

Product Information

Construct:	ALK2(R206H)(GST-147-509(end))
Mutation:	R206H
Concentration:	1.04 mg/ml
Species:	Human
Formulated In:	20 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 2 mM GSH
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_001105.5
MW:	68 kDa
Purity:	90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

