

---

Product Manual

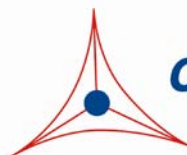
# 96-well Checkpoint Kinase Activity Assay Kit

## Catalog Number

STA-414	96 assays
STA-414-5	5 x 96 assays

**FOR RESEARCH USE ONLY**  
Not for use in diagnostic procedures

---



**CELL BIOLABS, INC.**  
*Creating Solutions for Life Science Research*

## **Introduction**

Cdc25C is a protein phosphatase responsible for dephosphorylating and activating Cdc2, a critical step in regulating mitosis in eukaryotic cells. Cdc25C is constitutively phosphorylated at Ser<sup>216</sup> throughout interphase by c-TAK1 (Cdc25C associated protein kinase), while phosphorylation at this site is DNA damage dependent at the G2/M checkpoint. Prior to mitosis, checkpoint kinases (CHK1 and CHK2) can be activated in response to DNA damage, phosphorylating Cdc25C at Ser<sup>216</sup>, ultimately leading to cell cycle arrest. Consequently, a cell is prevented from entering mitosis and passing this damage to daughter cells. Mutations to these checkpoint kinases lead to decreased DNA repair and increased susceptibility to cancer.

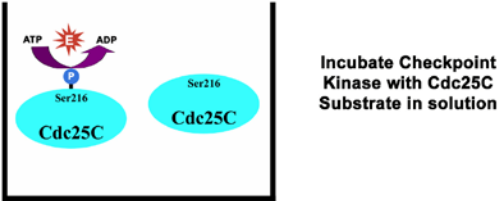
Cell Biolabs' 96-well Checkpoint Kinase Activity Assay Kit is an enzyme immunoassay developed for detection of the specific phosphorylation of Cdc25C at Ser<sup>216</sup> by Checkpoint Kinases (CHK1 and CHK2). First, the substrate is incubated with checkpoint kinase sample in solution (such as purified kinase, cell or tissue lysate). After the kinase reaction, the solution is transferred to a Kinase Substrate Capture Plate, where the phosphorylated Cdc25C is finally detected by an Anti-Phospho-Cdc25C (Ser<sup>216</sup>) Antibody (see Assay Principle).

Cell Biolabs' 96-well Checkpoint Kinase Activity Assay Kit provides a non-isotopic, sensitive and specific method to monitor checkpoint kinase activity using its physiological substrate; it can also be used in screening checkpoint kinase inhibitors. The kit has detection sensitivity limit of ~300 pg of active CHK1. A recombinant active CHK1 is also provided as a positive control. Each kit provides sufficient quantities to perform up to 96 assays.

## **Related Products**

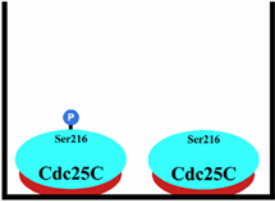
1. STA-320: OxiSelect™ Oxidative DNA Damage ELISA Kit (8-OHdG Quantitation)
2. STA-321: OxiSelect™ DNA Double-Strand Break (DSB) Staining Kit
3. STA-322: OxiSelect™ UV-induced DNA Damage ELISA Kit (CPD Quantitation)
4. STA-323: OxiSelect™ UV-induced DNA Damage ELISA Kit (6-4PP Quantitation)
5. STA-324: OxiSelect™ Oxidative DNA Damage Quantitation Kit (AP sites)
6. STA-326: OxiSelect™ Cellular UV-induced DNA Damage ELISA Kit (CPD)
7. STA-327: OxiSelect™ Cellular UV-induced DNA Damage Staining Kit (CPD)
8. STA-328: OxiSelect™ Cellular UV-induced DNA Damage ELISA Kit (6-4PP)
9. STA-329: OxiSelect™ Cellular UV-induced DNA Damage Staining Kit (6-4PP)
10. STA-413: Checkpoint Kinase Activity Immunoblot Kit
11. STA-415: ROCK Activity Immunoblot Kit
12. STA-416: 96-well ROCK Activity Assay Kit

# Assay Principle



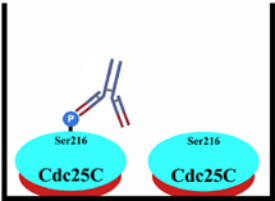
Incubate Checkpoint Kinase with Cdc25C Substrate in solution

