

PhosphoSens-Lysate Assay Protocol & Validation Data for DNA-PK:

Determination of DNA-PK Activity in Crude Lysates from HCT-116 Cells Using the Selective AQT0440 Sensor Peptide Substrate

HGNC Name: PRKDC (DNA-PK)

Long Names: DNA-dependent protein kinase catalytic subunit

DNA-PK PhosphoSens-Lysate Kinase Discovery Kit



MATERIALS INCLUDED

TABLE 1

Components	Description	100 Assay Kit Volumes
PhosphoSens Kinase-Selective Lysate Substrate, AQT0440, 1mM	DNA-PK-selective sensor peptide substrate for assaying kinase activity in complex biological samples	45 µL
ATP Solution, 100 mM	100 mM ATP in nuclease-free water	30 µL
DTT Solution, 1 M	1 M DTT in nuclease-free water	5 µL
Enzyme Reaction Buffer (ERB), 10X	500 mM HEPES, pH 7.5, 0.1% Brij-35, 100 mM MgCl ₂	300 µL
Enzyme Dilution Buffer (EDB), 5X Base	20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 1 mg/mL Bovine Serum Albumin (BSA)	800 µL
EGTA Solution, 550 mM	550 mM EGTA in water (pH adjusted with NaOH)	30 µL
PhosphoPreserve Cell Extraction Buffer, 1X Base*	Please reach out to support@assayquant.com to inquire	1,000 µL
PhosphoPreserve Phosphatase Inhibitor Cocktail - 1, 100X	Please reach out to support@assayquant.com to inquire	20 µL
PhosphoPreserve Phosphatase Inhibitor Cocktail - 3, 60X	Please reach out to support@assayquant.com to inquire	20 µL
PhosphoPreserve Protease Inhibitor Cocktail, 60X	Please reach out to support@assayquant.com to inquire	20 µL
PhosphoSens PhosphoControl Peptide, AQT0442, 1 mM	Fully phosphorylated version of the DNA-PK-selective sensor peptide substrate, AQT0440.	10 µL
HCC1143 Lysate Control	Crude lysate of HCC1143 cells in PhosphoPreserve CEB, supplemented with DTT, and PhosphoPreserve Protease and Phosphatase inhibitor cocktails.	10 µL
DNA-PK Inhibitor: AZD-7648, 10 mM	AZD-7648 dissolved in 100% DMSO	5 µL

MATERIALS NOT INCLUDED

1. Ultra-pure deionized water
2. Dimethyl Sulfoxide (DMSO), (CH₃)₂SO
3. Precision pipettes capable of dispensing down to 0.5 µL and pipette tips. Having both single and multichannel pipettes is helpful
4. Plasticware: Low Protein Binding Microcentrifuge Tubes (0.5 and 1.5 mL), and materials to make your own cell or tissue lysates for a titration.
5. Centrifuge capable of spinning plates at 200xg and microfuge tubes at 10,000xg (standard microcentrifuge)
6. Fluorescence microplate reader capable of reading kinetically with filter setup for excitation (360 nm) and emission (485 nm) wavelengths. Alternatively, an instrument with a monochromator can be used to set the excitation (360 nm) and emission (485 nm) wavelengths, although this can reduce the assay sensitivity.
7. 96-well plate: We have validated Corning, half-area 96-well, white flat-bottom polystyrene NBS microplates (Cat. # 3642)
8. Plate Seal and Sealing Paddle: We have validated Dot Scientific, SealPlate 50um Polyester Sealing Film, Non Sterile (Cat. #T480).

*PhosphoPreserve Cell Extraction Buffer, 1X Base, requires protease and phosphatase inhibitors to preserve native kinase activity (prevent protein degradation and dephosphorylation). See notes in point 7 on the right side of this slide.

Lysate Assay Conditions and Reaction Setup

Assay Conditions:

54 mM HEPES, pH 7.5

1 mM ATP

1.2 mM DTT

0.012% Brij-35

1% glycerol

0.2 mg/mL BSA

0.54 mM EGTA

1 µg/mL Activated Calf Thymus DNA (used with recombinant DNA-PK only, not necessary with HCC1143 lysate)

10 mM MgCl₂

15 µM Sensor Peptide substrate, AQT0440 or Phosphopeptide control, AQT0442

Kinase Enzyme:

- 20 units/well DNA-PK Recombinant Enzyme (not provided with kit, but can be ordered from Promega (V5811, 2,500 units))
- **For simple lysate assay** – 2.2 µg/well HCC1143 lysates
- ***For Lysate titration** – 0, 313, 625, 1,250, 2,500, 5,000, 10,000, 20,000 ng/well lysate. Samples are prepared by serial dilution of the crude lysate.

