

Product Information

Construct:	IL6R (20-365-Fc(IgG1))
Concentration:	3.46 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110mM NaCl, 2.2 mM KCl, and 20% glycerol.
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_000565
MW:	65 kDa +glycans
Purity:	≥90%
Notes:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Control Data

4-20% SDS-Page Coomassie Staining

