

## Product Information

<b>Description:</b>	Recombinant human anti-TFR antibody designed towards human TFR (transferrin receptor). This anti-TFR antibody is a purified recombinant antibody. This antibody contains an Avi-Tag™ on the C-terminus of the Heavy Chain.
<b>Background:</b>	TFR (transferrin receptor protein-1), also known as TfR1 or CD71, is a cell surface receptor that binds holo-transferrin (Tf) with high affinity, forming an iron-Tf-TfR complex that is endocytosed. TFR1 is crucial for cellular health, as iron acts as a co-factor for many enzymes involved in metabolism and ROS (reactive oxygen species) formation. Deregulated iron levels, and consequently ROS formation, correlate with multiple disease states including Parkinson's Disease, Amyotrophic Lateral Sclerosis (ALS), Huntington's Disease, Multiple Sclerosis (MS), and tumorigenesis. TFR1 is upregulated on the surface of tumor cells and endothelial cells in the brain, presenting TFR1 as a promising candidate for targeted therapy in leukemia, lymphoma and neurological diseases. Studies of TFR1 function and expression patterns will broaden our knowledge and support possible applications in therapy.
<b>Species:</b>	Human
<b>Isotype:</b>	IgG1
<b>Clonality:</b>	Monoclonal
<b>Concentration:</b>	0.86 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: 51 kDa + glycans; Light Chain: 24 kDa
<b>Glycosylation:</b>	This antibody has 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>Applications:</b>	Useful for SDS-PAGE.

## Quality Control Data

## 4-20% SDS-PAGE Coomassie Staining

