

Product Information

Description:	Recombinant human Angiopoietin-2 fusion protein. This protein contains an N-terminal FLAG-tag followed by the oligomerization domain of COMP (a.a 28-72 of mouse COMP (Cartilage Oligomeric Matrix Protein, also designated thrombospondin 5); fused to amino acids 258-496(end) of human Angiopoietin-2.
Construct:	COMP(mouse)-ANGPT2(human) (FLAG-(28-72)-(258-496(end))).
Concentration:	0.41 mg/ml
Species:	ANGPT2: Human; COMP: 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:	ANGPT2: NM_001147; COMP: NM_016685
MW:	33 kDa
Purity:	84%

Quality Control Data

4-20% SDS-Page Coomassie Staining

