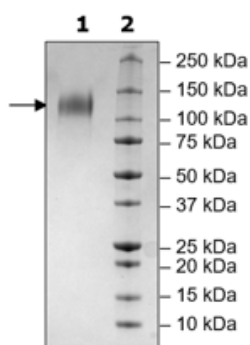


Product Information

Construct:	CD34 (32-290-Avi-His)
Concentration:	1.10 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, 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_001773
MW:	31 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Aggregation:	<10%

Quality Control Data

4-20% SDS-Page Coomassie Staining



Gel Filtration Trace