1-2 Hrs at 30 °C



Cdc25C is captured by the Substrate Capture Plate

1 Hour at Room Temp



Wash and detect with Anti-Phospho-Cdc25C (Ser216) Antibody

1 Hour at Room Temp



Wash and add HRP Conjugated Secondary Antibody

1 Hour at Room Temp



Wash and add Substrate Solution for 5-20 Min, then read OD

## **Kit Components**

### **Box 1 (shipped at room temperature)**

1. Checkpoint Kinase Substrate Capture Plate (Part No 241401): One strip well 96-well plate.
2. 10X Checkpoint Kinase Substrate (Part No 241402): One vial – 1.5 mL containing recombinant Cdc25C.
3. 10X Kinase Buffer (Part No. 241602): One bottle – 20 mL of 250 mM Tris, pH 7.5, 100 mM MgCl<sub>2</sub>, 50 mM Glycerol-2-Phosphate, 1 mM Na<sub>3</sub>VO<sub>4</sub>.
4. ATP Solution (Part No. 241604): One vial – 100 µL of 100 mM ATP.
5. Anti-Phospho-Cdc25C (Ser216) Antibody (Part No. 241403): One vial – 20 µL.
6. Secondary Antibody, HRP Conjugate (Part No. 10902): One vial – 20 µL.
7. Assay Diluent (Part No. 310804): One 50 mL bottle.
8. 10X Wash Buffer (Part No. 310806): One 100 mL bottle.
9. Substrate Solution (Part No. 310807): One 12 mL amber bottle.
10. Stop Solution (Part. No. 310808): One 12 mL bottle.

### **Box 2 (shipped on blue ice packs)**

1. Active CHK1 (Part No. 241303): One vial – 20 µL containing 50 ng active CHK1 in 25 mM Tris, pH 7.5, 10 mM MgCl<sub>2</sub>, 5 mM Glycerol-2-Phosphate, 0.1 mM Na<sub>3</sub>VO<sub>4</sub>, 10% Glycerol, 0.1% BSA.

## **Materials Not Supplied**

1. Checkpoint kinase sample (purified kinase, cell or tissue lysate)
2. Lysis Buffer: 50 mM Tris-HCl, pH 7.5, 150 mM NaCl, 1 mM 2-glycerophosphate, 1 % Triton X-100 or 1 % Nonidet P-40, 1 mM EDTA, 1 mM EGTA, 1 mM Na<sub>3</sub>VO<sub>4</sub> and Proteinase inhibitors.
3. 1M DTT
4. 30°C incubator or water bath
5. 10 µL to 1000 µL adjustable single channel micropipettes with disposable tips
6. 50 µL to 300 µL adjustable multichannel micropipette with disposable tips
7. Multichannel micropipette reservoir\
8. Microplate reader capable of reading at 450 nm (620 nm as optional reference wave length)

## **Storage**

Store active CHK1 at -80°C, ATP Solution at -20°C and all other kit components at 4°C until their expiration dates. Avoid multiple freeze/thaw cycles.

## **Preparation of Reagents**

- 1X Kinase Buffer: Dilute the 10X Kinase Buffer Concentrate to 1X with deionized water. Stir to homogeneity.  
*Note: Do not dilute the entire bottle of 10X Kinase Buffer. The kinase reaction requires 10X Kinase Buffer concentrate. Prepare only the minimum volume of 1X for the experiment.*
- 10X Kinase Reaction Buffer/DTT/ATP: Just prior to usage, add DTT to a final concentration of 10 mM and ATP to a final concentration of 2 mM to the 10X Kinase Buffer. For Example, add 10  $\mu$ L of 1M DTT (not provided) and 20  $\mu$ L of 100 mM ATP solution to 970  $\mu$ L of 10X Kinase Buffer. 10X Kinase Reaction Buffer containing DTT and ATP may be stored at 4°C for short term (1-2 weeks).
- 1X Wash Buffer: Dilute the 10X Wash Buffer Concentrate to 1X with deionized water. Stir to homogeneity.
- Anti-Phospho-Cdc25C (Ser<sup>216</sup>) Antibody and HRP-Conjugated Secondary Antibody: Immediately before use dilute the Anti-Phospho-Cdc25C (Ser<sup>216</sup>) Antibody 1:1000 and HRP-Conjugated Secondary Antibody 1:1000 with Assay Diluent. Do not store diluted solutions.

