

Data Sheet

STK19, GST-Tag

Human, Recombinant, N-terminal GST-tag

Catalog #: 100910

Lot #: 200923

Conc.: 0.05 mg/ml

Formulated in: 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.

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

References:

1. Yin, C. *et al.*, *Cell*. **176(5)**: 1113-1127, 2019.
2. Qian, L. *et al.*, *Clin. Cancer Res.* **26(13)**: 3420-3430, 2020.

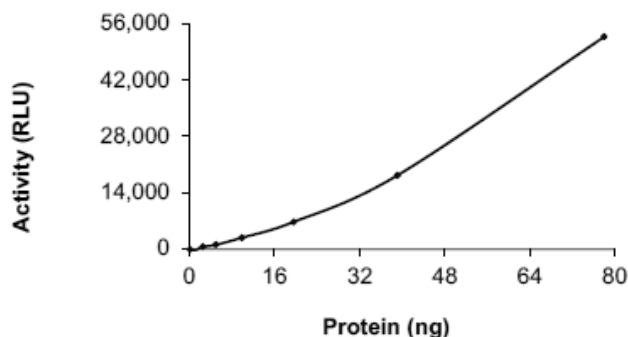
Description: Human Serine/threonine-protein kinase 19 (STK19), also known as G11 and RP1, GenBank Accession No. NM_004197, a.a. 1-368(full length), with N-terminal GST-tag, expressed in Sf9 insect cells via a baculovirus expression system. MW=70 kDa.

Specific Activity: 1.70 pmol/min/ μg
 Assay Conditions: STK19 was diluted with kinase dilution buffer and 1 mg/ml CREBtide (KRREILSRRPYSYR) and assayed according to the ADP-Glo™ Kinase Activity Assay Kit (Promega, #V9101).

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

Quality Assurance

Specific Activity



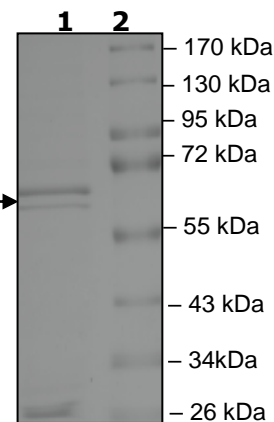
4-20% SDS-PAGE Coomassie staining

Lane 1:
STK19

Lane 2:
Protein Marker

Purity: >75%

MW: 70 kDa



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com