

Product Information

Construct:	14-3-3 zeta (His-Full Length)
Concentration:	0.20 mg/ml
Species:	Human
Formulated In:	50 mM sodium phosphate, pH 7.0, 300 mM NaCl, 150 mM imidazole, 0.1 mM PMSF, 0.25 mM DTT, 25% glycerol
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_003406
MW:	31 kDa
Purity:	90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

