

Product Information

Construct:	Peroxiredoxin I (K197Q)(His-1-199(end))
Mutation:	K197Q
Concentration:	0.84 mg/ml
Species:	Human
Formulated In:	25 mM Tris-HCl, pH 8.0, 100 mM NaCl, 0.05% Tween-20, 50% Glycerol, 3 mM DTT.
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_002574
MW:	23 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

