

Product Information

Construct:	Angiopoietin 2 (19-496-His) (Rabbit)
Concentration:	0.43 mg/ml
Species:	Rabbit
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:	XM_002720896.2
MW:	58 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	79%

Quality Control Data

4-20% SDS-Page Coomassie Staining

