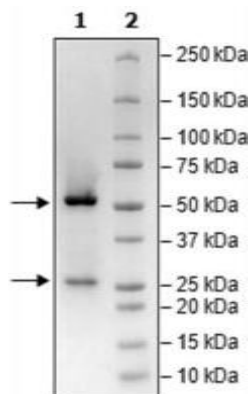


Product Information

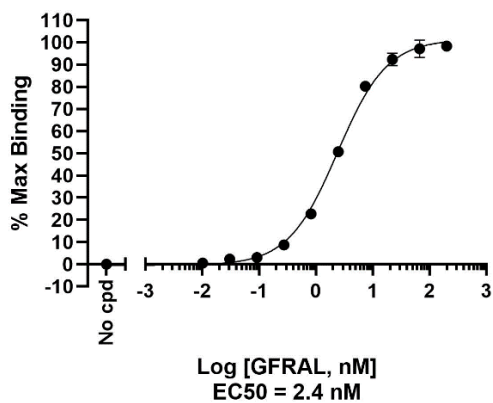
Description:	Recombinant anti-GFRAL (GDNF Family Receptor alpha Like) monoclonal antibody which recognizes GFRAL. This recombinant antibody has been tested for specific binding to purified recombinant human GFRAL (BPS Bioscience #101013) in an ELISA binding assay.
Species	Human
Concentration:	0.63 mg/ml
Clonality:	Monoclonal
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	Heavy Chain: 50 kDa; Light Chain: 24 kDa
Stability:	At least 12 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	An ELISA binding assay was performed with this antibody and its target protein, GFRAL (BPS Bioscience #101013). The antibody was coated onto an assay plate overnight at 2 µg/ml at 4°. On the next day, the plate was washed on blocked. Then, serial dilutions of the target protein were added (0 to 200 nM) for 1 hour at room temperature. After washing, HRP-Streptavidin was incubated for 30 minutes at room temperature. Following another wash step, colorimetric HRP substrate was added and quenched with 1N HCL. Then we measured absorbance on a plate reader at 450 nm. EC50 = 2.4 nM
Applications:	Useful for studying the binding and neutralization of GFRAL in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



GFRAL:Anti-GFRAL Binding Assay



Anti-GFRAL Neutralization Assay

