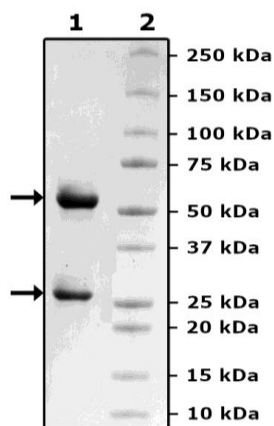


Product Information

Description:	Human recombinant anti-TFR (Transferrin Receptor) monoclonal antibody, which recognizes TFR. The purified antibody was enzymatically biotinylated using Avi-tag™ technology. This affinity purified antibody has been tested for specific binding to purified human TFR1 (BPS Bioscience #101287) in an ELISA binding assay.
Species	Human
Clonality:	Monoclonal
Concentration:	0.64 mg/ml
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 + glycans
Glycosylation:	This antibody runs at a higher MW by SDS-PAGE due to glycosylation.
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, TFR1 (BPS Bioscience #101287). The target protein was coated onto an assay plate overnight at 2 µg/ml at 4°. The next day, the plate was washed on blocked. Then, serial dilutions of the antibody 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 = 0.61 nM
Applications:	Useful for studying the binding of TFR in ELISA and in cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



TFR1:Anti-TFR (Biotin) Binding Assay

