

# **PhosphoSens® Cell Lysate Activity Assay Format**

## **DNA-PK Assay Validation Using the AQ0440 Selective Sensor Peptide**

HGNC Name: PRKDC (DNA-PK)

Long Names: DNA-dependent protein kinase catalytic subunit

INTRODUCING A NEW INITIATIVE FROM OUR DIRECT-TO-PHASE II NIH/NCI SBIR AWARD (\$2M):

# PhosphoSens-Lysate

 PhosphoSens-Lysate assays are here!

## Experience the Benefits of Running Kinase Activity Assays in Complex Samples

Study Kinase Enzyme Activity Where Biology Happens

Simply Lyse Cells & Read  
No Fractionation Needed



**PhosphoSens-Lysate Assays**  
transform kinase activity  
measurement by enabling  
continuous, real-time kinetic analysis  
directly in cell and tissue lysates. This  
innovative approach provides key  
insights into your target's native  
environment throughout the drug  
discovery pipeline.

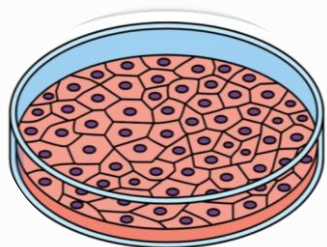
**NEW PRODUCTS** – Available Now for ERK1/2, AKT1/2/3, GSK3 $\alpha/\beta$ , JNK1-3, and DNA-PK!

**COMING SOON** – From our active R&D pipeline of selective sensor peptide substrates for:  
p38A/B/G/D, CDK1-3/5, CDK4/6, MEK1/2, SGK1/2/3, and PIM1/3.

# PhosphoSens-Lysate Assays

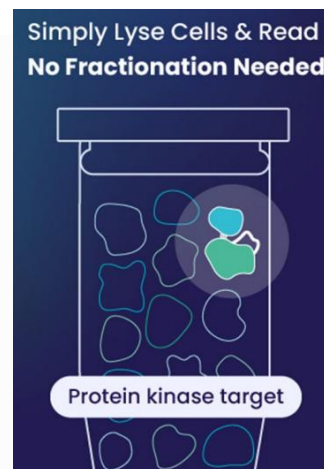
## A Simple & Powerful Solution to a Complex Problem

1. Grow Cells  
(+/- Pathway Activation)  
or Access Tissues



2. Harvest

ADD **PhosphoPreserve**  
Lysis Buffer + Protease &  
Phosphatase Inhibitors



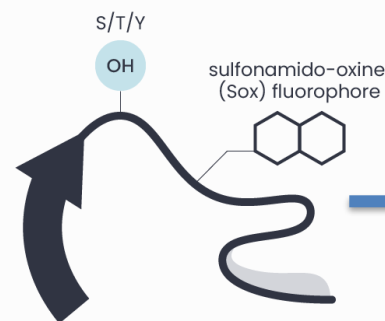
## Study Kinase Enzyme Activity Where Biology Happens

We have combined innovative PhosphoSens detection technology, invented at MIT, with high-throughput peptide synthesis methods to design sensor peptide substrates that are highly selective for the target of interest. This development integrates the advantages of the PhosphoSens platform (activity-based, direct, homogeneous, continuous/kinetic, quantitative) with the enabling capability of measuring endogenous target kinase activity in unfractionated cell or tissue lysates.

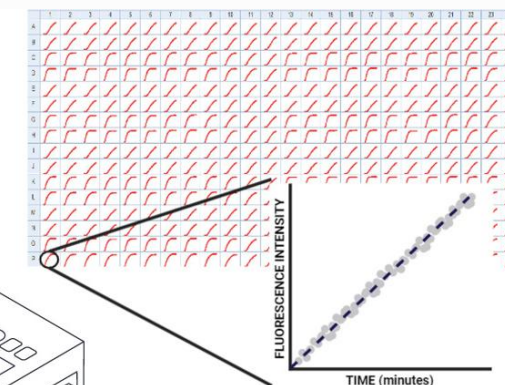
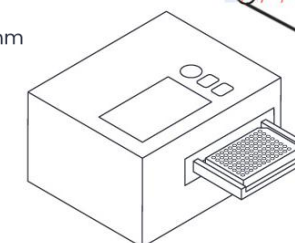
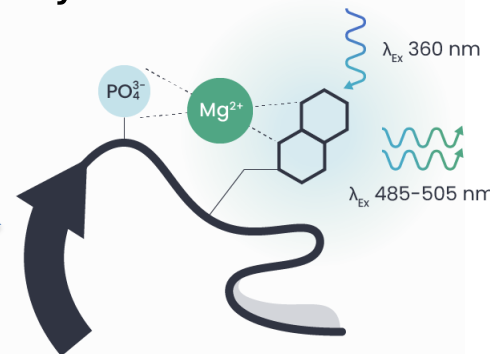


3. Prepare Reaction Mix  
(PhosphoSens Sensor Peptide  
Substrate, ATP, Mg<sup>2+</sup>)

Easy-to-use Kits &  
Bulk Sensor Peptide  
(Native substrates, stable,  
scalable, excellent lot consistency)

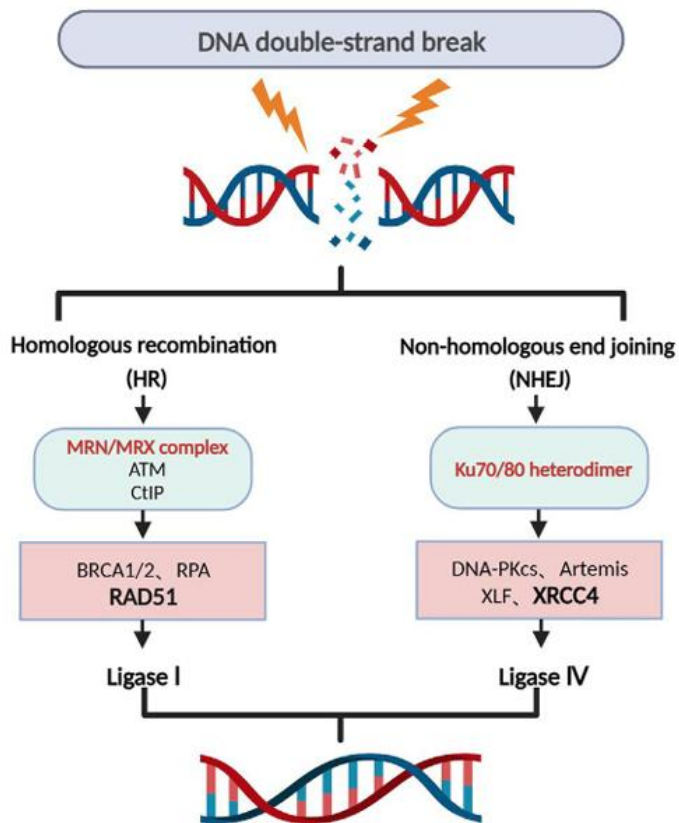


ADD Lysate & Read

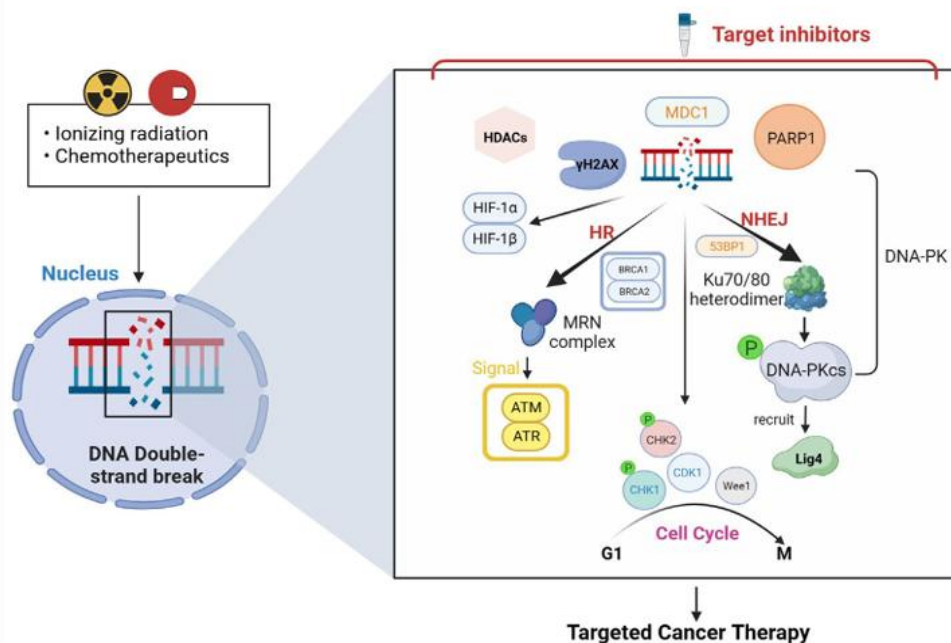




# DNA-PK Plays a Critical Role in the Repair of Double-strand DNA Breaks from Ionizing Radiation Used in Cancer Radiotherapy



## Cancer



**Fig. 1 of Jinpeng Tan, et al. 2023:** Double-strand break (DSB), a significant DNA damage brought on by ionizing radiation, acts as an initiating signal in tumor radiotherapy, causing cancer cells death. The two primary pathways for DNA DSB repair in mammalian cells are nonhomologous end joining (NHEJ) and homologous recombination (HR), which cooperate and compete with one another to achieve effective repair.

**Fig. 9 of Jinpeng Tan, et al. 2023:** HR and NHEJ are the two primary repair procedures in the DSB damage repair pathway. DNA double-strand breaks can be harmed by radiation or toxins. Pathway proteins that are important in DSB damage repair include the MRN complex, BRCA1/2, ATM/ATR, Ku70/80 heterodimer, DNA-PKcs, Lig4, MDC1, 53BP1, HDACs, HIF-1, PARP1, and H2AX. The creation of inhibitors to enable tailored therapy for the relevant tumors might conceivably use these proteins as targets.

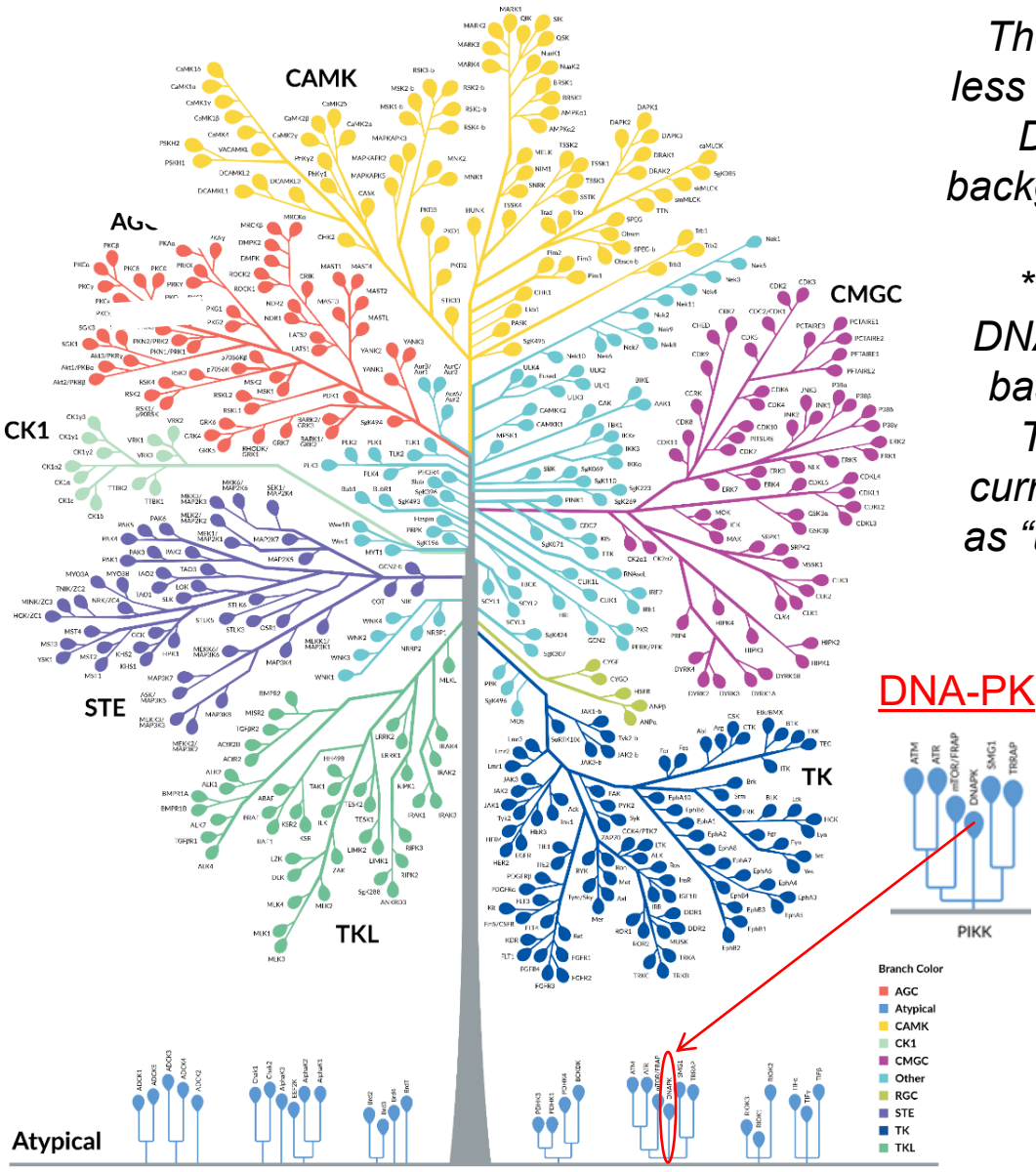
*Chrome-extension://efaidnbmninnibpcapjpcglclefindmkaj/  
<https://pmc.ncbi.nlm.nih.gov/articles/PMC10556206/pdf/MCO2-4-e388.pdf>*

