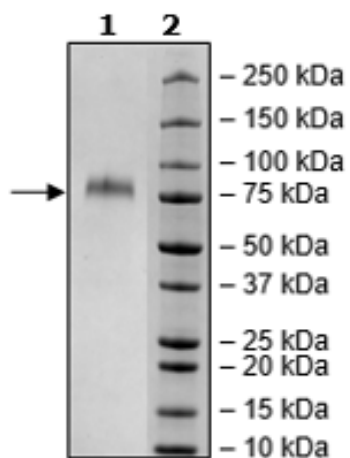


## Product Information

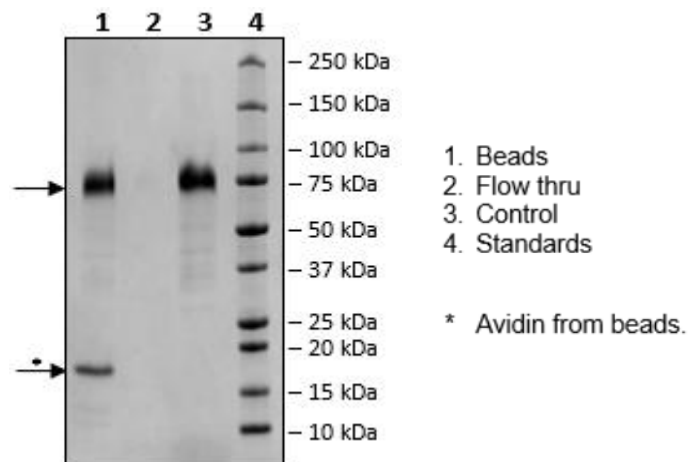
<b>Description:</b>	Recombinant human GFRAL (GDNF family receptor alpha like) encompassing amino acids 19-351(end). This protein is fused in C-terminal with the Fc fragment of human IgG1 followed by a C-terminal Avi-Tag™. The recombinant protein was enzymatically biotinylated using the Avi-Tag™ and affinity purified.
<b>Species</b>	Human
<b>Construct:</b>	GFRAL (19-351(end)-Fc(IgG1)-Avi)-(Biotin)
<b>Concentration:</b>	0.64 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>	67 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_207410.2
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
<b>Stability:</b>	At least 6 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>	<p>An ELISA binding assay was performed with GFRAL and GDF15. GDF15 was coated onto an assay plate overnight at 4 µg/ml at 4°. On the next day, the plate was washed on blocked. Then, serial dilutions of GFRAL 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. Absorbance was measured on a plate reader at 450 nm.</p> <p>A neutralization binding assay was performed with Anti-GFRAL Antibody (BPS Bioscience #101351) to test its ability to inhibit the binding of GFRAL (BPS Bioscience #101013) and GDF15. GDF15 was coated onto an assay plate overnight at 4 µg/ml at 4°. On the next day, the plate was washed on blocked. Then, serial dilutions of anti-GFRAL were added (0 to 200 nM) and allowed to preincubate for 30 minutes at room temperature before 1.5 ng/µl of GFRAL was added and continued to incubate for 1 hour. 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. Absorbance was measured on a plate reader at 450 nm.</p>
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

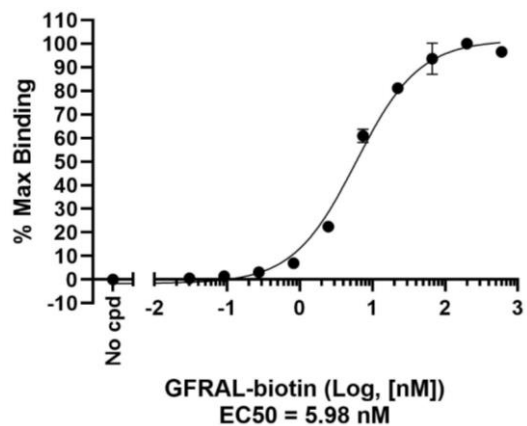
### 4-20% SDS-PAGE Coomassie Staining



### Biotin-Avidin Pulldown



### GDF15: GFRAL-Biotin Binding Assay



### Anti-GFRAL Neutralization Assay

