

Data Sheet

Anti-VEGF Neutralizing Antibody

Recombinant, Human, Monoclonal

Catalog #: 79478-2

Lot #: 180220

Conc.: 1.61 mg/ml

Host Species | Isotype:

Human | IgG1

Formulated in: 8mM phosphate, 110mM NaCl, 2.2mM KCl, pH7.4, and 20% glycerol

Stability: Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles. Store diluted antibody with carrier protein (BSA 0.1 – 0.5%) at 4°C, stable for 2 weeks.

References:

1. Rosenfeld, P. et al., *Healio*. 2005; 36(4): 331-335.
2. Avery, R., et al., *Ophthalmology*. 2006; 113(3): 363-372.

Description: Anti-VEGF Neutralizing Recombinant Human antibody. This antibody has not been tested for cross-reactivity with other species.

Purification: Protein A affinity purification from HEK293 supernatants.

Application: VEGF is a cell surface protein which signals for the growth of blood vessels. Anti-VEGF monoclonal antibodies are used to block tumor angiogenesis.

Specific Activity

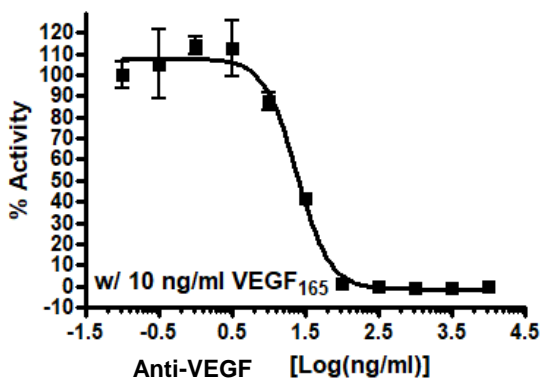


Fig.1 The Anti-VEGF Antibody was evaluated using KDR-NFAT-Luc (HEK293) Cell Line (BPS# 79387). VEGF165 (10 ng/ml at final concentration) was incubated with various concentrations of Anti-VEGF Antibody in the assay medium (DMEM + 1% BSA) for 45 min at room temperature and added to the cells. After 4 hour incubation at 37°C under 5% CO₂, ONE-Step Luciferase Assay System (BPS# 60690) was used to measure the luciferase activity.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com