*DNA-PK, ATM, ATR, CDK4/6, CDK1, WEE1, and CHK1/2 are all high priority targets for lysate assays*

# High Selectivity for DNA-PK with the AQT0440 Sensor Peptide

## Top 25 Hits with AQT0440 in Kinome Profiling with 411 Wild-type Kinases

Kinase target	Enzyme conc.	Average (RFU/pmol/min)	Rank	Selectivity Ratio	% Activity
DNA-PK	10	2176	1	1	100
CAMK1B	1.0	140	2	16	6.42
YES	1.0	125	3	17	5.73
CK1G1	0.80	92	4	24	4.22
CK1G2	0.80	71	5	31	3.24
CK1G3	0.80	64	6	34	2.95
TSSK2	2.0	53	7	41	2.45
SMG1	8.3	42	8	51	1.95
RSK4	1.0	40	9	54	1.85
CAMK1G	2.7	37	10	59	1.69
SRC	0.32	35	11	63	1.59
FAM20C	10	27	12	81	1.23
MARK4	1.0	23	13	94	1.07
SRPK2	2.0	23	14	96	1.04
TBK1	0.50	23	15	96	1.04
ATR-ATRIP	10	20	16	111	0.90
FGR	5.0	18	17	123	0.82
ZAP70	0.60	16	18	139	0.72
IKKB	8.0	15	19	148	0.67
ATM	10	14	20	151	0.66
CDK15/B1	4.1	13	21	162	0.62
GLK/KHS2	2.0	13	22	163	0.61
P70S6KB	4.6	11	23	200	0.50
TLK1	15	9.32	24	234	0.43
CDK9/T2	1.8	9.30	25	234	0.43

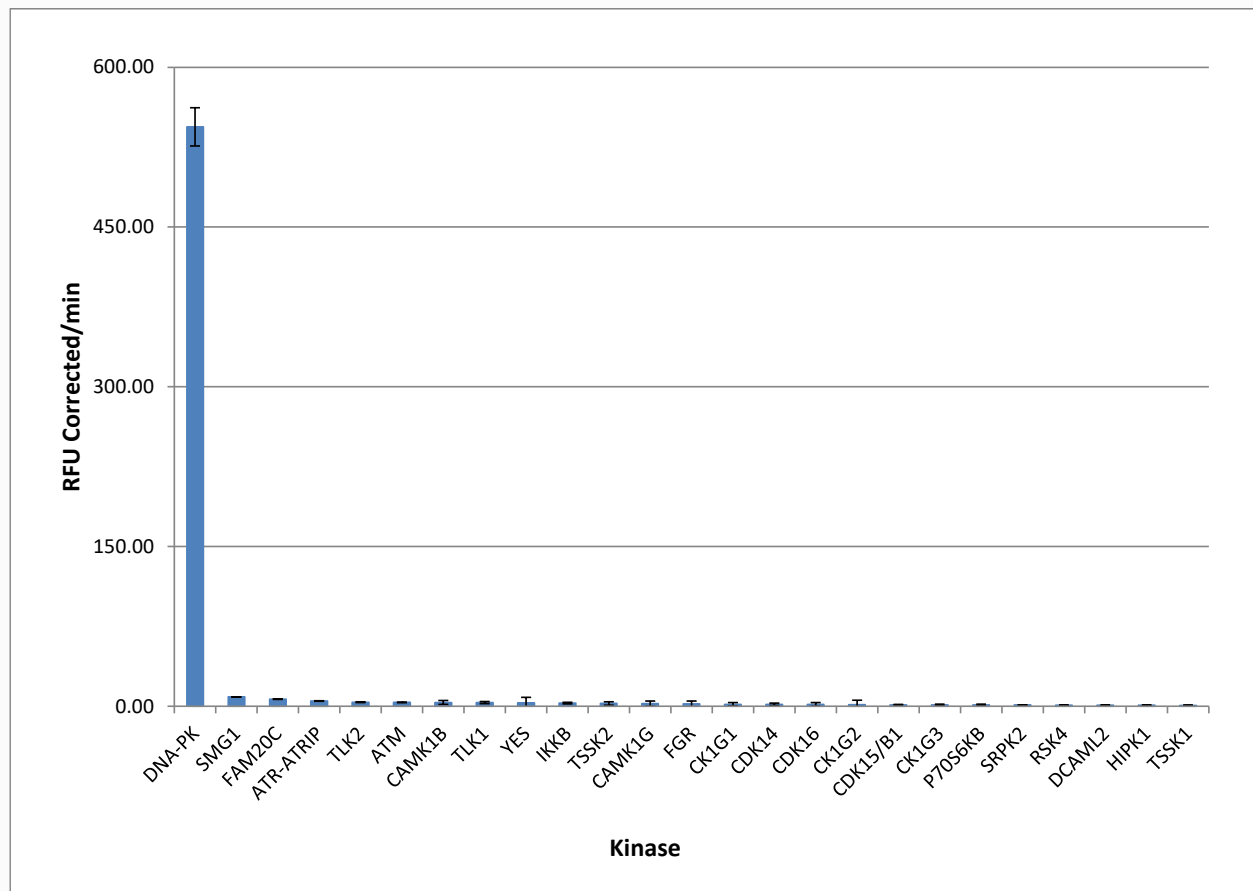


The off-target kinases have less than 6.5% of the activity of DNA-PK, which is in the background range (see slide 6).

\*The nM concentration of DNA-PK has been estimated based on a previous batch. The amount of DNA-PK currently provided is reported as “units” rather than “mg/mL” by the vendor.

# Selectivity of AQT0440 Retested Manually with the Top 25 Kinases from AQT's Kinome Profiling

Top 25 Kinases from the Profiling Run of 411 Kinases



*High-throughput  
Milestone Optimization  
led to the identification  
of the AQT0440 sensor  
peptide substrate.*

*The top 25 off-target  
kinases were reassessed  
manually and showed  
only 1.7% or less of the  
DNA-PK signal.*

*AQT0440 shows exquisite selectivity for the target kinase, a critical requirement for accurately measuring the activity of DNA-PK in crude lysates*

# Outline for this Study

## PhosphoSens-Lysate Assay Validation

### Lysate Source:

- ❖ HCT116 Cells +/- 0.50 to 1.0 µg/mL Neocarzinostatin (MedChemExpress, HY-111183; resuspended in 100% DMSO) for 1 hour or 100 µM Cobalt Chloride (Alfa Aesar; resuspended in water) for 48 hours

### Reference Compound Information:

- ❖ DNA-PK Inhibitor AZD-7648 (Cayman Chemical Co., 28598; resuspended in 100% DMSO)

### Experimental Validation at AssayQuant:

- ❖ Treated HCT116 cell lysate titration
- ❖ AQT0440 substrate  $K_m$  determination
- ❖ DMSO Tolerance Test
- ❖ Reference Compound  $IC_{50}$  Determination
- ❖ Full-length DNA-PK with Ku domain (Promega, V5811), to serve as a positive control
- ❖ Phosphopeptide control (AQT0442)
- ❖ Assessment of DNA-PK Activity in Multiple Cell Lines +/- Neocarzinostatin or Cobalt Treatment



# Preparation of Crude Cell Lysates from HCT116 Cells Treated ±Neocarzinostatin or Cobalt to Promote Double-Strand DNA Breaks

- 1) HCT116 Cells (passage 3) were plated in T-175 flasks and grown to 75% confluency over 48 hours at 37°C in McCoy's 5A Medium (ATCC, 30-2007) with 10% FBS (ATCC, 30-2020) and 1% PenStrep (30-2300) in an atmosphere of 5% CO<sub>2</sub>. Cells were then serum-starved in culture medium with 0.1% FBS for 24 hours and then incubated for 1 hour with or without 0.50 to 1.0 µg/mL Neocarzinostatin (MedChemExpress, HY-111183; resuspended in 100% DMSO) to promote double-strand DNA breaks. Alternatively, after 48 hours of normal cell growth, cells were incubated with 100 µM Cobalt Chloride (Alfa Aesar, 10692; resuspended in water) for 48 hours to generate reactive oxygen species (ROS), which also promote double-strand DNA breaks. Cells were placed on ice, then washed with cold PBS and lysed with 500 µL of cold Cell Extraction Buffer (CEB) as follows:

## Cell Extraction Buffer (CEB) with Protease and Phosphatase Inhibitors

- 50 mM HEPES, pH 7.4
- 150 mM NaCl
- 2 mM EGTA
- 1 mM DTT
- 1% Triton X-100
- 0.2 mM Na<sub>4</sub>P<sub>2</sub>O<sub>7</sub>
- 100 µM Na<sub>3</sub>VO<sub>4</sub>
- 0.2 mM NaF
- 50 mM β-glycerophosphate
- Protease Inhibitor Cocktail diluted 60-fold into lysis buffer
- Phosphatase Inhibitor Cocktail diluted 60-fold into lysis buffer

- 2) Cells were triturated to ensure solubilization and then lysates were centrifuged for 5 minutes at 12,000 x g to pellet cellular debris. Lysate supernatants were removed, aliquoted, supplemented with 10% glycerol, and used immediately or frozen at –80 °C, with each aliquot **used only once (one freeze/thaw cycle)**. Long-term storage of DNA-PK Lysates is not recommended, *as a loss in activity was observed over time.*

**Note:** The DNA-PK complex is a holoenzyme composed of the Ku heterodimer (Ku70 and Ku80) and a DNA-dependent Protein Kinase catalytic subunit (DNA-PKcs), a serine/threonine protein kinase. Ku binds to the ends of DNA double-strand breaks and recruits and activates DNA-PKcs to carry out its function in DNA repair. Because of the large size of this complex, measuring endogenous DNA-PK activity in lysates provides many advantages.