Notes:

1. Total protein concentration for cell lysates was determined using a modified Bradford assay (Cat # 5000006, BIO-RAD).
2. **Cell Extraction Buffer (CEB):** 50 mM HEPES, pH 7.5, 150 mM NaCl, 2 mM EGTA, 1% Triton X-100, **supplemented with DTT, protease inhibitors, and phosphatase inhibitors 1 & 3** as indicated in slide 3 and as prepared on slides 5 and 11 **just before use**.
3. **Final 1X Enzyme Dilution Buffer (EDB):** 50 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 1 mg/ml Bovine Serum Albumin (BSA), supplemented with 0.2 mM EGTA and 1 mM DTT just before use.
4. Reactions were run in 25 µL final volume in Corning, low volume 384-well, white flat bottom polystyrene NBS microplates (Cat. #3824) after sealing using optically-clear adhesive film (TopSealA-Plus plate seal, PerkinElmer [Cat. #6050185] or Dot Scientific [Cat. #T480]) in a Biotek Synergy Neo2 microplate reader with filter setup for excitation (360 nm) and emission (485 nm) wavelengths.
 - Alternatively, reactions can be run in 50 µL in Corning, half-area 96-well, white flat-bottom polystyrene NBS microplates (Cat. # 3642).

****This test is performed with your cell lysates and not the samples provided. Lysates should be 1 mg/mL or higher total protein concentration. In the example experiment we ran and detailed, we started at 20,000 ng/well. This should be determined empirically and will vary depending on the cell line and the treatment conditions.***

Reagent Preparation for a Low Volume 384-well Format (25 μ L final reaction volume)*

*If you are working in a Corning 96-well plate (Cat. # 3642), multiply the volume of components by 2 for a final reaction volume of 50 μ L per well

- Using the stock solutions provided with the kit (Table 1), prepare the reagents shown in Table 2 and Table 3 (final concentrations shown in parentheses).
 - 10 mM ATP**: Make 90 μ L of 10 mM ATP by adding 9 μ L of 100 mM ATP to 81 μ L of ultrapure deionized water.
 - 10 μ g/mL Activated Calf Thymus DNA (DNA)***: Make 90 μ L of 10 μ g/mL DNA by adding 9 μ L of 100 μ g/mL to 81 μ L of ultrapure deionized water. ***not provided, but can be purchased from Cytiva if running a recombinant DNA-PK Assay**
 - 10 mM DTT**: Make 300 μ L of 10 mM DTT by adding 3 μ L of 1M (1000 mM) DTT to 297 μ L of ultrapure deionized water.
 - 5.5 mM EGTA**: Make 300 μ L of 5.5 mM EGTA by adding 3 μ L of 550 mM EGTA to 297 μ L of ultrapure deionized water.
 - 150 μ M Sensor Peptide**: Make 90 μ L of 150 μ M AQT0440 by adding 13.5 μ L of 1 mM AQT0440 to 76.5 μ L of ultrapure deionized water. Make 14 μ L of 150 μ M AQT0442 by adding 2.1 μ L of 1 mM AQT0442 to 11.9 μ L of ultrapure deionized water.
 - 50 μ M DNA-PK Inhibitor AZD-7648 (50X)**: Make 100 μ L of 50 μ M AZD-7648 by adding 0.5 μ L of 10 mM stock into 99.5 μ L DMSO.
- Prepare final 1X *PhosphoPreserve* Cell Extraction Buffer (CEB) by adding 1 μ L of 1M DTT, 16.7 μ L of the provided *PhosphoPreserve* Inhibitor Cocktail for Proteases (60X), 16.7 μ L of the provided *PhosphoPreserve* Inhibitor Cocktail for Phosphatases-3 (60X), and 10 μ L of the provided AQT Inhibitor Cocktail for Phosphatases-1 (100X) to the 1000 μ L of *PhosphoPreserve* Cell Extraction Buffer (CEB), Base. Keep on ice prior to extracting cells.
- Prepare '1.28X Master Mix' by combining volumes of the components listed in Tables 2 and 3. The volumes for a single well and 32 wells (Table 2) or 5 wells (Table 3) are shown.
- Prepare 624 μ L (sufficient for 32 wells but test requires only 28 wells and includes dead volume) of 1.28X Master mix per Table 2 and 97.5 μ L of 1.28X Master mix per Table 3.
- When adjusting the volume for a different number of wells, ensure that you include an additional 8% dead volume above the actual volume required.

