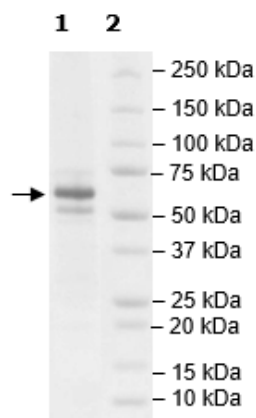


## Product Information

<b>Construct:</b>	CERK (1-531(end)-FLAG)
<b>Concentration:</b>	0.26 mg/ml
<b>Species:</b>	Rat
<b>Formulated In:</b>	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 100 µg/ml FLAG peptide
<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_00113486
<b>MW:</b>	61 kDa
<b>Purity:</b>	77%
<b>Specific Activity:</b>	14 pmol/min/µg
<b>Assay Conditions:</b>	CERK activity was measured at 37°C for 25 min in a buffer (20 mM HEPES pH7.2, 80 mM KCl, 3 mM CaCl <sub>2</sub> , 10 mM MgCl <sub>2</sub> , 2 mM DTT, 5 µM ATP) using a µ-Octylglucoside mixed micellar assay with 100 µM Ceramide (BPS Bioscience #40618). ATP reduction was detected using Kinase-Glo® Luminescent Kinase Assay Platform (Promega, Madison, WI).
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



CERK Activity

