

Product Information

Construct:	ANGPT2 (mouse) (19-496-Avi-His)
Concentration:	0.66 mg/ml
Species:	Mouse
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_007426.4
MW:	58 kDa +glycans
Glycosylation:	This protein runs at higher M.W. by SDS-PAGE due to glycosylation.
Purity:	83%

Quality Control Data

4-20% SDS-Page Coomassie Staining