# Lysate Titration with the AQT0440 Sensor Peptide



## Reaction Conditions and Set Up

### Reaction 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 calf thymus DNA

10 mM MgCl<sub>2</sub>

15 µM AQT0440 sensor peptide substrate

0, 0.020, 0.039, 0.078, 0.16, 0.31, 0.63, 1.3, 2.5, 5.0, 10, and 20 µg/well **crude cell lysate** from HCT116 cells treated with 0.5 µg/mL Neocarzinostatin

### Reaction Set Up:

20 µL Reaction Mix with AQT0440, ATP, & DTT

Seal plate and incubate at 30 °C for 15 minutes to equilibrate

5 µL Enzyme dilution buffer (EDB) with DNA-PK Lysate Buffer (1x) or DNA-PK Lysate in Lysate Buffer (5x in EDB)

25 µL Final reaction volume

Reaction was run at 30 °C for 240 minutes in either Corning, low volume 384-well, white flat-bottom polystyrene NBS microplates (Cat. #3824) at 20 or 25 µL final well volume or in in PerkinElmer, ProxiPlate-384 Plus, white shallow well microplates (Cat. #6008280) at 20 µL final well volume after sealing using optically-clear adhesive film (TopSealA-Plus plate seal, PerkinElmer [Cat. #6050185]) in a Biotek Synergy Neo 2 microplate reader with excitation (360 nm) and emission (485 nm) wavelengths.

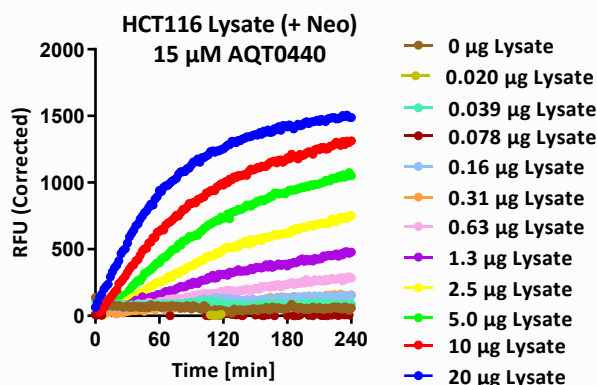
### Notes:

Enzyme Dilution Buffer (EDB): 20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 0.5 mM EGTA, 1 mM DTT, 1 mg/ml Bovine Serum Albumin.

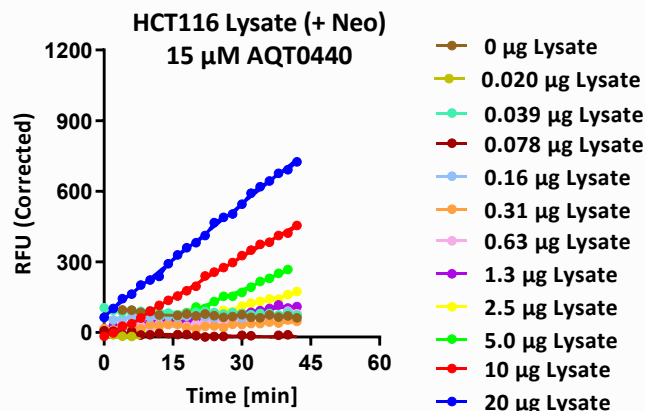
# Lysate Titration with the AQT0440 Sensor Peptide

## Progress Curves

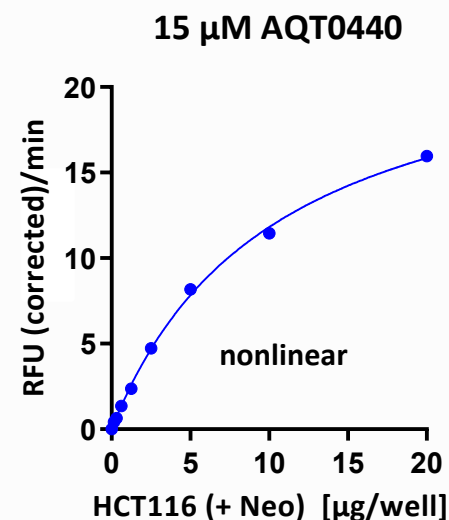
### Complete Progress Curves



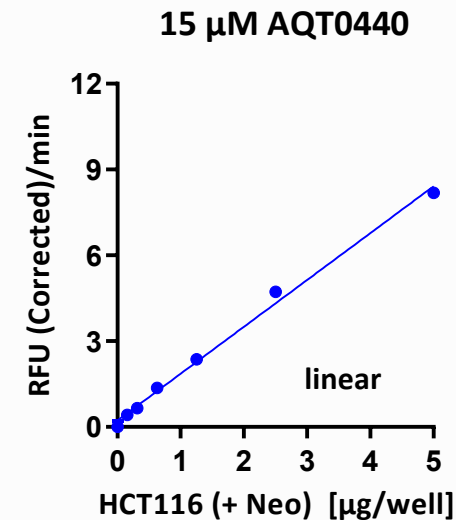
### Linear Region of Progress Curves



### Nonlinear Reaction



### Linear Reaction



The DNA-PK Lysate Assay is linear from 0.078 – 5.0  $\mu$ g/well of lysate (64-fold)

# DNA-PK Sensor Peptide $K_m$ Determination

## Reaction Conditions and Set Up

### Reaction 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 calf thymus DNA

10 mM  $MgCl_2$

0, 0.20, 0.39, 0.78, 1.6, 3.1, 6.3, 13, 25, 50, and 100 µM

**AQT0440 sensor peptide substrate**

4.0 µg/well NIH-3T3 crude cell lysate (from cells treated with 0.5 µg/mL Neocarzinostatin)

### Reaction Set Up:

2.5 µL 10X AQT0440 Substrate dilutions

17.5 µL Reaction Mix with ATP & DTT

Seal plate and incubate at 30 °C for 15 minutes to equilibrate

5 µL Enzyme dilution buffer (EDB) with lysate buffer(1x) or Lysate in lysis buffer (5x in EDB)

25 µL Final reaction volume

Reaction was run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat-bottom polystyrene NBS microplates (Cat. #3824) at 20 or 25 µL final well volume or in in PerkinElmer, ProxiPlate-384 Plus, white shallow well microplates (Cat. #6008280) at 20 µL final well volume after sealing using optically-clear adhesive film (TopSealA-Plus plate seal, PerkinElmer [Cat. #6050185]) in a Biotek Synergy Neo 2 microplate reader with excitation (360 nm) and emission (485 nm) wavelengths.

### Notes:

Enzyme Dilution Buffer (EDB): 20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 0.5 mM EGTA, 1 mM DTT, 1 mg/ml Bovine Serum Albumin.

# Sensor Peptide $K_m$ Determination for AQT0440

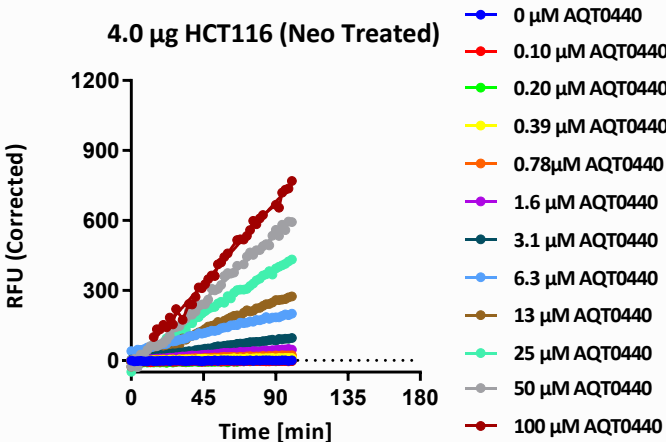
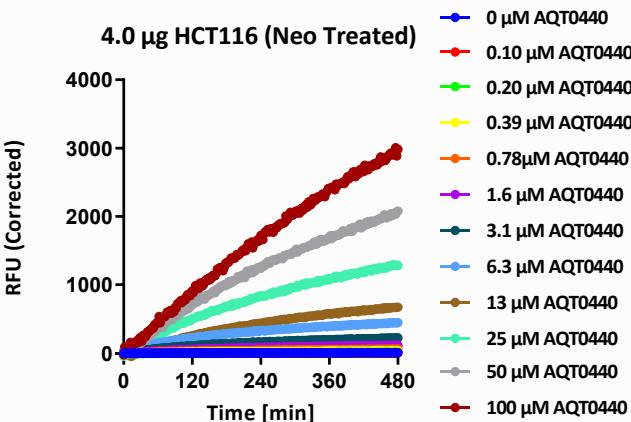


## Titration Curves and $K_m$ Plot and Table

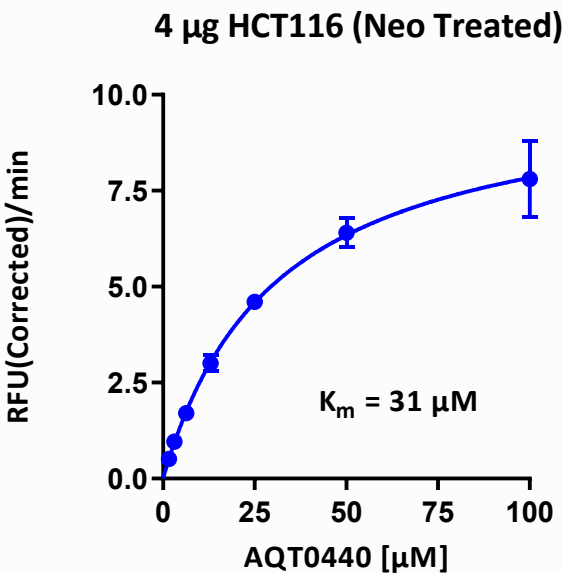
### Sensor Peptide Titration Progress Curves

Complete Time Course

Linear Region



### Sensor Peptide $K_m$ Plot



Michaelis-Menten	
Best-fit values	
Vmax	10.26
Km	30.91
95% CI (profile likelihood)	
Vmax	10.03 to 10.49
Km	29.28 to 32.63
Goodness of Fit	
Degrees of Freedom	5
R squared	0.9998
Sum of Squares	0.008537
Sy.x	0.04132
Constraints	
Km	Km > 0
Number of points	
# of X values	7
# Y values analyzed	7

The  $K_m$  value for AQT0440 is 31 µM.



# DMSO Tolerance Test with the AQT0440 Sensor Peptide

## Reaction Conditions and Set Up



### Reaction 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 calf thymus DNA

10 mM MgCl<sub>2</sub>

15 µM AQT0440 sensor peptide substrate

4.0 µg/well HCT116 crude cell lysate (from cells treated with  
0.5 µg/mL Neocarzinostatin)

**0-10% DMSO**

### Reaction Set Up:

2.5 µL 10X DMSO Titration

17.5 µL Reaction Mix with CSx Substrate, ATP & DTT

Seal plate and incubate at 30 °C for 15 minutes to equilibrate

5 µL Enzyme dilution buffer (EDB) (1x) or Lysate (5x in EDB)

25 µL Final reaction volume

Reaction was run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat-bottom polystyrene NBS microplates (Cat. #3824) at 20 or 25 µL final well volume or in in PerkinElmer, ProxiPlate-384 Plus, white shallow well microplates (Cat. #6008280) at 20 µL final well volume after sealing using optically-clear adhesive film (TopSealA-Plus plate seal, PerkinElmer [Cat. #6050185]) in a Biotek Synergy Neo 2 microplate reader with excitation (360 nm) and emission (485 nm) wavelengths.

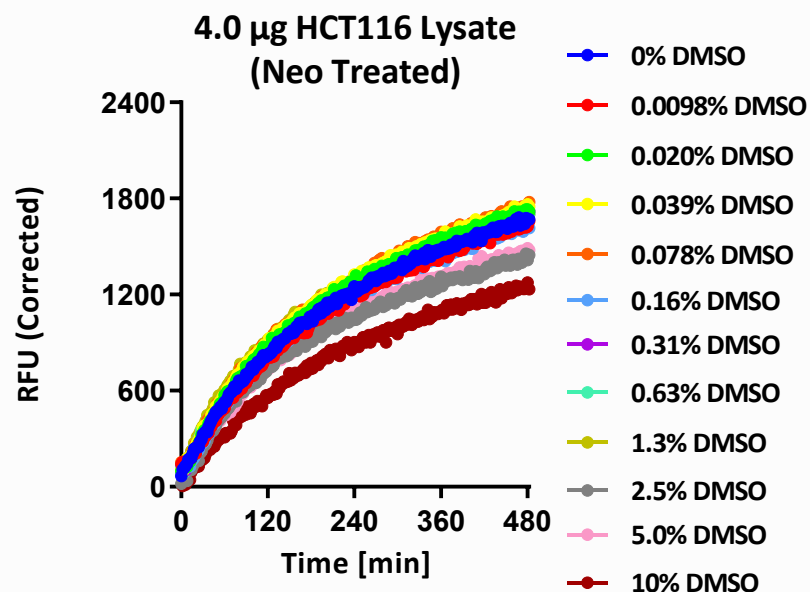
### Notes:

Enzyme Dilution Buffer (EDB): 20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 0.5 mM EGTA, 1 mM DTT, 1 mg/ml Bovine Serum Albumin.

