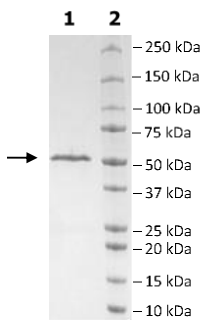


Product Information

Construct: FAK (GST-411-686)
Concentration: 0.61 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
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_153831
MW: 57 kDa
Purity: ≥90%
Specific Activity: 80 pmol/min/μg
Assay Conditions: Assay was done according to the FAK Kinase Assay Kit (BPS Bioscience #40722) with FAK titrated from 320 ng/rxn – 5 ng/rxn with 20 μM ATP instead of 10 μM ATP.
Applications: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

