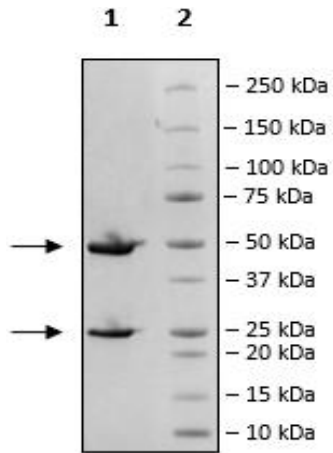


## Product Information

<b>Description:</b>	This anti-GPRC5D (G protein-coupled receptor class C group 5 member D) antibody is a purified, recombinant monoclonal antibody that recognizes human GPRC5D. The heavy chain construct includes a C-terminal Fc domain. This antibody has been tested for specific binding to human GPRC5D by flow cytometry.
<b>Background:</b>	GPRC5D (G protein-coupled receptor class C group 5 member D) is an orphan receptor of little-known function belonging to the large family of G protein-coupled receptors. The protein is highly expressed in malignant plasma cells such multiple myeloma and is now considered a therapeutic target for antibody-based treatment of multiple myeloma. For example, CAR-T cells and GPCR5DxCD3 bispecific antibodies resulted in promising outcomes at levels well tolerated by the patients.
<b>Species:</b>	Human
<b>Isotype:</b>	IgG4
<b>Clonality:</b>	Monoclonal
<b>Concentration:</b>	0.32 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>	Heavy Chain: 49 kDa + glycans; Light Chain: 24 kDa
<b>Glycosylation:</b>	This antibody runs at a higher MW by SDS-PAGE due to glycosylation.
<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>	One million of GPRC5D CHO cells (BPS Bioscience #78337) (green) and CHO-K1 cells (blue) were incubated with 1 µg of Anti-GPRC5D Antibody (BPS Bioscience #102000) for 30 minutes on ice. Cells were washed three times, then stained with Goat anti-Human IgG Fc Secondary Antibody, PE, eBioscience™ (Invitrogen #12-4998-82), washed three times, and analyzed by flow cytometry. The Y-axis represents the cell count. The X-axis indicates PE intensity.
<b>Applications:</b>	Useful for the flow cytometry analysis of human GPRC5D.

## Quality Control Data

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



### Flow Cytometry

