



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

Data Sheet

CHK2

human, recombinant, N-terminal HIS tag
Catalog #: 40040

Formulated in: 50mM Tris-HCl, pH 7.5,
150mM NaCl, 40% glycerol and 3mM DTT.

Stability: >12 months at -80°C

References:

1. Reinhardt HC and Yaffe MB. Curr Opin Cell Biol. 2009 Feb 19.
2. Stracker TH and Petrini JH. Cell Cycle. 2008 Dec;7(23):3618-21.

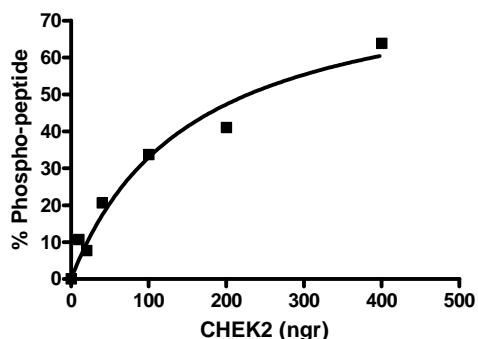
Description: Human Check point kinase 2 (GenBank Accession No. NM_007194), full length, with N-terminal HIS-tag. MW=91kDa. Expressed in a Baculovirus infected Sf9 cell expression system.

Specific Activity: 6.9 pmoles/min/μg. Assay conditions: CHEK2 was incubated with a substrate for 1h at RT in 1Xkinase buffer supplemented with ATP. Developer solution was added to reaction and reaction was stopped after 1h of incubation at RT.

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

Quality Assurance

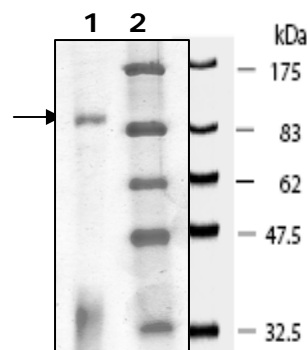
CHK2 Specific Activity



10% SDS-PAGE Coomassie staining

Lane 1:
2 μg CHK2
Lane 2:
Protein Marker
BioLabs (#P7708L)

MW: 91 kDa.
Purity: 50%



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

Please visit our website at: www.bpsbioscience.com