

Product Information

Construct:	Tau-381 (Full Length)
Concentration:	0.20 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 0.25 mM DTT, 0.1 mM PMSF, 25% glycerol
Expression System:	<i>E. coli</i>
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Uniprot:	P10636-4
MW:	54 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

