

Product Information

| | |
|---------------------------|---|
| Construct: | MKK6 (GST-2-334(end)) |
| Concentration: | 0.90 mg/ml |
| Species: | Human |
| Formulated In: | 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, 0.67 mM glutathione |
| 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_002758.3 |
| MW: | 64 kDa |
| Purity: | ≥90% |

Quality Control Data

4-20% SDS-Page Coomassie Staining

