

Data Sheet

RIPK1, GST-Tag (Mouse)

Mouse, Recombinant, N-terminal GST-Tag

Catalog #: 100340

Lot #: 190807-G2 **Conc.:** 1.6 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 122 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 3mM DTT and 20% glycerol.

Stability: At least 6 months at -80°C .
 Avoid freeze/thaw cycles.

References:

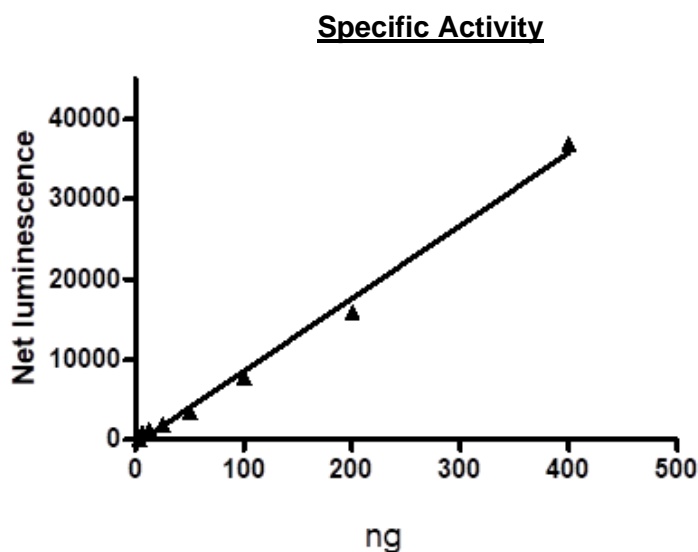
1. Hsu, H. et al., Immunity 4: 387-396, 1996.
2. Kim, J Y. et al., J Cell Sci. 2011 Feb 15;124(Pt 4):647-56.

Description: Mouse RIPK1, also known as RCK and MAP3K19, GenBank Accession No. NM_009068, a.a. 1-327, with N-terminal GST-tag, expressed in a Baculovirus-infected Sf9 cell expression system. MW = 64 kDa.

Assay Conditions: Assay was done in a Kinase buffer (BPS Cat. #79334) using MBP (0.2 mg/ml) as a substrate with 10 μM ATP at 30°C for 50 minutes. The amount of ATP transferred was calculated using ADP-Glo reagent (Promega).

Application: Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance



4-20% SDS-PAGE
Coomassie staining

Lane 1:
 3 μg RIPK1
Lane 2:
 Protein Marker

MW: 64 kDa

Purity: $\geq 90\%$

