

Product Information

Construct:	ATR (GST-2245-end)
Concentration:	0.05 mg/ml
Species:	Human
Formulated In:	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 0.1 mM PMSF, 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_001184
MW:	70 kDa
Purity:	75%

Quality Control Data

4-20% SDS-Page Coomassie Staining