Table 2 - Selective Sensor Peptide Substrate AQT0440

Components for 1.28X Master Mix:	For 1 Well:	For 32 Wells:
Enzyme Reaction Buffer (10X)	2.5 μ L	80 μ L
ATP (10 mM)	2.5 μ L	80 μ L
DTT solution (10 mM)	2.5 μ L	80 μ L
EGTA Solution (5.5 mM)	2.5 μ L	80 μ L
DNA (10 μ g/mL)* replace with water in lysate assay	2.5 μ L	80 μ L
Selective Sensor Peptide Substrate AQT0440 (150 μ M)	2.5 μ L	80 μ L
Ultrapure deionized water	4.5 μ L	144 μ L
Total volume	19.5 μL	624 μL

Table 3 - Phospho-Peptide Control AQT0442

Components for 1.28X Master Mix:	For 1 Well:	For 5 Wells:
Enzyme Reaction Buffer (10X)	2.5 μ L	12.5 μ L
ATP (10 mM)	2.5 μ L	12.5 μ L
DTT solution (10 mM)	2.5 μ L	12.5 μ L
DNA (10 μ g/mL)* replace with water in lysate assay	2.5 μ L	12.5 μ L
EGTA Solution (5.5 mM)	2.5 μ L	12.5 μ L
Sensor Phosphopeptide Control AQT0442 (150 μ M)	2.5 μ L	12.5 μ L
Ultrapure deionized water	4.5 μ L	35.0 μ L
Total volume	19.5 μL	85 μL

Step-by-Step Guide to Performing a Lysate Activity Assay in a 384-well Plate

A plate map for a simple lysate assay is shown on the next slide, which serves as a guide for making additions to the plate as outlined below.

1. Prepare 1X EDB using the 5X stock of EDB, Base provided, and supplement with DTT and EGTA. For example, to make 5000 μL of 1X EDB, add 1000 μL 5X EDB Base along with 5 μL of 1M DTT, 5 μL of 550 mM EGTA, and 3990 μL of ultrapure deionized water to create the final composition shown on slide 5. **Keep on ice.**
2. Prepare 250 μL BLANK by combining 249.571 μL 1X EDB and 0.429 μL of 1X Final PhosphoPreserve Cell Extraction Buffer .
3. *Prepare 40 μL of 20 units/ μL recombinant DNA-PK (5X) by adding 38 μL of 1X EDB to the 2.0 μL vial of the 80 units/ μL DNA-PK stock. **Keep on ice until needed.** **skip if not running a recombinant DNA-PK Assay with Promega DNA-PK*
4. Prepare 35 μL of 0.44 mg/mL lysate (5X) by adding 2.46 μL of the 6.25 mg/mL cobalt treated HCC1143 lysate stock to 32.54 μL of 1X EDB. **Keep on ice until needed.**
5. Transfer 20 μL of lysate to a new tube and heat at 95 $^{\circ}\text{C}$ for 5 minutes to serve as a Heat Inactivated (HI) negative control. Remove the tube and cool to room temperature.
6. Add 0.5 μL of 50 μM DNA-PK Inhibitor to all wells in columns 2, 5 and 7 (as shown on the plate map on the next slide). Add 0.5 μL of DMSO to all wells without the tool compound in columns 1, 3, 4, 6 and 8 (as shown on the plate map on the next slide).
7. Add 19.5 μL of 1.28X Master Mix for AQT0440 from Table 2 to all the wells except the wells labeled 'AQT0442'. For the wells labeled 'AQT0442', add 19.5 μL of 1.28X Master Mix for AQT0442 from Table 3.
8. Seal the plate using the plate seal supplied and press down with the supplied paddle. Incubate at 30 $^{\circ}\text{C}$ for 15 minutes to equilibrate the plate and Master Mix. This can be done by placing the plate inside a plate reader set at 30 $^{\circ}\text{C}$. This step is important to prevent temperature changes that can create anomalies in the data at the beginning of the reaction.
Note: The additions in step 8 and step 9 should be performed quickly since the reaction will start with these additions.
9. Source of Kinase Enzyme or control (BLANK or EDB, for background determination and the AQT0442 sensor phosphopeptide positive control):
 - Add 5 μL BLANK (prepared in step 2 above) to the wells labeled 'BLANK' or 1X EDB to wells labeled '1X EDB'
 - Add 5 μL 1X EDB to wells labeled 'AQT0442'
 - Add 5 μL of the HCC1143 lysate to the corresponding wells labeled '2.2 μg Lysate'
 - *Add 5 μL of 20 units/ μL Recombinant DNA-PK (5X) to the wells labeled '20 units DNA-PK'. **skip if not running a recombinant DNA-PK Assay*
10. Re-seal the plate, centrifuge at 200xg for one minute, and place the plate in the microplate reader set at 30 $^{\circ}\text{C}$.
11. Read the plate in kinetic mode with continuous fluorescence intensity detection (Ex/Em 360/485 nm) every 2 minutes for 1-4 hours. The frequency of the readings and the overall duration can be adjusted as needed.