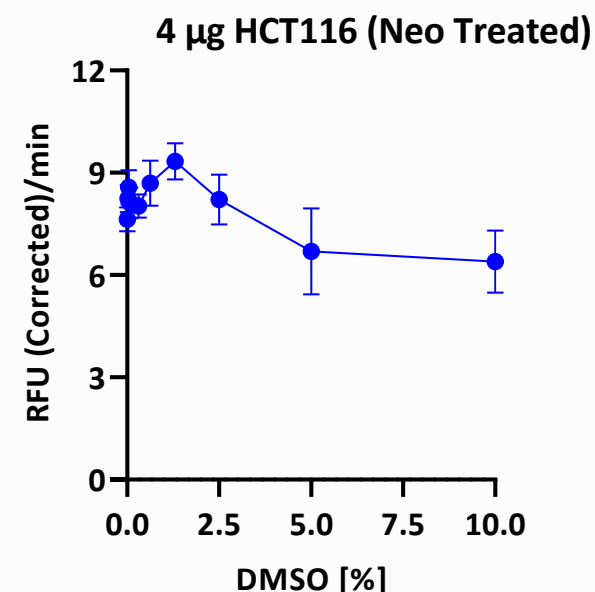
# DMSO Tolerance Test with the AQT0440 Sensor Peptide

## Titration Curves and Inhibition Plot

### Complete Progress Curves



### Reaction Rate vs [DMSO] Plot



*No significant loss in enzyme activity was observed up to 2.5% DMSO. There was a 13% loss in signal at 5% DMSO, and a 25% loss in signal at 10% DMSO. We assessed compound potency using 2% DMSO final.*

# IC<sub>50</sub> Determination with the AQT0440 Sensor Peptide



## Reaction Conditions and Set Up

### Reaction Conditions:

54 mM HEPES, pH 7.5  
1.0 mM ATP  
1.2 mM DTT  
0.012% Brij-35  
1% glycerol  
0.2 mg/ml BSA  
0.54 mM EGTA  
1 µg/mL calf thymus DNA (for recombinant DNA-PK only)  
10 mM MgCl<sub>2</sub>  
15 µM AQT0440 sensor peptide substrate

### **Compounds:**

- AZD-7648 was titrated with 3-fold dilutions in 100% DMSO at 50X the final concentrations and then diluted 50-fold into the assay for final concentrations from 0-10 µM in 2% DMSO.

### **DNA-PK Enzyme:**

- **Cell Lysate for IC<sub>50</sub> determination:**
  - 2.0 µg/well untreated HCT116 crude cell lysate
- **Recombinant DNA-PK for AQT1196 phosphorylation:**
  - 20 Units/well Full-length DNA-PK protein with the Ku domain (Promega, V5811)

### Reaction Set Up:

0.5 µL 50X AZD-648 diluted in 100% DMSO or DMSO alone  
19.5 µL Reaction Mix with CSx Substrate, ATP & DTT  
Seal plate and incubate at 30 °C for 15 minutes to equilibrate  
5 µL Enzyme dilution buffer (EDB) (1x), Lysate Buffer or DNA-PK in Lysate Buffer (5x in EDB)  
25 µL Final reaction volume

Reaction was run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat-bottom polystyrene NBS microplates (Cat. #3824) at 20 or 25 µL final well volume or in PerkinElmer, ProxiPlate-384 Plus, white shallow well microplates (Cat. #6008280) at 20 µL final well volume after sealing using optically-clear adhesive film (TopSealA-Plus plate seal, PerkinElmer [Cat. #6050185]) in a Biotek Synergy Neo 2 microplate reader with excitation (360 nm) and emission (485 nm) wavelengths.

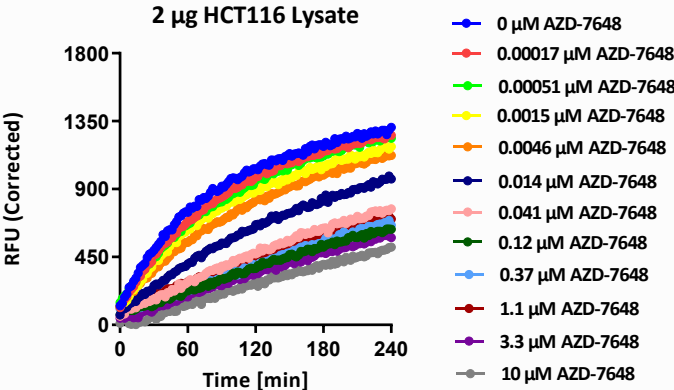
### Notes:

1X Enzyme Dilution Buffer (EDB) is used to dilute enzyme and for the blank. Composition is 20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 0.5 mM EGTA, 1 mM DTT, 1 mg/ml Bovine Serum Albumin.

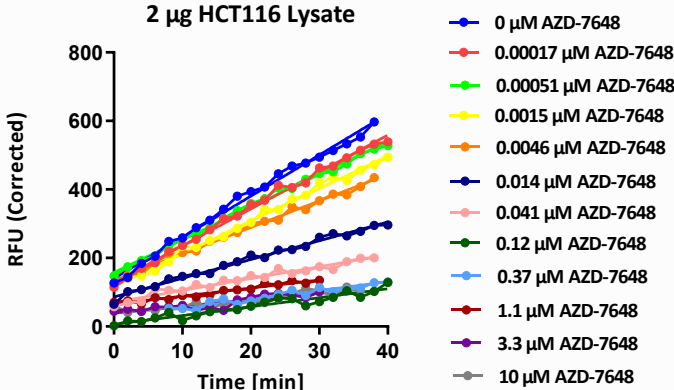
# IC<sub>50</sub> Determination with Untreated HCT116 Lysate using AZD-7648 and the AQT0440 Sensor Peptide



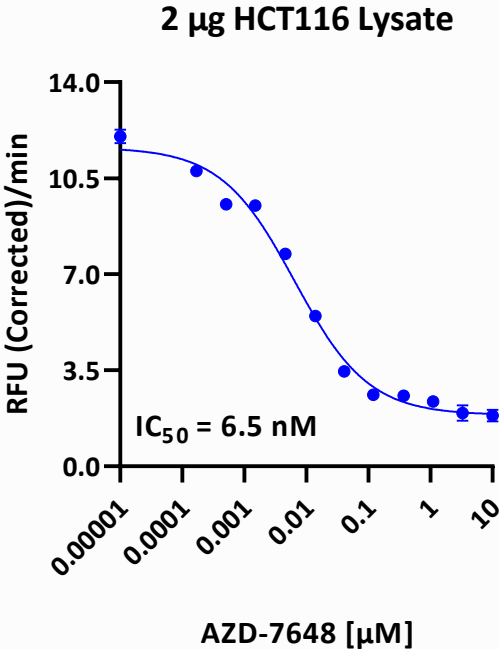
## Full Progress Curves



## Progress Curves for Linear Region



## IC<sub>50</sub> Curve



[Inhibitor] vs. response -- Variable slope (four parameters)

Best-fit values	
Bottom	1.865
Top	11.62
IC50	0.006506
HillSlope	-0.7331
logIC50	-2.187
Span	9.760
95% CI (profile likelihood)	
Bottom	1.116 to 2.455
Top	10.68 to 12.90
IC50	0.003635 to 0.01069
HillSlope	-1.105 to -0.5001
logIC50	-2.440 to -1.971
Goodness of Fit	
Degrees of Freedom	8
R squared	0.9906
Sum of Squares	1.534
Sy.x	0.4379
Constraints	
IC50	IC50 > 0
Number of points	
# of X values	12
# Y values analyzed	12

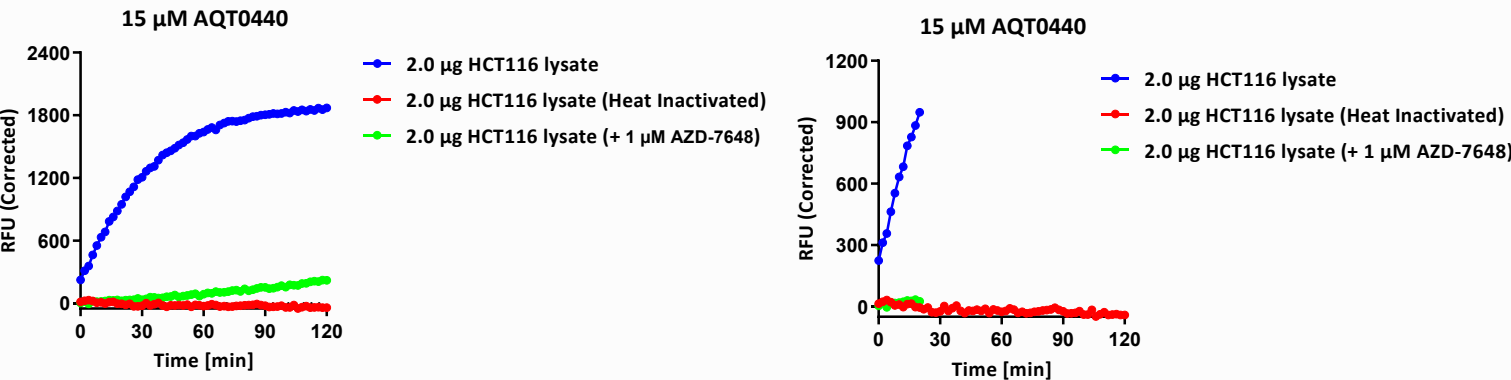
The IC<sub>50</sub> value for AZD-7648 in HCT116 Cell Lysate is 6.5 nM.



# DNA-PK Lysate Activity Assay Using the AQT0440 Sensor Peptide



## 1) Full Time Course (0–120 min.) & Linear Range (0–40 min.) with 2.0 µg/well

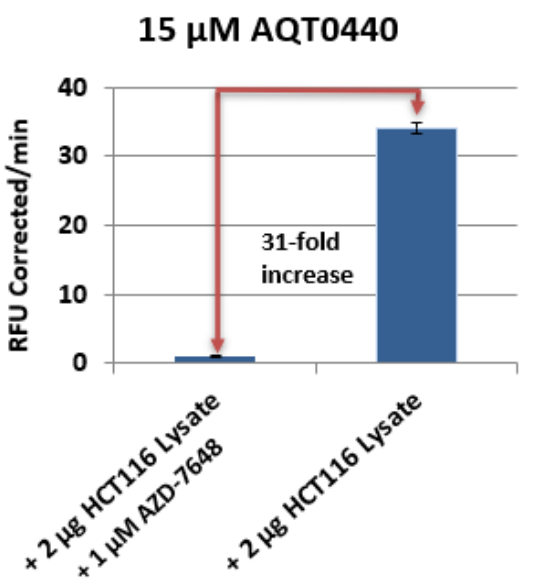


## 2) Reaction Rates to Assess Change

	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

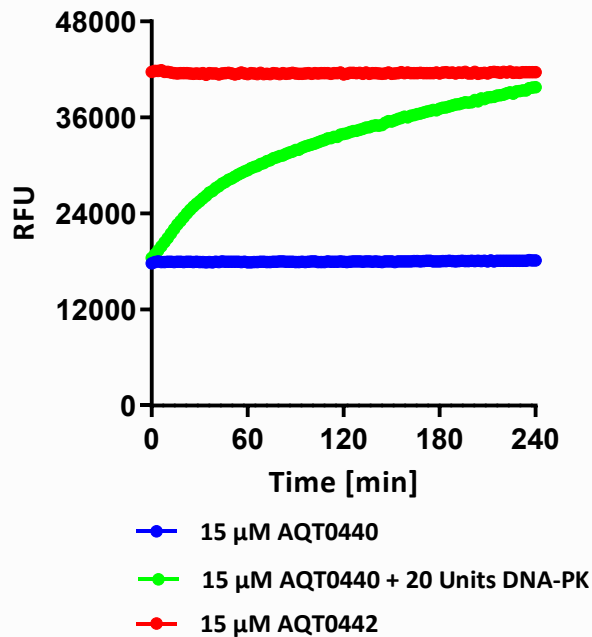
**Crude lysate samples:** The AQT0440 sensor peptide (15 µM) was used to generate **1)** Full progress curve time course and the linear range. The reaction rates (RFU Corrected values [Total – Background]/min. +/- standard deviations) are the slope of the linear region of each progress curve, which are presented in **2)**, highlighting complete elimination of signal with heat inactivation and by AZD-7648, a DNA-PK inhibitor.