## **Assay Protocol**

### **I. Kinase Reaction (samples should be assayed in duplicate)**

- 1a. For immunoprecipitations with anti-checkpoint kinase antibody: Kinase is first immunoprecipitated from cell or tissue lysate with an anti-checkpoint kinase antibody and Protein A/G bead slurry. Immediately before the kinase assay, wash the bead slurry once with 1X Kinase Buffer. Remove all supernatant and add 120  $\mu$ L of 1X Kinase Buffer. Next, add 15  $\mu$ L of 10X Kinase Reaction Buffer/DTT/ATP, mixing well (see Preparation of Reagents Section). Initiate the kinase reaction by adding 15  $\mu$ L of 10X Checkpoint Kinase Substrate.
- 1b. Purified Kinase or Cell Lysate: Purified kinase or cell lysate sample can be used directly in the kinase assay or further diluted with 1X Kinase Buffer. Add 120  $\mu$ L of checkpoint kinase sample to a microcentrifuge tube or microtiter plate. Next, add 15  $\mu$ L of 10X Kinase Reaction Buffer/DTT/ATP, mixing well (see Preparation of Reagents Section). Initiate the kinase reaction by adding 15  $\mu$ L of 10X Checkpoint Kinase Substrate.
2. (optional) Add 4  $\mu$ L of the provided active CHK1 and 116  $\mu$ L of 1X Kinase Buffer to a microcentrifuge tube or microtiter plate. Next, add 15  $\mu$ L of 10X Kinase Reaction Buffer/DTT/ATP, mixing well (see Preparation of Reagents Section). Initiate the kinase reaction by adding 15  $\mu$ L of 10X Checkpoint Kinase Substrate.
3. Incubate the tubes or plate at 30°C for 1-2 hours with gentle agitation.
4. For immunoprecipitated samples (step 1a), centrifuge tubes for 10 seconds at 12,000 x g.

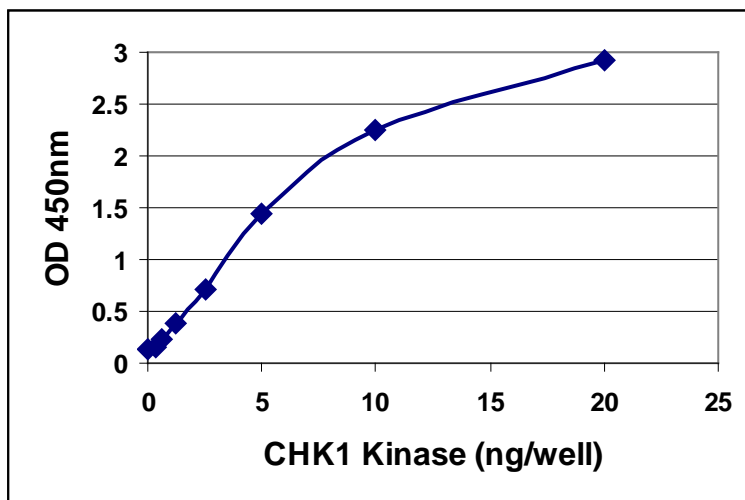
### **II. Transfer and Detection**

1. Transfer 100  $\mu$ L of kinase reaction sample to a well of the Checkpoint Kinase Substrate Capture Plate.
2. Cover with a plate cover and incubate at room temperature for 1 hour on an orbital shaker.

