

Product Information

Construct:	TACI (23-166-Fc(IgG1)-Avi)
Concentration:	0.79 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM NaCl, 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_012452
MW:	45 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

