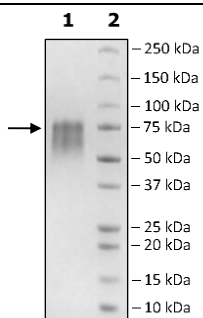


Product Information

Construct:	CD55 (35-353-Avi-His)
Concentration:	1.48 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_000574
MW:	38 kDa +glycans
Purity:	≥90%
Aggregation:	<10%
Notes:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Gel Filtration Trace

