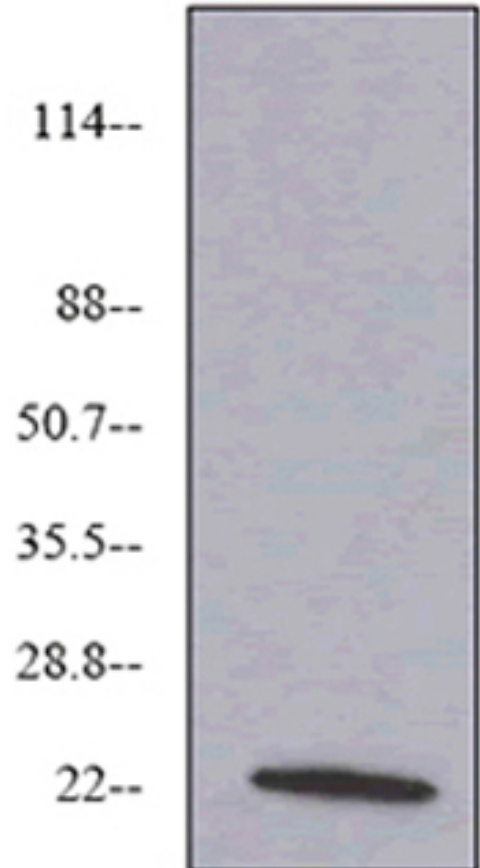
ADI-905-164-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



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



Western blot analysis of VEGF probed with VEGF, mAb (6B7) (0.1 μ 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 .