

Product Information

Construct:	MAPKAPK2 (GST-Th-1-400(end))
Concentration:	1.09 mg/ml
Species:	Human
Formulated In:	40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 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_032960
MW:	72 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

