

Product Information

Construct:	BRPF3 (GST-576-701)
Concentration:	4.34 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20 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_015695
MW:	41 kDa
Purity:	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

