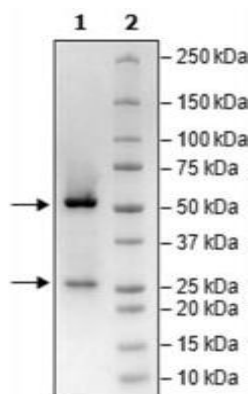


## Product Information

<b>Description:</b>	Recombinant anti-GFRAL (GDNF Family Receptor alpha Like) 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.
<b>Species</b>	Human
<b>Concentration:</b>	0.63 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	High Chain: 50 kDa; Low Chain: 24 kDa
<b>Stability:</b>	At least 12 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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.
<b>Assay Conditions:</b>	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
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



GFRAL:Anti-GFRAL Binding Assay

