



11526 Sorrento Valley Rd SteA2  
San Diego Ca 92121  
Tel: 1.858.202.1401  
1.858.829.3082  
Fax: 1.858.259.5734  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## Data Sheet

### CHK1, Active

human, recombinant, N-terminal GST tag  
Catalog #: 40039

Formulated in: 50mM Tris-HCl, pH 7.5,  
150mM NaCl, 0.25mM DTT, 0.1mM EGTA,  
0.1mM EDTA, 0.1mM PMSF, 25% glycerol.

Stability: >6 months at -80°C

References:

1. Walworth, N. et al. Nature. 1993 May 27;363(6427):368-71.
2. Walworth, NC. et al. Science. 1996 Jan 19;271(5247):353-6.

Description:

Recombinant full-length human CHK1 was expressed by baculovirus in Sf9 insect cells using a N-terminal GST tag. The gene accession number is NM\_001274.

Specific Activity:

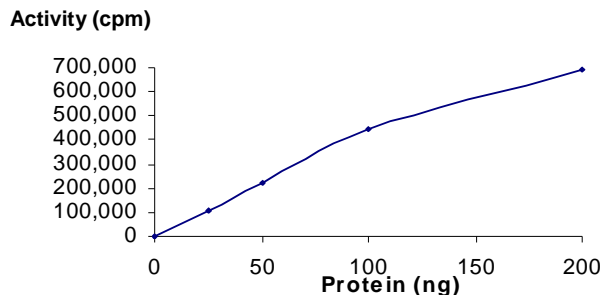
214 nmol/min/mg

Application:

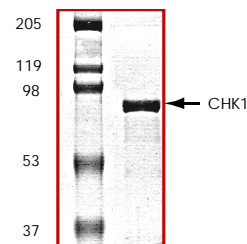
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### Quality Assurance

CHK1 Activity  
214 nmol/min/mg



10% SDS-PAGE  
Coomassie staining



MW: 82 kDa.

Purity: 95%

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** or **1.858.829.3082**, Fax **1.858.259.5734**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)