**Note:** The amount of activation depends on several factors, including cell type and passage number, serum deprivation pretreatment used to make cells quiescent, and the nature, concentration, and duration of the exposure to DNA-damaging agents. With HCT116 Cell Lysates, we saw little to no effect on activity with 0.5 µg/mL Neocarzinostatin treatment for 1 hour. 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.



# DNA-PK Recombinant Activity Assay with AQT0440 Sensor Peptide and Phosphocontrol AQT0442

## Full Time Course (0–4 hours)

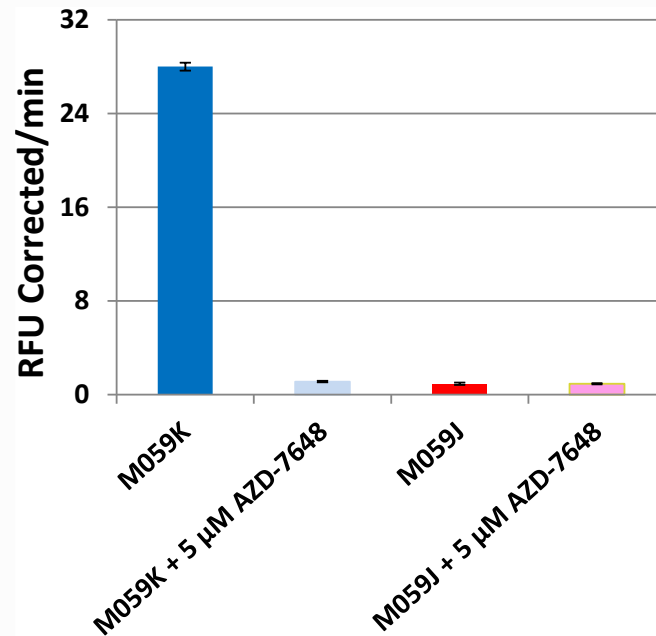


Time course for phosphorylation of the AQT0440 sensor peptide substrate (15 μM) ± DNA-PK recombinant enzyme, compared to the signal with the AQT0442 phosphopeptide control (15 μM): Full-length DNA-PK protein with the Ku domain (20 Units, Promega, V5811) fully phosphorylated the AQT0440 sensor peptide substrate by 4 hours, 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). It was necessary to add calf thymus DNA to this reaction since recombinant DNA-PK was used rather than a cell lysate. The signal with AQT0442 is used to convert RFU (Corrected) values to nmoles of phosphopeptide product/minute.

AQT0442 is a chemically synthesized phosphopeptide control for AQT0440

# DNA-PK Lysate Activity Assay Using AQT0440 with a Paired Set of M059 Glioblastoma Cell Lines

## 2 $\mu$ g M059K/M059J Lysates/well with AQT0440 Sensor Peptide

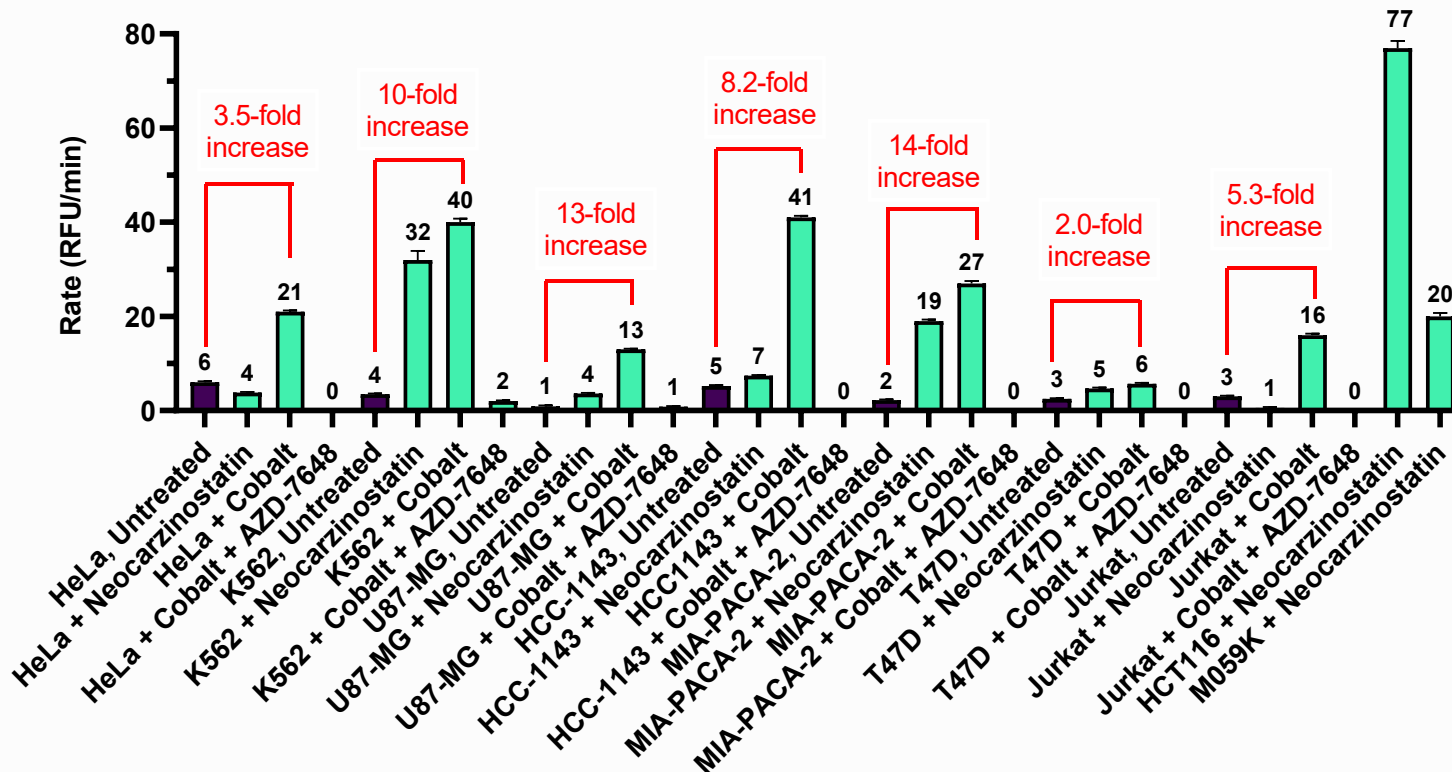


**Background:** M059K and M059J were originally isolated concurrently from the same tumor specimen from a 33-year-old male patient with untreated glioblastoma. M059K cells express normal levels of DNA-dependent protein kinase (DNA-PK), while M059J cells lack DNA-PK protein and activity. M059K cells are approximately 30-fold less sensitive to ionizing radiation than M059J cells due to the ability of DNA-PK to promote double-strand DNA break repair. M059K cells are also less sensitive than M059J cells to the cytotoxic effects of bleomycin, N, N-bis(2-chloroethyl)-N-nitrosourea and nitrogen mustard. M059K cells are DNA double-strand break repair proficient, while M059J cells are deficient in the repair of DNA double-strand breaks. The method described on slide 8 was used to prepare cell lysates for the cell lines in the histogram. Standard lysate assay conditions on slide 15 were used to run the assay. Reaction rates (RFU Corrected/min) were determined from the slopes using the linear portion of each progress curve. Values are the average of triplicate reactions  $\pm$  standard deviation. In M059K cell lysates vs M059J cell lysates, there was a 25-fold increase in the rate. This activity was inhibited by 5  $\mu$ M AZD-4648, a selective DNA-PK inhibitor.

*Demonstrates detection of endogenous DNA-PK activity in M059K cells and full inhibition with the DNA-PK inhibitor AZD-7648. As expected, no DNA-PK activity was observed with the M059J Control cell line.*

# DNA-PK Lysate Activity Assay Using the AQT0440 Sensor Peptide Across a Variety of Cell Types

## 2 µg Lysate/well with AQT0440 Sensor Peptide



### Description of cell lines where lysates had the highest DNA-PK activity (and/or highest increases with treatment):

- K562 cells** have a hematopoietic morphology and were isolated from the bone marrow of a 53-year-old female with chronic myeloid leukemia (CML) at blast crisis.
- U87-MG cells** have an epithelial morphology and were isolated from malignant gliomas from a male patient, likely with Glioblastoma.
- HCC1143 cells** are a triple negative (ER-/PR-/HER2-) mammary gland breast cancer line isolated from a 54-year-old female patient with ductal carcinoma TMN Stage IIA, grade 3.
- MIA-PaCa-2 cells** are an adherent epithelial cell line isolated from the pancreas of a 65-year-old male with carcinoma.
- HCT116 cells** are an adherent cell line isolated from the colon of a patient with colon cancer. There is a mutation in codon 13 of the ras-proto-oncogene. This cell line is near-diploid and has a relatively stable genetic profile.

**Method** - described on slide 8 was used to prepare cell lysates for the cell lines in the histogram. Standard lysate assay conditions on slide 15 were used to run the assay. Reaction rates (RFU Corrected/min) were determined from the slopes using the linear portion of each progress curve. Values are the average of duplicate reactions +/- standard error.

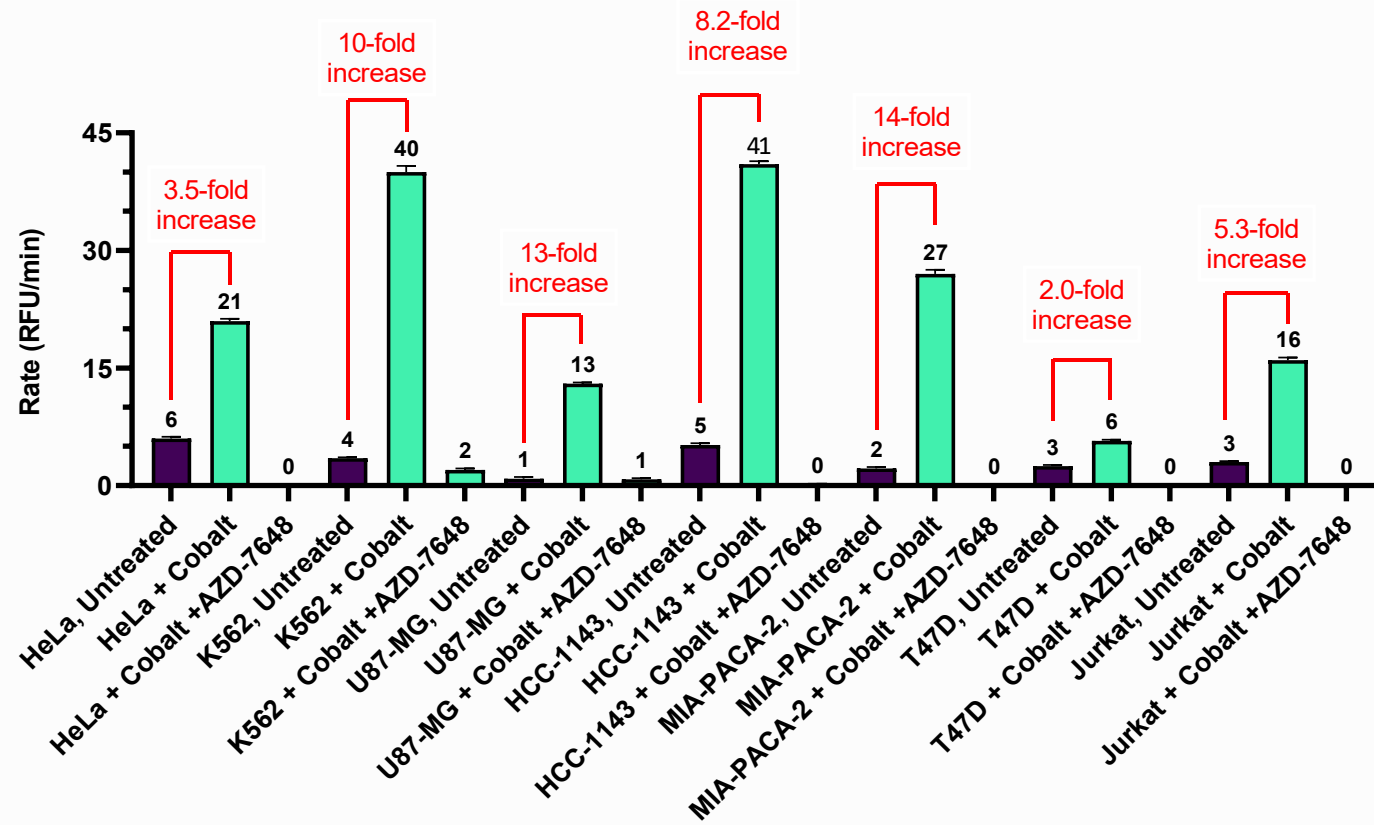
**Results** – Cobalt was more effective than Neocarzinostatin in increasing reaction rates, likely because it is more effective in causing double-strand DNA breaks. No or very low DNA-PK activity was observed with A549 and MCF-7 cells, and no significant increase in rate was observed with either treatment (data not shown).

