

PRODUCT INFORMATION



Chk1 (human, recombinant)

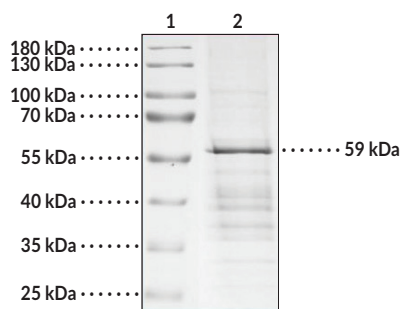
Item No. 43751

Overview and Properties

Synonyms:	Cell Cycle Checkpoint Kinase 1, Checkpoint Kinase 1, CHK1 Checkpoint Homolog, Serine/threonine-protein Kinase Chk1
Source:	Active recombinant human N-terminal His-tagged Chk1 expressed in insect cells
Amino Acids:	1-476 (full length)
Uniprot No.:	O14757
Molecular Weight:	59 kDa
Storage:	-80°C (as supplied)
Stability:	≥1 year
Purity:	≥80% estimated by SDS-PAGE
Supplied in:	Sterile 50 mM sodium phosphate, pH 7.5, with 300 mM sodium chloride, 50-200 mM imidazole, 0-0.1 mM PMSF, 0.2-1.0 mM DTT, and 10-25% glycerol
Bioactivity:	See figure for details
Specific Activity:	batch specific

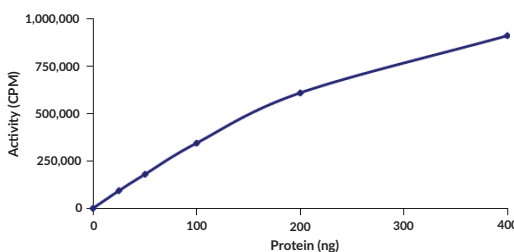
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: Chk1

SDS-PAGE Analysis of Chk1. This protein has a calculated molecular weight of 59 kDa.



The specific activity of Chk1 was determined to be 165 nmol/min/mg as per activity assay protocol.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/09/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Checkpoint kinase 1 (Chk1) is a serine/threonine kinase and downstream effector of ataxia-telangiectasia and RAD3-related protein/kinase (ATR) in the cellular DNA damage response.^{1,2} It is composed of an N-terminal kinase domain, a flexible linker region, a serine/glutamine cluster, and a C-terminal regulatory domain, which contains two conserved motifs, CM1 and CM2, that mediate kinase autoinhibition.^{2,3} Chk1 is ubiquitously expressed and localized primarily to the nucleus.^{3,4} In response to single-stranded DNA (ssDNA) generated from replication stress, DNA crosslinking, or other damage, ATR phosphorylates Chk1 at serine 345 and 317 inducing its activation and cell cycle arrest at the S or G₂/M phase.⁴ Loss of one copy of *Chk1* induces erythropoietic defects, anemia, and sudden death in mice, while *Chk1* knockout is embryonically lethal. Pharmacologic inhibition of Chk1 increases DNA damage and enhances antitumor efficacy of chemotherapeutic agents in various mouse xenograft models.⁵ Cayman's Chk1 (human, recombinant) protein can be used for enzyme activity assay and Western blot (WB) applications. This protein has a calculated molecular weight of 59 kDa.

References

1. Qiu, Z., Oleinick, N.L., and Zhang, J. ATR/CHK1 inhibitors and cancer therapy. *Radiother. Oncol.* **126**(3), 450-464 (2018).
2. Zhang, Y. and Hunter, T. Roles of Chk1 in cell biology and cancer therapy. *Int. J. Cancer* **134**(5), 1013-1023 (2014).
3. Day, M., Parry-Morris, S., Houghton-Gisby, J., et al. Structural basis for recruitment of the CHK1 DNA damage kinase by the CLASPIN scaffold protein. *Structure* **29**(6), 531-539 (2021).
4. Neizer-Ashun, F. and Bhattacharya, R. Reality CHEK: Understanding the biology and clinical potential of CHK1. *Cancer Lett.* **497**, 202-211 (2021).
5. Barnard, D., Diaz, H.B., Burke, T., et al. LY2603618, a selective CHK1 inhibitor, enhances the anti-tumor effect of gemcitabine in xenograft tumor models. *Invest New Drugs* **34**(1), 49-60 (2016).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM