

Product Information

Construct:	IGF1R (His-960-1367(end))
Concentration:	0.34 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 150 mM Imidazole, 2.2 mM KCl, 0.04% Tween-20, 0.20% Triton X-100, 20% glycerol
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_000875
MW:	47 kDa
Purity:	80%

Quality Control Data

4-20% SDS-Page Coomassie Staining