*Demonstrates detection of high DNA-PK activity in HCT116 (Colorectal), HCC1143 (triple negative breast), K562 (CML), and MIA-PACA-2 (Pancreatic) cell lysates after treatment with Cobalt or Neocarzinostatin and inhibition with the selective DNA-PK inhibitor AZD-7648.*



# DNA-PK Lysate Activity Assay Using the AQT0440 Sensor Peptide Across a Variety of Cell Types Following Treatment with Cobalt

## 2 µg Lysate/well with AQT0440 Sensor Peptide



### Cell Lysates with Cobalt Treatment ± AZD-7648

HCC1143, K562, U87-MG, and MIA-PACA-2 cells are described on slide 20. Lysates from these cells had the largest increase in DNA-PK activity with Cobalt treatment vs the untreated controls with increases in rate of 8-fold for HCC1143, 10-fold for K562, 13-fold for U87-MG, and 14-fold for MIA-PACA-2. AZD-7648 inhibited all DNA-PK activity in these lysates for 30 minutes. Cobalt treatment results in an increase in reactive oxygen species, which can cause DNA double-strand breaks that activates DNA-PK catalytic kinase activity that is measured selectively with the AQT0440 sensor peptide.

*Demonstrates activation of DNA-PK activity in all cells (2-14-fold) by cobalt treatment with inhibition by the selective DNA-PK inhibitor AZD-7648.*

# Summary

- ❖ The PhosphoSens-Lysate Assay for DNA-PK using the AQT0440 selective sensor peptide demonstrates a robust, sensitive, and physiologically relevant assay that selectively measures endogenous DNA-PK activity with all the cellular components and signaling complexes using only 2 µg of crude cell lysate per test.
- ❖ Results include:
  - Treatment with Neocarzinostatin or Cobalt resulted in up to 10-fold and 14-fold increase in DNA-PK activity, respectively, with activation observed across multiple cell lines. Both chemicals result in DNA double-strand breaks. The highest DNA-PK activity was with HCT116 cells treated with Neocarzinostatin, resulting in a rate that was almost 2-fold higher than any other cell line. This lysate was used for further validation studies.
  - DNA-PK activity with lysates from Neocarzinostatin-treated HCT116 cells was linear from 0.078 to 5.0 µg/well, a 64-fold linear range.
  - The DNA-PK lysate activity was inhibited by 97% with 1 µM of AZD-7648 DNA-PK inhibitor reference compound.
  - The IC<sub>50</sub> value for AZD-7648 with lysate from untreated HCT116 cells was 6.5 nM.
  - A K<sub>m</sub> of 31 µM was determined for AQT0440 with the Neocarzinostatin-treated HCT116 cell lysate.
- ❖ These DNA-PK activity measurements are direct, highly quantitative, and in an easy-to-use format. This allows functional assessment of the native DNA-PK/Ku activity in complex samples using crude cell lysates or tissue homogenates, thereby providing a more physiological and economical approach to study DNA-PK.

# Company & Technology Supporting Slides

# AQT Senior Scientific Team

*Experienced & Diverse Team from Leading Institutions*



**Dr. Bill Radany**  
CEO & Head of  
Business Development



**Dr. Erik Schaefer**  
Co-Founder, CSO



**Dr. Barbara Imperiali**  
Co-Founder &  
CTO, MIT Faculty



**Dr. Earl May**  
Senior Director, Discovery  
Technologies (DT)



**Dr. Eric Berg**  
Director of Peptide  
Operations



**Dr. Daniel Urul**  
Senior Scientist, DT



**Dr. Venky Nemmara**  
Senior Scientist, DT

- ❖ We apply our >200 years of collective experience in kinase biology, drug or assay development, chemical-biology and automation to solve critical problems in target biology and drug development.
- ❖ AQT services allow us to evaluate drugs in diverse ways, delivering high-quality data combined with strong communication to enable discovery



**Susan Cornell-Kennon**  
Senior Scientist, DT

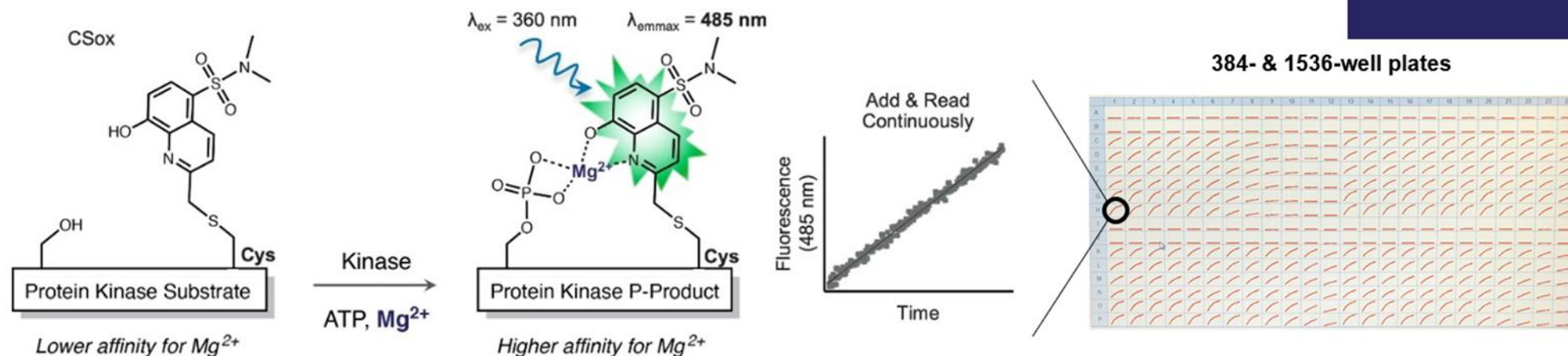


**Dr. Satish Pimrale**  
Business  
Development Leader





# Sensor Peptide Substrates for Continuous (Kinetic) Monitoring of Protein Kinase Activity



- ❖ Uses chelation-enhanced fluorescence via sulfonamido-oxine (Sox) chromophore, invented by Barbara Imperiali (MIT).
- ❖ 10 patents with 1 pending, exclusively licensed from MIT. Rich trade secret portfolio
- ❖ Sox is small, minimally hydrophobic, and neutral
- ❖ Assay is direct, homogeneous, and kinetic (continuous)
- ❖ Can use a wide range of ATP concentrations (Low  $\mu\text{M}$ , ATP  $K_m$  or mM [physiological])
- ❖ Runs on commonly available readers with any plate type (96-, 384- or 1536-well)

