

VEGFA monoclonal antibody (R012)

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

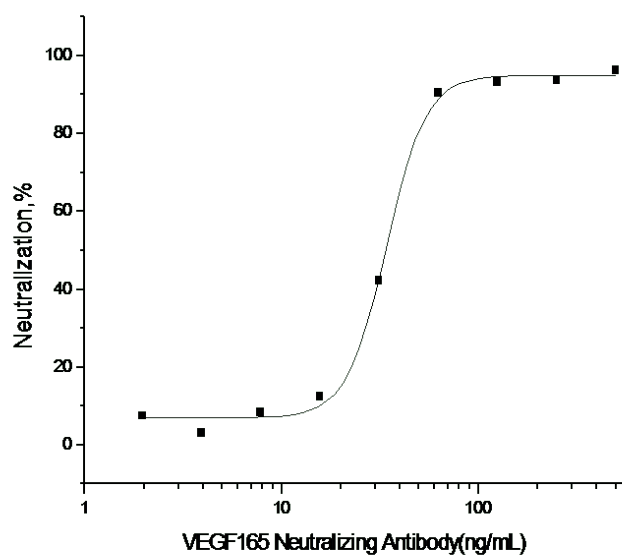
Ordering Information

[Order Online »](#)

ENZ-ABS971-0200	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Measured by its ability to neutralize VEGF165 induced proliferation of human umbilical vein endothelial cells (HUVEC). The Neutralization Titer (IC50) is typically 15-75ng/mL in the presence of 10 ng/mL Recombinant Human VEGF165.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name VEGF165, Vascular endothelial growth factor A, Vascular permeability factor, VPF

Application Neutralization

Clone R012

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant human / Cynomolgus VEGF165/VEGFA protein.

Isotype IgG

Purity Detail Protein A purified.

Species Reactivity Cynomolgus, Human

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

.

Last modified: September 17, 2024



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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

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