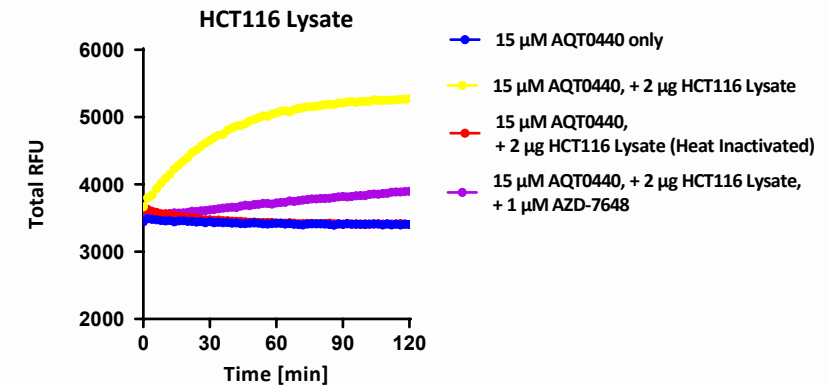
Plate Additions and Plate Map For a Simple Lysate Activity Assay

Component	Volume to add to wells
Tool Compound or 100% DMSO	0.5 μ L
Master mix	19.5 μ L
Lysate, DNA-PK, BLANK, or 1X EDB	5 μ L

Total volume 25 μ L

	Plate map for Lysate activity assay							
	HCC1143 Cell Lysate			With recombinant enzyme* optional*		EMPTY		
	No tool compound	With tool compound (AZD-7648)	Heat inactivated lysate	No tool compound	With tool compound (AZD-7648)	No tool compound	With tool compound (AZD-7648)	Phosphopeptide control
	1	2	3	4	5	6	7	8
A	BLANK	BLANK	BLANK	1X EDB	1X EDB	EMPTY	EMPTY	AQT0442
B	BLANK	BLANK	BLANK	1X EDB	1X EDB	EMPTY	EMPTY	AQT0442
C	2.2 ug Lysate	2.2 ug Lysate	2.2 ug HI lysate	20 Units DNA-PK	20 Units DNA-PK	EMPTY	EMPTY	AQT0442
D	2.2 ug Lysate	2.2 ug Lysate	2.2 ug HI lysate	20 Units DNA-PK	20 Units DNA-PK	EMPTY	EMPTY	AQT0442

Progress Curves for Total Fluorescence with AQT0440



AQT0442 data can be found on slide 10

This test uses 32 wells (28 wells for the AQT0440 sensor peptide substrate and 4 wells for the AQT0442 phosphopeptide control). All conditions are tested in duplicate.

Data Analysis

1. Collect the data from the microplate reader. This is a complete time course (Progress Curve) for every well with values in Relative Fluorescence Units (RFU) for each time point for Total (for each experimental condition) and “EDB/blank” wells.
2. Take the average of duplicate “EDB/blank” wells for each condition at each time point. Subtract the average EDB/blank values from the corresponding Total RFU of individual wells for each condition at each time point to obtain the background-corrected RFU values. For example, take the average of A1 and B1, and subtract the value from the total RFU determined for individual wells C1 and D1 at each time point. You can then either plot these RFU (Corrected) values separately to assess individual wells or take the average of the RFU (Corrected) values at each time point and plot this data.
3. It is highly recommended to run the “EDB/blank” wells at each compound concentration to correct for tool compound autofluorescence, if any. Since this is a kinetic assay format, the background with compounds will not change over time and can be subtracted from the total.
4. From the plot of the RFU (Corrected) values, determine the slope from the points in the linear region. This is the “initial reaction rate” in RFU (Corrected)/min. We recommend using ~30 minutes of the linear region of the progress curve to determine the rate. This can be performed in Excel, Excel-Fit, GraphPad Prism, the software provided with your microplate reader, or any other suitable software package, such as DynaFit, GeneData Screener, KinTek, Mathematica, MATLAB, or SigmaPlot.
5. Compare the RFU (Corrected)/min values for the samples to evaluate the activity of the kinase in each sample.
6. Refer to slide 10 for representative validation data for this simple lysate assay.

