

Product Information

Construct:	Shh (25-198-His) (Mouse)
Concentration:	0.27 mg/ml
Species:	Mouse
Formulated In:	8 mM Phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% 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
Genbank Accession:	NM_009170
MW:	21 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

