

Data Sheet

STK33, GST-Taq

Human, Recombinant, N-terminal GST-tag
Catalog #: 40093
Lot#: 170929-G2 **Conc.:** 0.28 mg/ml

Formulated in: 40mM Tris-HCl pH 8.0, 110mM NaCl, 2.2mM KCl, 0.04% Tween-20, 3mM DTT and 20% glycerol.

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles. Protein may be diluted to ≥ 100 µg/ml in PBS + glycerol and stored at -80°C.

References:

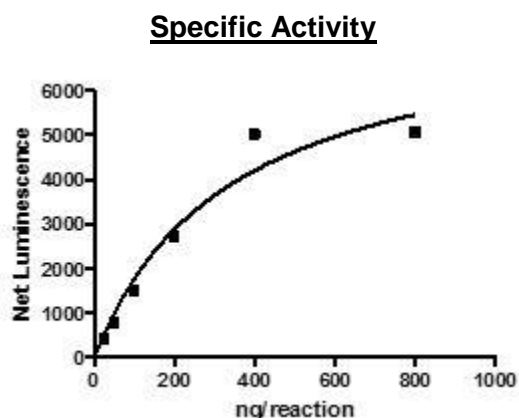
1. Mujica, A.O., *et al.*, 2001. *Gene* **280**: 175-181.
2. Brauksiepe, B., *et al.*, 2008. *Biochemistry* **9**: 25-37.

Description: STK33 (GenBank Accession No. NM_030906.2), amino acids 2-end, with N-terminal GST-tag, MW = 85 kDa, expressed in a Baculovirus-infected Sf9 cell expression system.

Specific Activity: ≥10 pmol/min/µg. Assay was done using ADP-Glo (Promega) reagents in Kinase buffer containing 2 mM DTT using MBP substrate (0.1 mg/ml) and 100 µM ATP. Reaction was done at 30°C for 40 min.

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

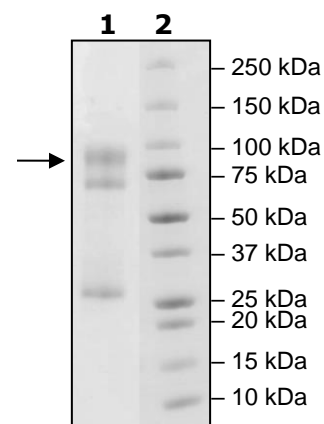
Quality Assurance



**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
2µg STK33
Lane 2:
Protein Marker

MW: 85 kDa
Purity: ≥46%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com