

## Recombinant SNK protein

**Catalog No:** 31194

**Expressed In:** Baculovirus

**Quantity:** 10 µg

**Source:** Human

**Buffer Contents:** 10 µg of Recombinant SNK protein in 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 4 mM reduced glutathione, 20% glycerol. Protein is supplied at 0.071 µg/µl.

**Protein Details:** Recombinant human SNK protein was produced using baculovirus infected Sf9 cells. The protein was made against amino acids M1-N685, accession number NM\_006622 and N-terminally fused to GST-HIS6-Thrombin site. Purified by GSH-agarose affinity purification.

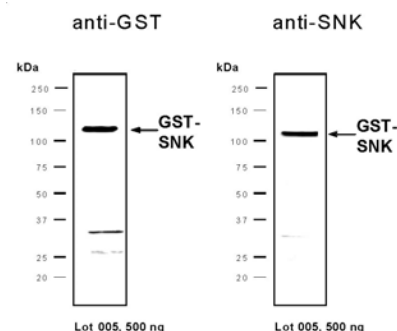
**Application Notes:** Recombinant SNK is suitable for kinase assays and Western blot. The molecular weight of the protein is ~107.435 kDa. The activity of the protein is ~ 2 pmol/µg min.

Recommended kinase reaction conditions: 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl<sub>2</sub>, 3 mM MnCl<sub>2</sub>, 3 µM Na-orthovanadate, 1.2 mM DTT, ATP (variable), 2.5 µg/50 µl PEG20.000, Substrate: Rb-CTF, 10 µg / 50 µl, Recombinant SNK: 200 ng/50 µl.

Kinase activity may vary depending on the substrate and reaction conditions used.

**Storage and Guarantee:** This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



### Recombinant SNK protein tested by SDS-PAGE gel.

0.5 µg of SNK protein was analyzed using SDS/PAGE followed by Western blot with either an anti-GSK or anti-SNK antibody.