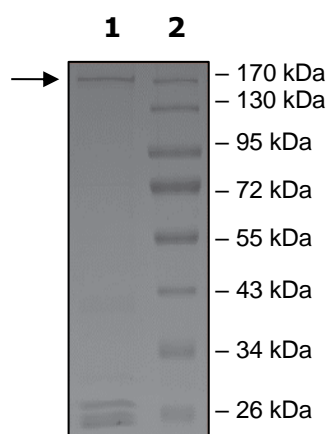


## Product Information

<b>Construct:</b>	LRRK2 (G2019S) (GST-968-end)
<b>Mutation:</b>	G2019S
<b>Concentration:</b>	0.10 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 0.1 mM PMSF, 25% glycerol.
<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_198578
<b>MW:</b>	210 kDa
<b>Purity:</b>	70%
<b>Specific Activity:</b>	4.2 pmol/min/μg
<b>Assay Conditions:</b>	Kinase activity was measured using ADP-Glo™ Kinase Assay kit (Promega; Cat# V9101) which quantifies the amount of ADP produced by the LRRK2 (G2019S) reaction. The ADP-Glo™ Reagent is added to terminate the reaction and quench the remaining ATP. The Kinase Detection Reagent is then added to convert ADP to ATP and to measure the newly converted ATP using a luciferase/luciferin reaction. Assay: Kinase activity was measured using substrate LRRKtide (RLGRDKYKTLRQIRQ) (stock 1 mg/ml). Increasing amounts of LRRK2 (G2019S) were incubated with a final concentration of 0.2 mg/ml substrate in 40 mM Tris-HCl, pH 7.4, 20 mM MgCl <sub>2</sub> , 0.1 mg/ml BSA, 50 μM fresh DTT, and 25 μM ATP. The total incubation time was 40 minutes at room temperature. The reaction was terminated by the addition of 5 μl of ADP-Glo™ Reagent and a subsequent incubation at room temperature for 40 minutes. Luminescence was measured by the addition of 10 μl Kinase Detection Reagent followed by incubation at room temperature for an additional 30 minutes. The blank was determined from a “no kinase” sample.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



Specific Activity