Data for a Simple Activity Assay Validation



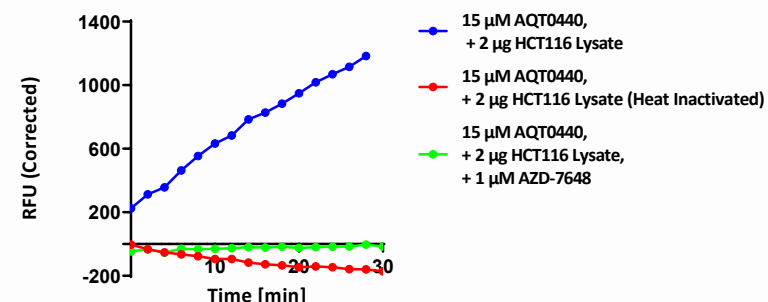
AssayQuant[®]
TECHNOLOGIES INC.

HCT116 Lysate or Recombinant DNA-PK with AQT0440 Sensor Peptide Substrate or AQT0442 Phospho-Control

PROGRESS CURVES

A. Crude Lysate Samples (2.0 µg/well)

1) Linear Time Course (0–30 min)

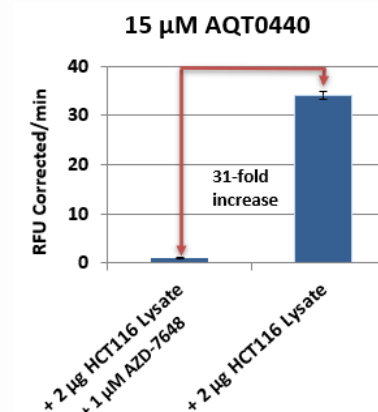


2) Reaction Rates to Assess Change

	Reaction Rate (RFU/min)	Change
2.0 µg HCT116 lysate (no treatment)	34 ± 0.82	
2.0 µg HCT116 lysate, Heat Inactivated	0 ± 0.34	100% inhibition DNA-PK Activity
2.0 µg HCT116 lysate, + 1.0 µM AZD-7648	1.1 ± 0.17	97% inhibition DNA-PK Activity

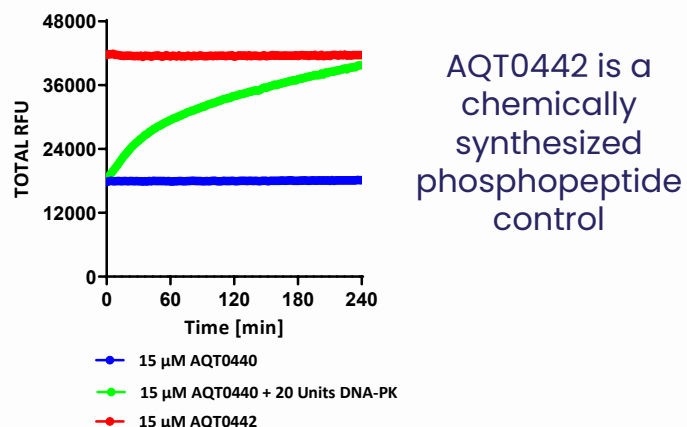
QUANTITATIVE ASSESSMENT

3) Histogram



B. Recombinant DNA-PK & AQT0442 Phospho-Control

1) Full Time Course (0–240 min.)



A. HCT116 Crude Lysate Sample: The AQT0440 sensor peptide was used to generate 1) A linear progress curve time course (0–30 min). The reaction rates (RFU Corrected values [Total – Background]/min. +/- standard deviations) are the slope of the linear region of each progress curve, which are presented in the table in 2) and as a histogram in 3), highlighting a 31-fold change in activity comparing the HCT116 Lysate to the HCT116 Lysate in the presence of AZD-7648. **Note:** The amount of activation depends on several factors, including cell type and passage number, the serum deprivation pretreatment used to make cells quiescent, as well as the nature, concentration, and duration of the activating stimulus or exposure to DNA damaging agent if applicable. These conditions can be varied to determine the effect on DNA-PK activity. The total amount of DNA-PK protein can be determined by Western Blotting or an ELISA; however, with the short stimulation times typically used, these levels are not expected to change.

