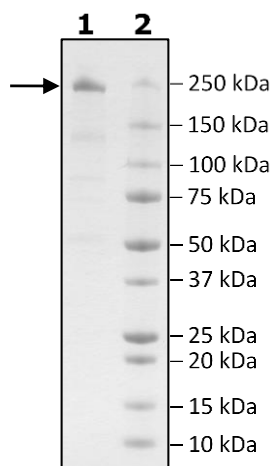


Product Information

Construct:	ACC1 (1-2383(end)-FLAG-His)
Concentration:	0.70 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 100 µg/ml FLAG peptide
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_198834
MW:	271 kDa
Purity:	74%
Specific Activity:	≥58 pmol/min/µg
Assay Conditions:	Assay was done according to the Acetyl-Coenzyme A Carboxylase 1 (ACC1) Assay Kit (BPS Bioscience #79315) with various amounts of ACC1.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