*Simple, Powerful & Flexible!*

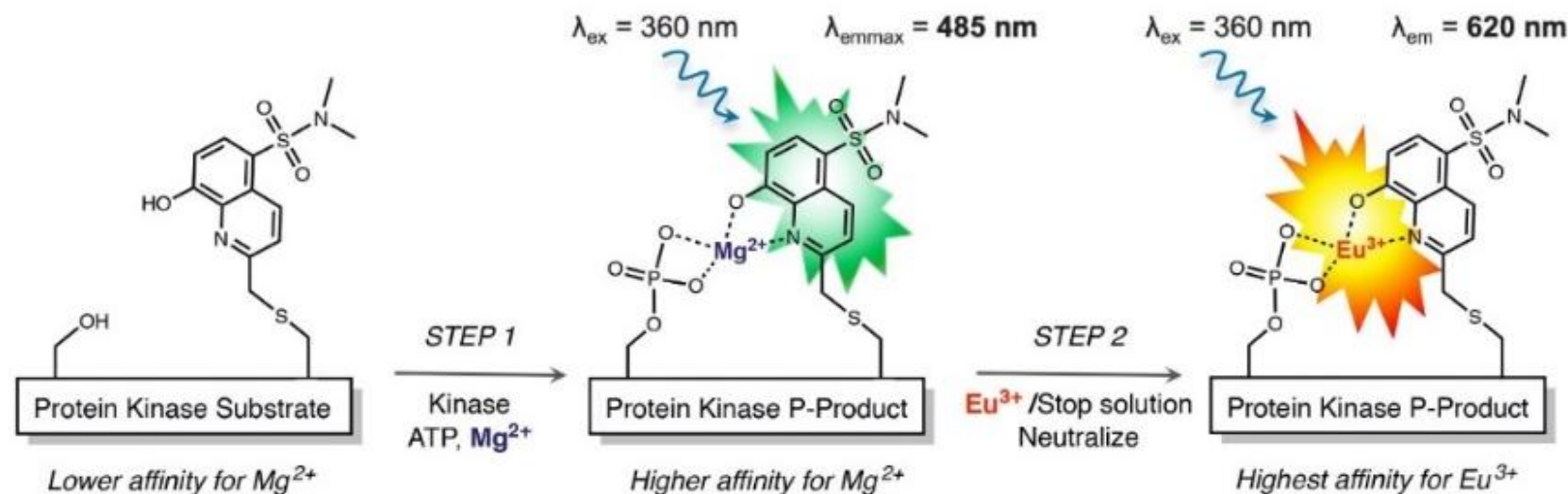
*a full progress  
curve in every well*

**Ideal for quantitative  
assessments including  
Profiling, Potency & MOA**

**PhosphoSens<sup>®</sup>**



# Replacing $\text{Mg}^{2+}$ with $\text{Eu}^{3+}$ Create **PhosphoSens<sup>®</sup> - Red**



**Ideal for:**

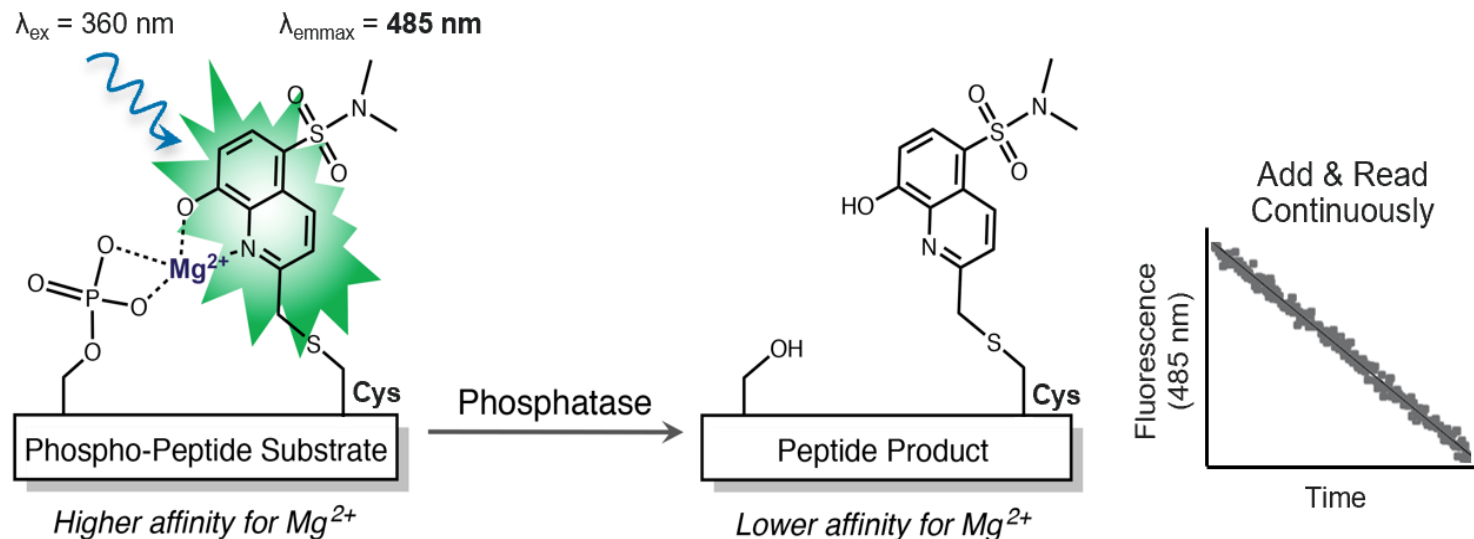
- **HTS**
- **SAR**

- ❖ Sox sensitizes Europium ( $\text{Eu}^{3+}$ ) luminescence for time-resolved fluorescence (TRF) with 620 nm emission and data acquisition delay, which eliminates compound interference
- ❖ Run as an Endpoint format; ideal for high-numbers of tests for HTS and SAR
- ❖ Patent approved in 2020

*Eliminates autofluorescence while using the same small fluorophore*

# Same Sox-based Detection Platform Enables Continuous Monitoring of **Protein**

**Pho**



- Same mechanism - ChEF ( $\text{Mg}^{2+}$ ) for kinetic assays and  $\text{Eu}^{3+}$ /TRF for Endpoint (Red), as for protein kinases
- Uses **CSox-based phosphopeptide substrates** derived from physiological targets to monitor loss of signal
- Covered by 10 patents (with 1 pending) and trade secret portfolio

*Increasing requests for tyrosine & serine/threonine phosphatases*

Kinetic and Endpoint

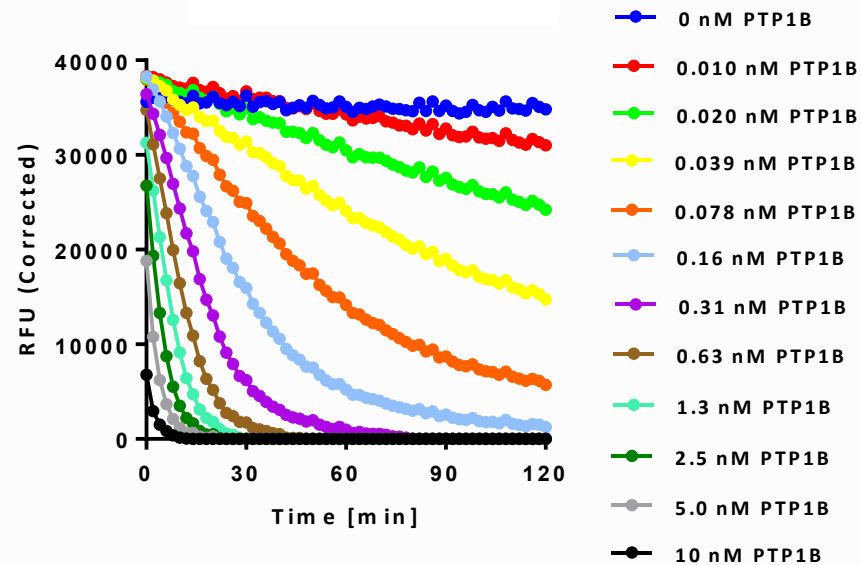
*One Technology,  
Two Fit-for-Purpose  
Formats*

Expanding menu  
(currently 34 of  
the  
190 Protein  
Phosphatases)

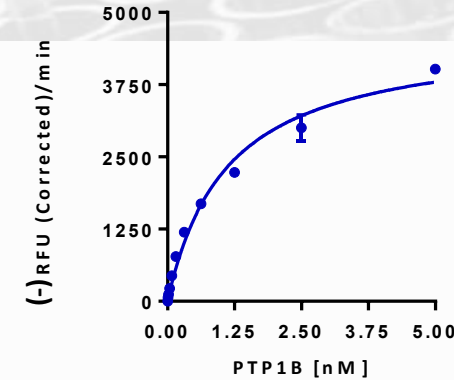
# Analysis of PTP1B & PTPN2 Tyrosine Phosphatases

## Optimized *PhosphoSens* Phosphopeptide Substrates

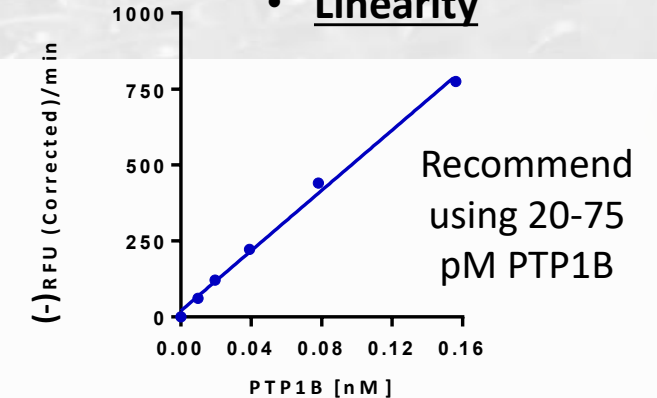
### Progress Curves & Linearity for PTP1B with AQT0266 Sensor Peptide, Net Signal



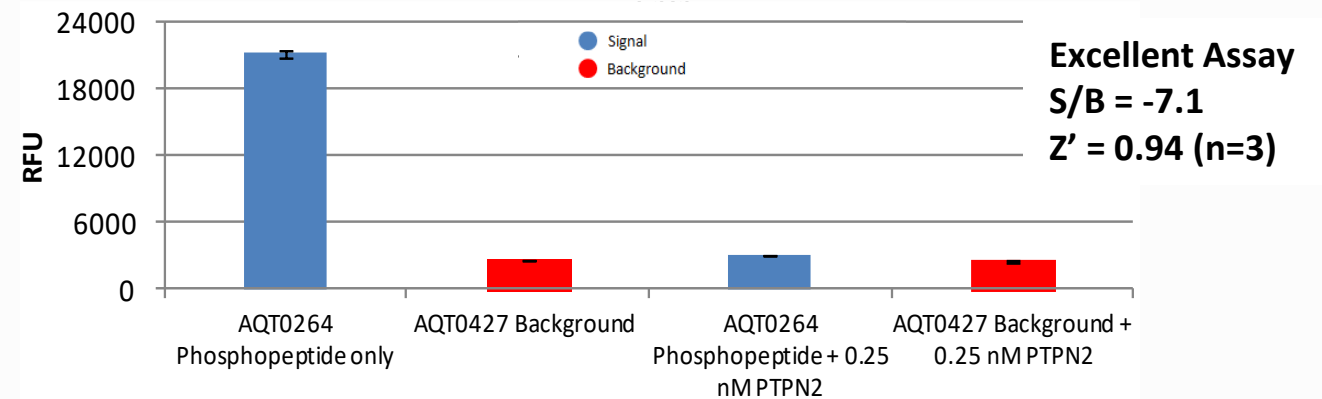
### Total Range



### Linearity



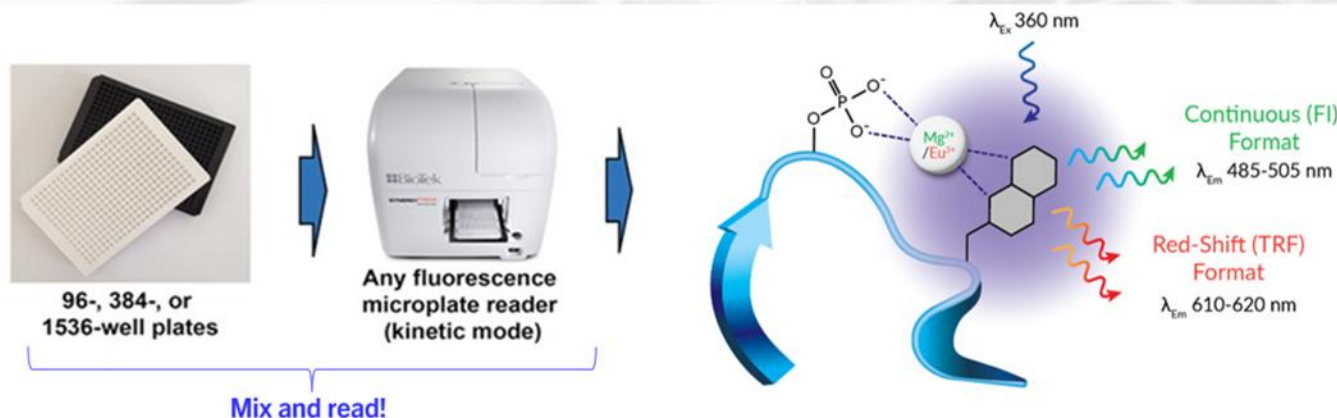
### Total & Background for PTPN2 with AQT0264 in PhosphoSens®-Red



*High-sensitivity & physiologically-relevant. Analysis of allosteric or substrate-competitive inhibitors*