B. Purified recombinant DNA-PK enzyme & AQT0442 Control: The DNA-PK protein complex (20 units) fully phosphorylated the AQT0442 sensor peptide substrate by 240 min., as shown by convergence with the signal obtained with the AQT0442 phosphopeptide positive control (a flat horizontal line defining the maximum RFU with this sensor peptide). The signal with DNA-PK enzyme was eliminated by adding 1 µM AZD-7648 inhibitor. The signal with AQT0442 can be used to convert RFU (Corrected) values to nmoles of Phosphate.

Reagent Preparation for a 384-well Format (25 µL final reaction volume) Lysate Titration*

*If you are working in a Corning 96-well plate (Cat. # 3642), multiply the volume of components by 2 for a final reaction volume of 50 µL. This test is performed with your cell lysates and not the samples provided. Lysates should be 1 mg/mL or higher total protein concentration

- Using the stock solutions provided with the kit (Table 1), prepare the reagents shown in Table 4 (final concentrations shown in parentheses).
 - 10 mM ATP:** Make 120 µL of 10 mM ATP by adding 12 µL of 100 mM ATP to 108 µL of ultrapure deionized water.
 - 10 mM DTT:** Make 300 µL of 10 mM DTT by adding 3 µL of 1M (1000 mM) DTT to 297 µL of ultrapure deionized water.
 - 5.5 mM EGTA:** Make 300 µL of 5.5 mM EGTA by adding 3 µL of 550 mM EGTA to 297 µL of ultrapure deionized water.
 - 150 µM Sensor Peptide:** Make 105 µL of 150 µM AQT1211 by adding 15.75 µL of 1 mM AQT0982 to 89.25 µL of ultrapure deionized water.
- Prepare final 1X *PhoshoPreserve* Cell Extraction Buffer (CEB) by adding 1 µL of 1M DTT, 16.7 µL of the provided *PhoshoPreserve* Inhibitor Cocktail for Proteases (60X), 16.7 µL of the provided *PhoshoPreserve* Inhibitor Cocktail for Phosphatases-3 (60X), and 10 µL of the provided *PhoshoPreserve* Inhibitor Cocktail for Phosphatases-1 (100X) to the 1000 µL of *PhoshoPreserve* Cell Extraction Buffer (CEB), Base. Keep on ice prior to extracting cells.
- Prepare lysate from cells, using sufficient cells to achieve a final concentration of 1 mg/mL of total protein determined using a modified Bradford assay (Cat # 5000006, BIO-RAD). See next slide for additional details.
- Prepare '1.25X Master Mix' by combining volumes of the components listed in Table 4. The volumes for a single well and 40 wells are shown.
- Prepare 800 µL (sufficient for 40 wells but this test requires only 36 wells and includes dead volume) of 1.25X Master mix per Table 4.
- When adjusting the volume for a different number of wells, ensure that you include an additional 8% dead volume above the actual volume required.

**Table 4 – Selective Sensor Peptide Substrate
AQT0440**

Components for 1.25X Master Mix:	For 1 Well:	For 40 Wells:
Enzyme Reaction Buffer (10X)	2.5 µL	100 µL
ATP (10 mM)	2.5 µL	100 µL
DTT solution (10 mM)	2.5 µL	100 µL
EGTA Solution (5.5 mM)	2.5 µL	100 µL
Selective Sensor Peptide Substrate AQT0440 (150 µM)	2.5 µL	100 µL
Ultrapure deionized water	7.5 µL	300 µL
Total volume	20 µL	800 µL

5.0 µL of crude lysate diluted in 1X EDB or BLANK (2:1 ratio of 1X EDB: 1X Final *PhoshoPreserve* Cell Extraction Buffer) alone for the blanks is added to each well (see slides 12 - 14 for detailed protocol).

Preparation of Crude Cell Lysates

Example: Preparing HCT116 Cell Lysates

Generate a lysate from cells using sufficient cells to achieve a final concentration of at least 1 mg/mL of total protein.

1. Cell Culture and Stimulation:

- a. 4.9×10^6 HCT116 cells were plated in 40 mL in a T-175 tissue culture-treated flask and incubated at 37 °C in McCoy's 5A Medium with 10% FBS and 1% PenStrep in an atmosphere of 5% CO₂.
- b. Cells at ~70% confluency (~ 16×10^6 cells) were washed with PBS, and the residual liquid was aspirated.
- c. Cells were then detached in 20 mL DPBS. Rotate the plate to cover the cells and use a cell scraper to ensure all cells are detached, followed by a pipette tip to wash the surface several times and break up any clumps of cells. Centrifuge the cells to a pellet, and aspirate the residual liquid.

