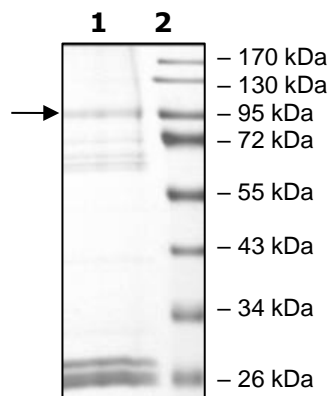


## Product Information

<b>Description:</b>	Recombinant human full-length RIPK3 (also known as receptor-interacting serine/threonine-protein kinase 3 or RIP3). This construct contains an N-terminal GST-tag. The recombinant protein was affinity purified and is active.
<b>Species:</b>	Human
<b>Construct:</b>	RIPK3 (GST-Full Length)
<b>Concentration:</b>	0.05 mg/ml
<b>Expression System:</b>	Sf9
<b>Purity:</b>	≥70%
<b>Format:</b>	Aqueous buffer solution.
<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>MW:</b>	96 kDa
<b>Genbank Accession:</b>	NM_006871
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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.
<b>Specific Activity:</b>	9.5 pmol/min/μg
<b>Assay Conditions:</b>	RIPK3 activity was measured by using MBP protein, diluted in 100 mM MOPS to a final concentration of 0.5 mg/ml, in the ADP Glo™ assay (Promega #V9101). Reaction was initiated by mixing increasing amounts of the RIPK3 with 25 μM ATP in 40 mM Tris-HCl, pH 7.4, 2.5 mM MgCl <sub>2</sub> , 0.1 mg/ml BSA prepared with 50 μM DTT, 2.5 mM co-factor MnCl <sub>2</sub> and substrate at a final concentration of 500 μg/ml final concentration. After a 40-minute incubation at Room Temperature, the reaction was terminated by addition of ADP-Glo™ Reagent, followed by a subsequent 40-minute incubation at room temperature. Kinase Detection Reagent was added, and the reaction was incubated for another 30 minutes at ambient temperature. Detection of luminescence was measured using the Luminescence Module Protocol on GloMax®-Multi Microplate Multimode Reader. The corrected activity (RLU) was calculated by removing the blank value for each sample. The Kinase Specific Activity was calculated as follows: $RLU / [(specific\ activity\ of\ ADP\ in\ RLU/pmol) * (Reaction\ time\ in\ min) * (Enzyme\ amount\ in\ \mu g\ or\ mg)]$ . The blank was determined from a “no kinase” sample by replacing the enzyme solution with an equal volume of Kinase Dilution Buffer X (1X).
<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

