

# PhosphoSens GSK3A/B Cell Lysate Kinase Assay Kit Protocol & Validation Data

---

Determination of GSK Activity in Crude Lysates from A375 Cells  
Using the Selective AQT1211 Sensor Peptide Substrate

HGNC Name: GSK3alpha and GSK3beta

Long Names: Glycogen synthase kinase-3 alpha & Glycogen synthase kinase-3 beta

# PhosphoSens GSK3A/B Cell Lysate Kinase Assay Kit (AQT1211-KL-100)

## MATERIALS INCLUDED

TABLE 1		
Components	Description	100 Assay Kit Volumes
<b>PhosphoSens Kinase-Selective Lysate Substrate, AQT1211, 1mM</b>	GSK3A/B-selective sensor peptide substrate for assaying kinase activity in complex biological samples	45 µL
<b>ATP Solution, 100 mM</b>	100 mM ATP in nuclease-free water	30 µL
<b>DTT Solution, 1 M</b>	1 M DTT in nuclease-free water	5 µL
<b>Enzyme Reaction Buffer (ERB), 10X</b>	500 mM HEPES, pH 7.5, 0.1% Brij-35, 100 mM MgCl <sub>2</sub>	300 µL
<b>Enzyme Dilution Buffer (EDB), 5X Base</b>	20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 1 mg/mL Bovine Serum Albumin (BSA)	800 µL
<b>EGTA Solution, 550 mM</b>	550 mM EGTA in water (pH adjusted with NaOH)	30 µL
<b>PhosphoPreserve Cell Extraction Buffer, 1X Base</b>	Please reach out to support@assayquant.com to inquire	1,000 µL
<b>PhosphoPreserve Protease Inhibitor Cocktail</b>	Please reach out to support@assayquant.com to inquire	20 µL
<b>PhosphoPreserve Phosphatase Inhibitor Cocktail</b>	Please reach out to support@assayquant.com to inquire	20 µL
<b>PhosphoSens PhosphoControl Peptide, AQT1158, 1 mM</b>	Fully phosphorylated version of the GSK3A/B-selective sensor peptide substrate, AQT1211.	10 µL
<b>A375 Lysate Control*</b>	Crude lysate of A375 cells in PhosphoPreserve CEB, supplemented with DTT, and PhosphoPreserve Protease and Phosphatase inhibitor cocktails.	25 µL
<b>GSK3B (Full Length, HIS Tag)*</b>	Recombinant Human GSK3B Protein (Full Length, HIS Tag), Active stored in 50 mM Tris-HCl, pH 7.5, 50-300mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25-1.0mM DTT, 0-0.1mM PMSF, 10-25% glycerol	2 µL
<b>LY2090314, 10 mM</b>	LY2090314 resuspended in 100% DMSO	5 µL

## MATERIALS NOT INCLUDED

1. Ultra-pure deionized water
2. Dimethyl Sulfoxide (DMSO), (CH<sub>3</sub>)<sub>2</sub>SO
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.

# 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

10 mM MgCl<sub>2</sub>

15 μM Sensor Peptide substrate, AQT1211 or Phosphopeptide control, AQT1213

## Kinase Enzyme:

- 0.5 nM GSK3b Recombinant Enzyme
- **For simple lysate assay** – 1.1 μg/well A375 lysates
- **\*For Lysate titration** – 0, 39, 78, 156, 313, 625, 1250, 2500 ng/well A375 lysates.  
Samples are prepared by serial dilution of the crude lysate.

***\*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 2500 ng/well. This should be determined empirically and will vary depending on the cell line and the treatment conditions.*

## Notes:

1. Total protein concentration for cell lysates was determined using a modified Bradford assay (Cat # 5000006, BIO-RAD).
2. **PhosphoPreserve Cell Extraction Buffer:** Should be **supplemented with included protease and phosphatase inhibitors** as prepared on slides 5 and 11 **just before use**.
3. **Final 1X Enzyme Dilution Buffer (EDB):** 20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 1 mg/ml Bovine Serum Albumin (BSA), and 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).