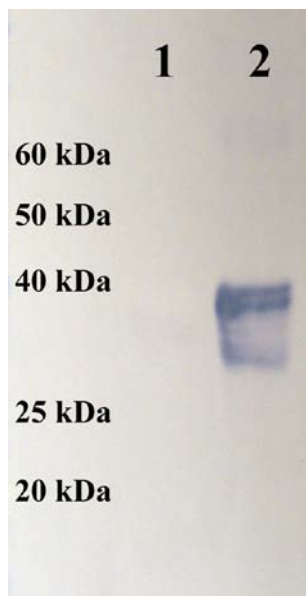
3. Remove plate cover and empty wells. Wash microwell strips once with 250  $\mu$ L 1X Wash Buffer. Empty the wells and tap microwell strips on absorbent pad or paper towel to remove excess 1X Wash Buffer.
4. Add 200  $\mu$ L of Assay Diluent to each well. Cover with a plate cover and incubate at room temperature for 30 minutes on an orbital shaker.
5. Remove plate cover and empty wells. Wash microwell strips 5 times with 250  $\mu$ L 1X Wash Buffer per well with thorough aspiration between each wash. After the last wash, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess 1X Wash Buffer.
6. Add 100  $\mu$ L of the diluted Anti-Phospho-Cdc25C (Ser<sup>216</sup>) Antibody to each well.
7. Cover with a plate cover and incubate at room temperature for 1 hour on an orbital shaker.
8. Remove plate cover and empty wells. Wash the strip wells 5 times according to step 5 above.
9. Add 100  $\mu$ L of the diluted Anti-Phospho-Cdc25C (Ser<sup>216</sup>) Antibody to each well.
10. Cover with a plate cover and incubate at room temperature for 1 hour on an orbital shaker.
11. Remove plate cover and empty wells. Wash the strip wells 5 times according to step 5 above.
12. Add 100  $\mu$ L of the diluted HRP-Conjugated Secondary Antibody to each well.
13. Cover with a plate cover and incubate at room temperature for 1 hour on an orbital shaker.
14. Remove plate cover and empty wells. Wash microwell strips 5 times according to step 5 above. Proceed immediately to the next step.
15. Warm Substrate Solution to room temperature. Add 100  $\mu$ L of Substrate Solution to each well, including the blank wells. Incubate at room temperature for 5-20 minutes on an orbital shaker.
16. Stop the enzyme reaction by adding 100  $\mu$ L of Stop Solution into each well, including the blank wells. Results should be read immediately (color will fade over time).
17. Read absorbance of each microwell on a spectrophotometer using 450 nm as the primary wave length.

## Example of Results

The following figure demonstrates typical results seen with Cell Biolabs' 96-well Checkpoint Kinase Activity Assay Kit. One should use the data below for reference only.



**Figure 1: CHK1 Kinase Activity Assay.** Active CHK1 was incubated for 2 hours at 30°C as described in the Assay Protocol.



**Figure 2: CHK1 Activity Immunoblot Assay.** 25  $\mu$ L of 1X Kinase Buffer containing 10 ng of active CHK1 was incubated with 50  $\mu$ L of 1X Kinase Buffer containing 0.2 mM ATP and 8  $\mu$ g of recombinant Cdc25C for 60 minutes at 30°C. Kinase reaction was stopped by adding 25  $\mu$ L of 4X SDS-PAGE Sample Buffer. **Lane 1:** Without kinase (negative control); **Lane 2:** with kinase. Phosphorylation of Cdc25C substrate was detected by anti-phospho-Cdc25C (Ser<sup>216</sup>) antibody as described in Assay Protocol for Cat. #STA-413.

## **References**

1. Zhou, B. B., and Elledge, S. J. (2000) *Nature* **408**, 433-439
2. Matsuoka, S., Huang, M., and Elledge, S. J. (1998) *Science* **282**, 1893-1897
3. Ahn, J., Urist, M., and Prives, C. (2004) *DNA Repair* **3**, 1039-1047
4. Ahn, J., and Prives, C. (2002) *J. Biol. Chem.* **277**, 48418-48426

## **Warranty**

These products are warranted to perform as described in their labeling and in Cell Biolabs literature when used in accordance with their instructions. THERE ARE NO WARRANTIES THAT EXTEND BEYOND THIS EXPRESSED WARRANTY AND CELL BIOLABS DISCLAIMS ANY IMPLIED WARRANTY OF MERCHANTABILITY OR WARRANTY OF FITNESS FOR PARTICULAR PURPOSE. CELL BIOLABS' sole obligation and purchaser's exclusive remedy for breach of this warranty shall be, at the option of CELL BIOLABS, to repair or replace the products. In no event shall CELL BIOLABS be liable for any proximate, incidental or consequential damages in connection with the products.

## **Contact Information**

Cell Biolabs, Inc.  
7758 Arjons Drive  
San Diego, CA 92126  
Worldwide: +1 858-271-6500  
USA Toll-Free: 1-888-CBL-0505  
E-mail: [tech@cellbiolabs.com](mailto:tech@cellbiolabs.com)  
[www.cellbiolabs.com](http://www.cellbiolabs.com)

©2011: Cell Biolabs, Inc. - All rights reserved. No part of these works may be reproduced in any form without permissions in writing.