

VEGF Recombinant monoclonal antibody (VG76e) (Rabbit IgGκ)

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

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

| | |
|---------------------------|---|
| Use/Stability | Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C. |
| Handling | Avoid freeze/thaw cycles. |
| Short Term Storage | +4°C |
| Long Term Storage | -20°C |
| Shipping | Blue Ice |

Regulatory Status RUO - Research Use Only

Product Details

| | |
|---------------------------------------|---|
| Alternative Name | Vascular endothelial growth factor, VEGF-A, Vascular permeability factor, VPF |
| Application | ELISA, IHC, WB |
| Clone | VG76e |
| Formulation | Liquid. In PBS containing 0.02% Proclin 300. |
| Host | Rabbit |
| Immunogen | Recombinant human VEGF189 expressed in E. coli. |
| Isotype | IgGk |
| Purity Detail | Protein A affinity purified. |
| Species Reactivity | Human, Mouse, Primate |
| Technical Info / Product Notes | This chimeric rabbit antibody was made using the variable domain sequences of the original Mouse IgG1 format, for improved compatibility with existing reagents, assays and techniques. |
| UniProt ID | P15692 |

Last modified: May 29, 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-
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)