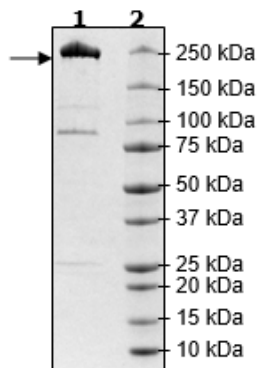


## Product Information

|                           |  |
|---------------------------|--|
| <b>Construct:</b>         | ACC2 (1-2458(end)-His)   |
| <b>Concentration:</b>     | 0.41 mg/ml   |
| <b>Species:</b>           | Human  |
| <b>Formulated In:</b>     | 40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, 100 mM imidazole   |
| <b>Expression System:</b> | Sf9  |
| <b>Format:</b>            | Aqueous buffer solution  |
| <b>Stability:</b>         | At least 6 months at -80°C. Avoid freeze/thaw cycles.  |
| <b>Storage:</b>           | -80°C  |
| <b>Genbank Accession:</b> | NM_001093  |
| <b>MW:</b>                | 277 kDa  |
| <b>Purity:</b>            | 86%  |
| <b>Specific Activity:</b> | 20 pmol/min/μg   |
| <b>Assay Conditions:</b>  | Assay was done using ADP-Glo (Promega) reagents with 20 μM ATP, 20 μM acetyl-CoA and 12 mM NaHCO <sub>3</sub> in 25 μL ACC assay buffer (30 mM HEPES pH=7.4, 2 mM MgCl <sub>2</sub> , 1 mM DTT, and 2 mM Sodium citrate). Reaction at room temperature for 40 min. |
| <b>Applications:</b>      | Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.  |

## Quality Control Data

4-20% SDS-Page Coomassie Staining



ACC2 Activity

