

Product Information

Description:	Recombinant human complex of FcRn also known as Neonatal Fc receptor, heterodimer of 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). These recombinant proteins are co-expressed in a HEK293 expression system and affinity purified.
Species:	Human
Construct:	FCGRT ((24-297)-His) / B2M (21-119)
Concentration:	0.97 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:	FCGRT: 32 kDa + glycans; B2M: 14 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	FCGRT: NM_001136019; B2M: NM_004048
Stability:	At least 6 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.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

