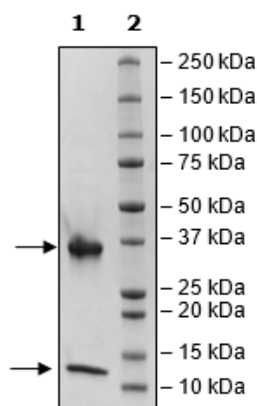


Product Information

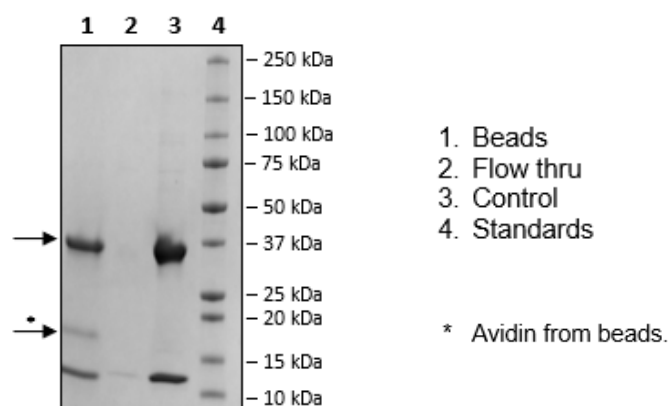
<b>Description:</b>	Recombinant human FcRn complex composed of FcRn (also known as Neonatal Fc receptor or FCGRT (Fc Fragment of IgG Receptor and Transporter)), encompassing amino acid 24-297 and B2M (beta 2 microglobulin), encompassing amino acids 21-119 respectively. FCGRT has a C-terminal His-tag (10xHis) followed by an Avi-tag. This protein is enzymatically biotinylated using Avi-Tag™ technology. These proteins are co-expressed in a HEK293 expression system and affinity purified. HiP™ indicates a high purity protein (≥90% pure) and less than 10% aggregation as measured by gel filtration.
<b>Species:</b>	Human
<b>Construct:</b>	FCGRT (24-297-His-Avi)-(Biotin) / B2M (21-119)
<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>	FCGRT: 34 kDa + glycans; B2M: 14 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Aggregation:</b>	<10%
<b>Genbank Accession:</b>	FCGRT: NM_001136019; B2M: NM_004048
<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.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown



Quality Control Data

Gel Filtration Curve

