

# VEGF Recombinant monoclonal antibody (VG76e) (Mouse IgG1κ)

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

## Ordering Information

[Order Online »](#)

|                 |       |
|-----------------|-------|
| ENZ-ABS659-0200 | 200µg |
|-----------------|-------|

## 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               | Mouse   |
| Immunogen          | Recombinant human VEGF189 expressed in E. coli.                               |
| Isotype            | IgG1k   |
| Purity Detail      | Protein A affinity purified.  |
| Species Reactivity | Human, Mouse, Primate   |
| UniProt ID         | P15692  |

### Worry-free Guarantee

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

.