

Data Sheet

Anti-Beta NGF (Beta Nerve Growth Factor) Neutralizing Antibody

Recombinant, Human, Monoclonal

Catalog #: 100436

Lot #: 190625 **Conc.:** 2.29 mg/ml

Host Species | Isotype:

Human | IgG1

Formulated in: 8 mM phosphate, 110 mM NaCl, 2.2 mM KCl, pH 7.4, and 20% glycerol

Stability: Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Norman, B.H. *et al.*, *J. Med. Chem.* 2017; **60(1)**: 66-88.
2. Lane, N.E., *et al.*, *N. Engl. J. Med.* 2010; **363(16)**: 1521-1531.

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

Purification: Protein A affinity purification from HEK293 supernatants.

Background: Beta NGF is a neurotrophic factor (NTF) that is involved in growth, proliferation and survival of neurons. Anti-NGF monoclonal antibodies are in clinical trials as a chronic pain control medicine.

Application: Useful for blocking beta NGF binding to its receptor in biochemical and cellular assays.

Specific Activity

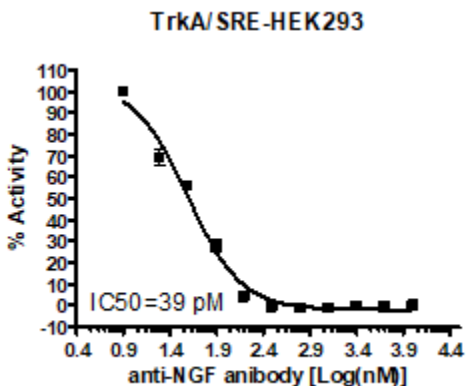


Fig.1 The Anti-NGF Antibody was evaluated using TrkA-SRE (Luc)- HEK293 cell line (BPS Bioscience #79798). NGF (500 pM at final concentration) was incubated with various concentrations of Anti-NGF Antibody in the assay medium (Opti-MEM) for 45 min. at room temperature and added to the cells. After 5 hour incubation at 37°C, ONE-Step™ Luciferase Assay System (BPS Bioscience #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 **1.858.202.1401** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com