2. Lysate Preparation:

- a. Cells were lysed in 600 µL of ice-cold 1X *PhosphoPreserve* CEB supplemented with DTT and protease and phosphatase inhibitors just before use per slide 11.
- b. Collect the lysate into a 1.5 mL microcentrifuge tube and then break up the DNA strands, if necessary, by passing through a 22-gauge needle 3 times or briefly sonicating on ice for 2 seconds on low power, followed by a 5 min spin at 10,000xg in a microcentrifuge at 4 °C. Remove and retain the supernatant and keep on ice. Determine the protein concentration and then immediately set up a lysate activity assay to determine the linear range. Alternatively, make aliquots and then snap freeze in liquid nitrogen or using dry ice in ethanol, and store at -80 °C.

This procedure yielded 800 µL per well of 4.5 mg/mL total protein. This can be scaled up or down with larger/smaller flasks or plates. The yield may vary slightly with cell size or passage number. The number of freeze-thaws should be minimized.

Step-by-Step Guide to Performing a Lysate Activity Assay Titration in a 384-well plate

A plate map for lysate dose-dependent assay is shown on the next slide, which serves as a guide for making additions to the plate as outlined below.

This test is performed with your cell lysates and not the samples provided. Lysates should be 1 mg/mL or higher total protein concentration

1. Transfer 20 μ L of 1.25X Master Mix per well to Rows A-C (triplicate samples) of the assay plate utilizing a multichannel pipette.
2. Seal the plate using the optically clear plate seal supplied and press down with the supplied paddle. Incubate at 30°C for 15 minutes to equilibrate the plate and Master Mix. This can be done by placing the plate inside a plate reader set at 30 °C. This preincubation is performed just prior to adding the 5 μ L of sample.
3. Prepare 1X EDB using the 5X stock of EDB, Base provided, and supplement with DTT and EGTA. For example, to make 5000 μ L of 1X EDB, add 1000 μ L 5X EDB Base along with 5 μ L of 1M DTT, 5 μ L of 550 mM EGTA, and 3990 μ L of ultrapure deionized water to create the final composition shown on slide 5. **Keep on ice.**
4. Prepare 60 μ L of 0.50 mg/mL of your stimulated lysate (for example, combine 30 μ L of a 1 mg/ml lysate stock with 30 μ L of 1X EDB; more concentrated lysate can be used but using lysate at less than 1 mg/mL is not recommended due to possible interference of lysate buffer in the assay when using higher amounts of lysate). **Keep on ice until needed.**
5. Prepare a 1:1 mixture of 1X EDB and CEB (50% EDB/50% lysate buffer) for diluting the lysate prepared above, by mixing 200 μ L of CEB with 200 μ L of 1X EDB (If the lysate is more concentrated, adjust this buffer to match the composition of the buffer used for the diluted lysate prepared in step 4).
6. Utilizing a separate lysate dilution plate, add 30 μ L of 50% EDB/50% lysate buffer per well to wells 1-11 in a single row. Add 30 μ L of the stimulated lysate prepared in step 4 to well 11. Add the remainder of the stimulated lysate to well 12.
7. Mix the contents of well 11 and transfer 30 μ L to well 10. Mix the contents of well 10 and transfer 30 μ L to well 9. Repeat this procedure down to well 2. Well 1 will be used for the “no enzyme” blank and receive only the 50% EDB/50% lysate buffer.
8. Transfer 5 μ L per well from the lysate dilution plate to rows A-C of the assay plate.
9. Centrifuge the plate at 200 x g for one minute, reseal, and place in the microplate reader set at 30 °C.
10. Read the plate in kinetic mode with continuous fluorescence intensity detection (Ex/Em 360/485 nm) every 2 minutes for 1-4 hours.

Note: The additions in step 8 should be performed quickly since the reaction will start with these additions.

Data analysis is performed as described on slide 9.

Plate Additions and Plate Map for a Lysate Titration

Component	Volume to add to wells
Master Mix	20 μ L
Lysate	5.0 μ L

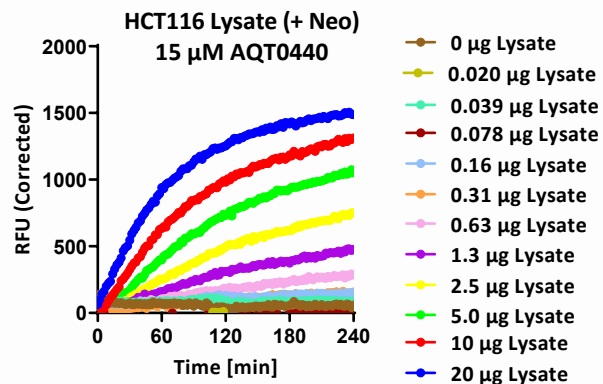
Total volume 25 μ L

Plate Map for Lysate Titration to Assess Dose Dependence													
		1	2	3	4	5	6	7	8	9	10	11	12
HCT116 Cell Lysates (ng)/well	A	0	2.4	4.9	9.8	20	39	78	156	313	625	1250	2500
	B	0	2.4	4.9	9.8	20	39	78	156	313	625	1250	2500
	C	0	2.4	4.9	9.8	20	39	78	156	313	625	1250	2500