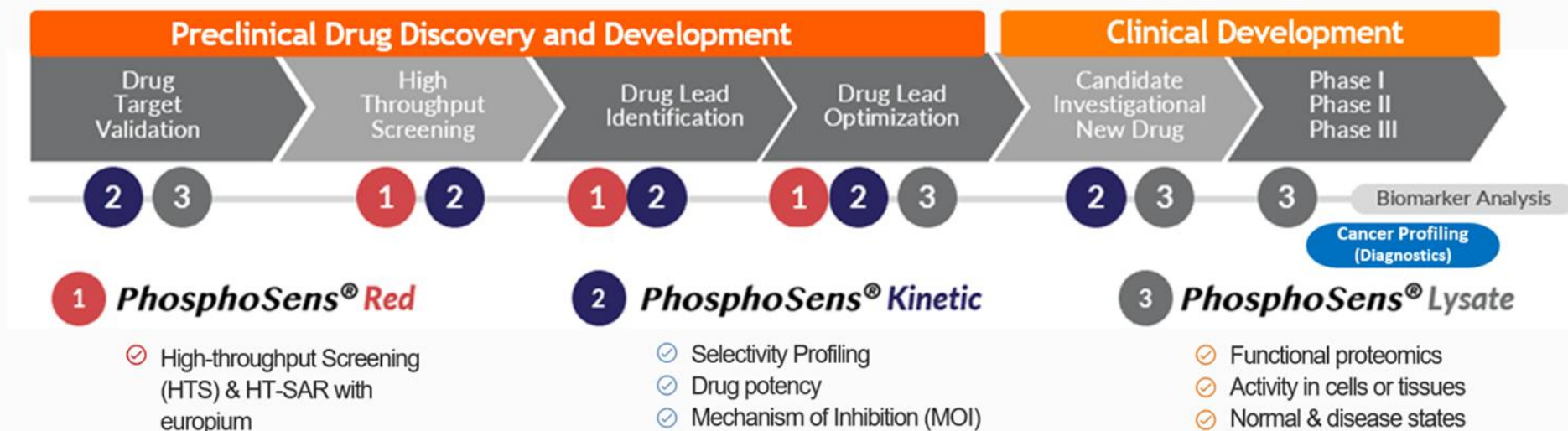
# PhosphoSens® Platform

Enabling Analysis Across the Drug Development Workflow



**One Technology,  
Two Fit-for-  
Purpose Formats**

**Kinetic and Endpoint**

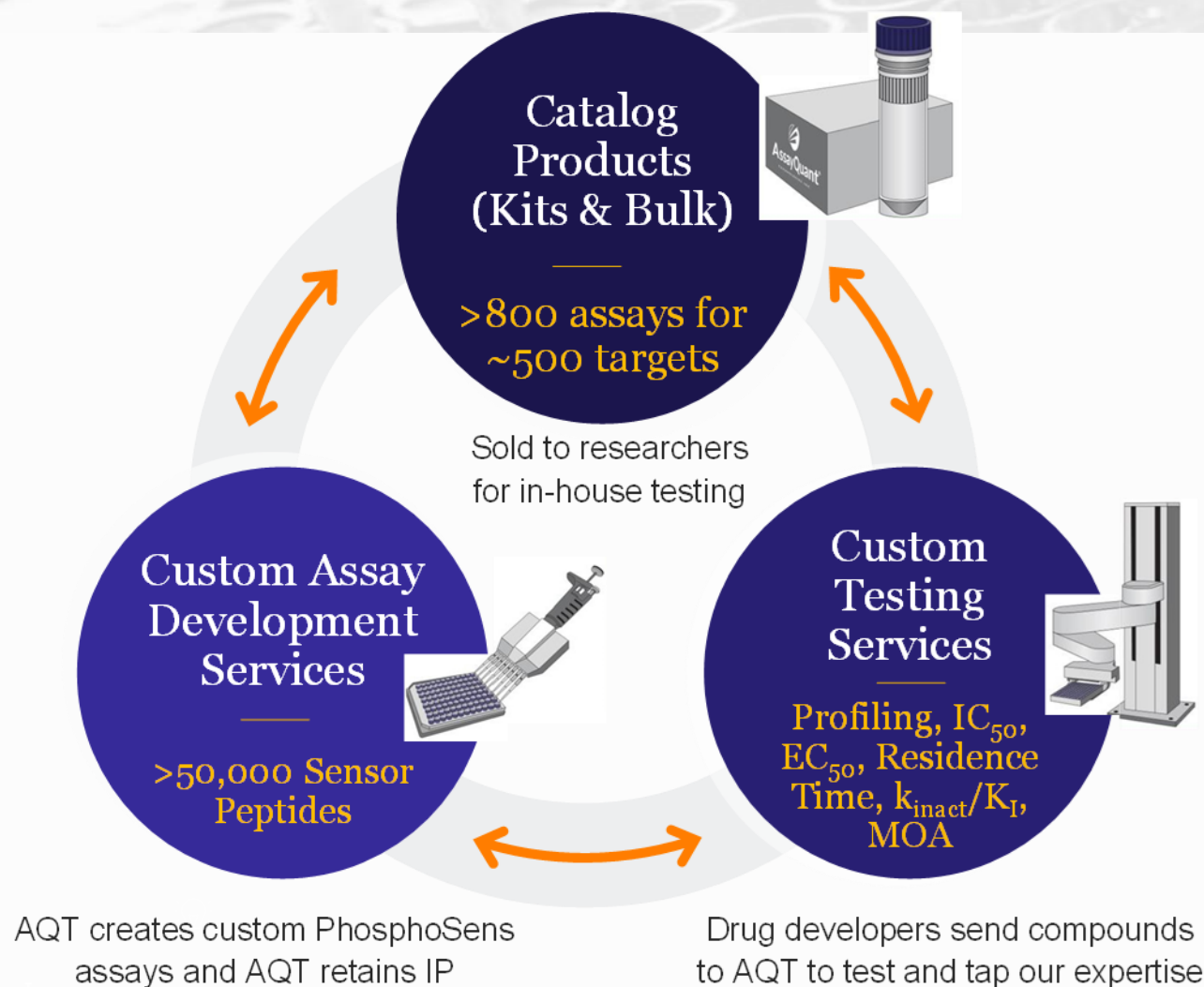


*Accelerates progress and improves outcomes*



# Integration of Catalog Products & Services

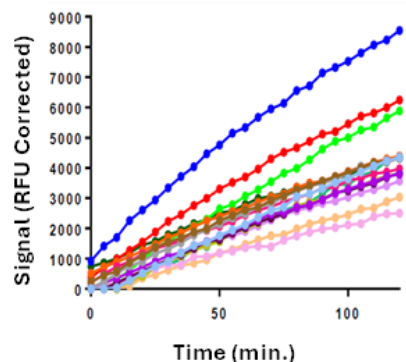
*Multiple entry & transition points to address your needs*



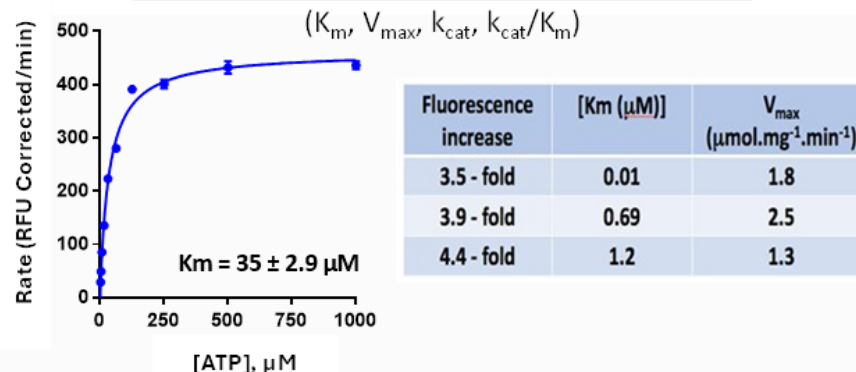


# PhosphoSens® Platform Continuous Workflow Applications

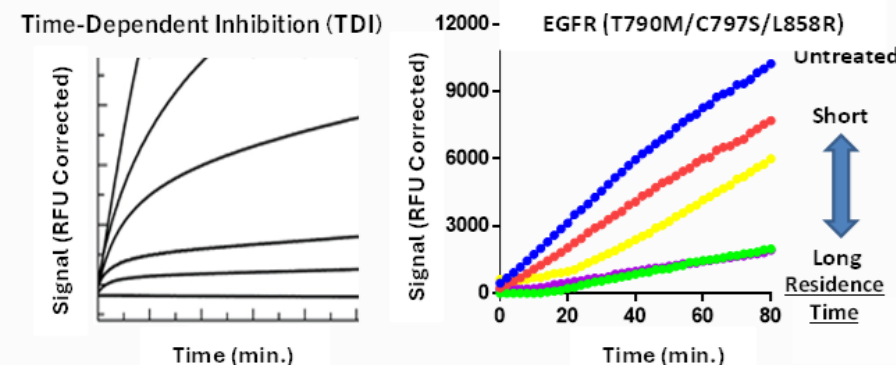
## Assay Development



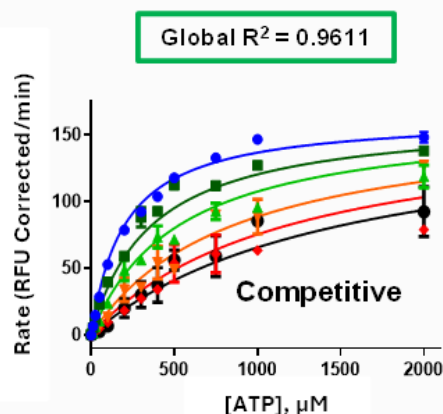
## Enzyme Kinetics with Recombinant Proteins



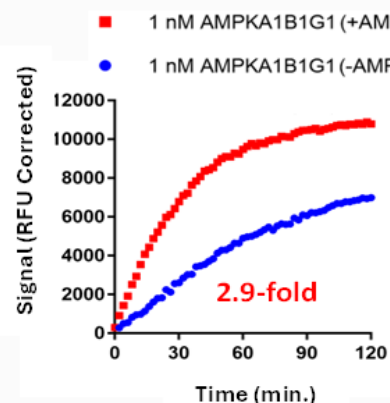
## Drug Potency ( $IC_{50}$ , $K_i$ , Residence Time, $k_{inact}/K_i$ , MOA, $EC_{50}$ )



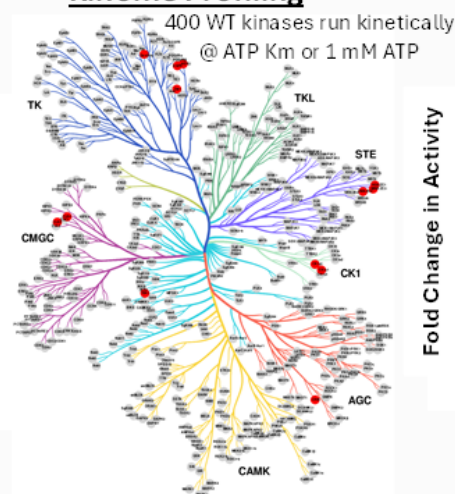
## Mechanism of Inhibition (MOI)



## Target Activation

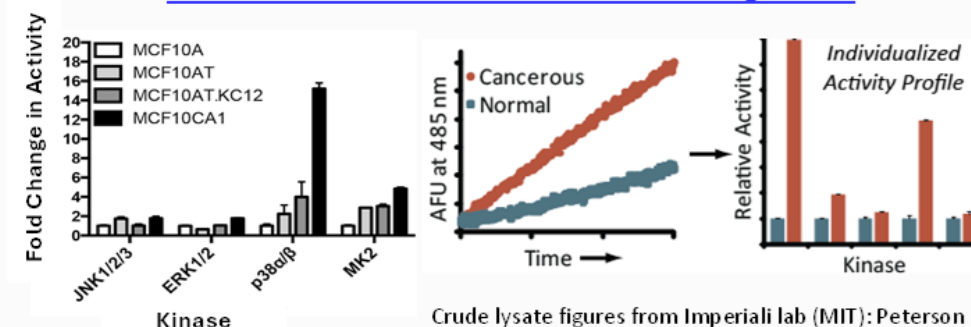


## Kinome Profiling



## Quantitative Enzyme Activity Measurements in Unfractionated Cell or Tissue Lysates

Direct-to-Phase II SBIR from NIH/NCI funded in August 2024



Crude lysate figures from Imperiali lab (MIT): Peterson *et al.* Biochemistry 2014, 53: 5771-5778; Stains *et al.* Chem Biol. 2012, 19(2): 210-217.

# PhosphoSens® Platform Advantages:

- ❖ **Sensitive** and **continuous (kinetic)** format based on Chelation-enhanced fluorescence.
- ❖ **Homogeneous (add & read)** and **direct** measure of enzyme activity resulting in a complete **progress curve in every well** for each condition.
- ❖ Sensor peptides or lipids based on **physiologically-relevant** substrates.
- ❖ Run under **optimal enzyme conditions**, with physiological  $Mg^{2+}$ ,  $Mn^{2+}$  &  $Ca^{2+}$  ions and **any ATP** concentration (1-2 mM = physiological; or at ATP  $K_m$ ) or sample type (lysates via selective sensor peptides or IP-kinase assays).
- ❖ Determination of **initial reaction rate** from **linear** portion of curve provides **high accuracy & precision** ( $Z' > 0.7$ ) and enables assessment of time-dependent inhibitor or activator potency ( $IC_{50}$ ,  $EC_{50}$ ,  $k_{inact}/K_i$ , residence time).
- ❖ **Corrects for any compound autofluorescence** – **Kinetic**: Background signal doesn't change over time; **Red**: Uses Europium ( $Eu^{3+}$ ) for time-resolved fluorescence (TRF) with a 100  $\mu$ sec delay & 620 nm emission.
- ❖ **Rapid & predictable development** of new assays is achieved even with difficult targets using >30,000 sensor peptide panel for protein kinases or phosphatases and any DAG species for DGK assays.
- ❖ **Rigorous manufacturing** process for PhosphoSens® sensors provides excellent lot to lot consistency.
- ❖ **Works across entire drug development workflow** – Improving efficiency and performance

*We apply our >200 years of collective experience in kinase biology, drug or assay development, and automation to address complex biology and evaluate drugs in diverse ways, delivering high-quality data to accelerate your programs.*