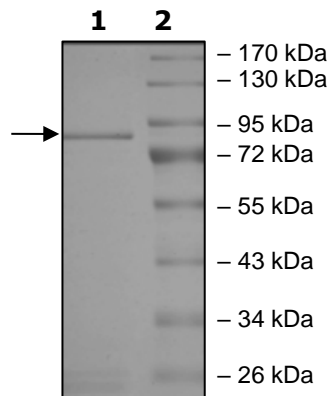


Product Information

Description:	Recombinant human ROS1 (G2101A) (ROS proto-oncogene 1, receptor tyrosine kinase), encompassing amino acids 1881-end. This construct contains an N-terminal GST-tag and was affinity purified.
Species:	Human
Construct:	ROS1 (G2101A) (GST-1881-end)
Mutation:	G2101A
Concentration:	0.10 mg/ml
Expression System:	Sf9
Purity:	90%
Format:	Aqueous buffer solution.
Formulated In:	50 mM Tris-HCl, pH 7.5, 300 mM NaCl, 10 mM glutathione, 0.1 mM EDTA, 0.25 mM DTT, 0.1 mM PMSF, 25% glycerol
MW:	82 kDa
Genbank Accession:	NM_002944
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Specific Activity:	131 pmol/min/μg
Assay Conditions:	<p>Kinase activity was measured using the ADP-Glo™ Kinase Assay Kit (Promega; Cat# V9101) which quantifies the amount of ADP produced. The ADP-Glo™ Reagent is added to terminate the reaction and deplete the remaining ATP. The Kinase Detection Reagent is then added to convert ADP to ATP and to measure the newly synthesized ATP using a luciferase reaction.</p> <p>ROS1 activity was measured by using IGFR1tide synthetic peptide (KKKSPGEYVNIEFG) diluted in water to a final concentration of 1 mg/ml. Reaction was initiated by mixing increasing amounts of the ROS1 with 25 μM ATP in 40 mM Tris-HCl, pH 7.4, 20 mM MgCl₂, 0.1 mg/ml BSA prepared with 250 μM DTT and substrate with a final concentration of 50 μg/ml.</p> <p>After a 40-minute incubation at 37°C, the reaction was terminated by addition of the AMP-Glo™ Reagent followed by a subsequent 40-minute incubation at room temperature. Kinase Detection Reagent was then added and incubated for another 30 minutes. Detection of luminescence was measured using the Luminescence Module Protocol on GloMax®-Multi Microplate reader. The corrected activity (RLU) was calculated by removing the blank value for each sample divided by the (specific activity of ADP in RLU/pmol)*(Reaction time in min)*(Enzyme amount in μg or mg). The blank was determined from a “no kinase” sample by replacing the enzyme working solution with an equal volume of Kinase Dilution Buffer IX (1X).</p>
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Specific Activity