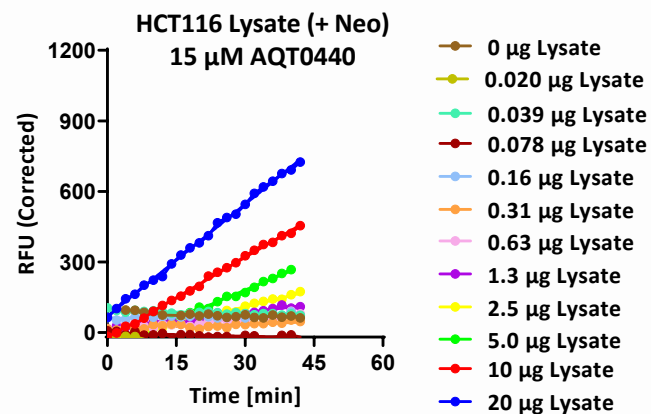
This test uses 36 wells and only the AQT0440 sensor peptide substrate and the lysate from HCT116 cells. Column 1 is the Blank. All conditions are tested in triplicate.

Lysate Titration for HCT116 Cells and DNA-PK Activity Measured with AQT0440

1. Full Time Course Progress Curves

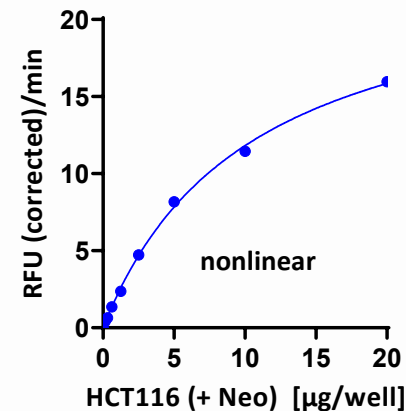


2. Linear Range of Progress Curves

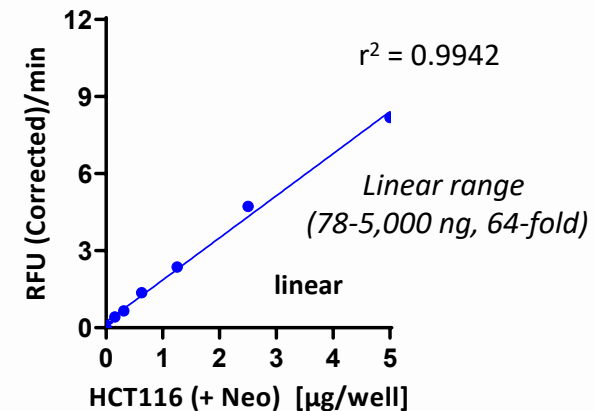


3. Reaction Rate vs Amount of Lysate per Well

A. All Lysate Amounts 15 μ M AQT0440



B. Lysate Amounts in Linear Range 15 μ M AQT0440



The AQT0440 sensor peptide was used at 15 μ M with an increasing amount of lysate from HCT116 cells treated for 1 hour with 0.50 μ g/mL Neocarzinostatin to induce DNA breaks. This experiment has also been run with untreated HCT116 Cell Lysate with similar results. RFU Corrected values (Total – Background) were determined for each condition. The results are presented for each amount of lysate for 1) Full time course of each progress curve (0-240 min.), and 2) Linear range of each progress curve, which was used to determine the slope for each amount of lysate. The results were then plotted as Reaction rates (RFU Corrected/min. +/- standard deviations) for all lysate amounts 3A), or those within the linear range as determined by an r^2 value > 0.99. Having the concentration of crude lysate samples at 1 mg/mL or higher, ensures that the amount of CEB in the reaction is minimized, even at the highest concentrations to avoid any inhibition of the kinase activity that can reduce the linear range.

The PhosphoSens-Lysate kinase activity assay provides a selective, highly quantitative, and accurate measure of kinase activity in a complex sample.