# Reagent Preparation for a Low Volume 384-well Format (25 $\mu$ 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  $\mu$ 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  $\mu$ L of 10 mM ATP by adding 9  $\mu$ L of 100 mM ATP to 81  $\mu$ L of ultrapure deionized water.
  - 10 mM DTT**: Make 300  $\mu$ L of 10 mM DTT by adding 3  $\mu$ L of 1M (1000 mM) DTT to 297  $\mu$ L of ultrapure deionized water.
  - 5.5 mM EGTA**: Make 300  $\mu$ L of 5.5 mM EGTA by adding 3  $\mu$ L of 550 mM EGTA to 297  $\mu$ L of ultrapure deionized water.
  - 150  $\mu$ M Sensor Peptide**: Make 90  $\mu$ L of 150  $\mu$ M [AQT1211](#) by adding 13.5  $\mu$ L of 1 mM [AQT1211](#) to 76.5  $\mu$ L of ultrapure deionized water. Make 14  $\mu$ L of 150  $\mu$ M [AQT1213](#) by adding 2.1  $\mu$ L of 1 mM [AQT1213](#) to 11.9  $\mu$ L of ultrapure deionized water.
  - 50  $\mu$ M GSK3A/B Inhibitor LY2090314 (50X)**: Make 100  $\mu$ L of 50  $\mu$ M LY2090314 by adding 0.5  $\mu$ L of 10 mM stock into 99.5  $\mu$ L DMSO.
- Prepare final 1X *PhosphoPreserve* Cell Extraction Buffer (CEB-B) by adding 1  $\mu$ L of 1M DTT, 16.7  $\mu$ L of the provided AQT Inhibitor Cocktail for Proteases (60X) and 16.7  $\mu$ L of the provided AQT Inhibitor Cocktail for Phosphatases (60X) to 1000  $\mu$ L of *PhosphoPreserve* Cell Extraction Buffer (CEB-B), 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  $\mu$ 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  $\mu$ 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 AQT1211

Components for 1.28X Master Mix:	For 1 Well:	For 32 Wells:
Enzyme Reaction Buffer (10X)	2.5 $\mu$ L	80 $\mu$ L
ATP (10 mM)	2.5 $\mu$ L	80 $\mu$ L
DTT solution (10 mM)	2.5 $\mu$ L	80 $\mu$ L
EGTA Solution (5.5 mM)	2.5 $\mu$ L	80 $\mu$ L
Selective Sensor Peptide Substrate <a href="#">AQT1211</a> (150 $\mu$ M)	2.5 $\mu$ L	80 $\mu$ L
Ultrapure deionized water	7.0 $\mu$ L	224 $\mu$ L
<b>Total volume</b>	<b>19.5 <math>\mu</math>L</b>	<b>624 <math>\mu</math>L</b>

## Table 3 - Phospho-Peptide Control AQT1213

Components for 1.28X Master Mix:	For 1 Well:	For 5 Wells:
Enzyme Reaction Buffer (10X)	2.5 $\mu$ L	12.5 $\mu$ L
ATP (10 mM)	2.5 $\mu$ L	12.5 $\mu$ L
DTT solution (10 mM)	2.5 $\mu$ L	12.5 $\mu$ L
EGTA Solution (5.5 mM)	2.5 $\mu$ L	12.5 $\mu$ L
Sensor Phosphopeptide Control <a href="#">AQT1213</a> (150 $\mu$ M)	2.5 $\mu$ L	12.5 $\mu$ L
Ultrapure deionized water	7.0 $\mu$ L	35.0 $\mu$ L
<b>Total volume</b>	<b>19.5 <math>\mu</math>L</b>	<b>97.5 <math>\mu</math>L</b>

# 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  $\mu\text{L}$  of 1X EDB, add 1000  $\mu\text{L}$  5X EDB Base along with 5  $\mu\text{L}$  of 1M DTT, 5  $\mu\text{L}$  of 550 mM EGTA, and 3990  $\mu\text{L}$  of ultrapure deionized water to create the final composition shown on slide 5. **Keep on ice.**
2. Prepare 250  $\mu\text{L}$  BLANK by combining 249.571  $\mu\text{L}$  1X EDB and 0.429  $\mu\text{L}$  of 1X Final PhosphoPreserve Cell Extraction Buffer .
3. Prepare 250  $\mu\text{L}$  of 2.5 nM recombinant GSK3beta (5X) by adding 0.429  $\mu\text{L}$  of the 1458 nM GSK3beta stock to 249.571  $\mu\text{L}$  of 1X EDB. **Keep on ice until needed.**
4. Prepare 35  $\mu\text{L}$  of 0.22 mg/mL lysate (5X) by adding 2.14  $\mu\text{L}$  of the 3.6 mg/mL untreated A375 lysate stock to 32.86  $\mu\text{L}$  of 1X EDB. **Keep on ice until needed.**
5. Transfer 20  $\mu\text{L}$  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  $\mu\text{L}$  of 50  $\mu\text{M}$  GSK3A/B Inhibitor to all wells in columns 2, 5 and 7 (as shown on the plate map on the next slide). Add 0.5  $\mu\text{L}$  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  $\mu\text{L}$  of 1.28X Master Mix for AQT1211 from Table 2 to all the wells except the wells labeled 'AQT1213'. For the wells labeled 'AQT1213', add 19.5  $\mu\text{L}$  of 1.28X Master Mix for AQT1213 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 AQT1213 sensor phosphopeptide positive control):
  - Add 5  $\mu\text{L}$  BLANK (prepared in step 2 above) to the wells labeled 'BLANK' or 1X EDB to wells labeled '1X EDB'
  - Add 5  $\mu\text{L}$  1X EDB to wells labeled 'AQT1213'
  - Add 5  $\mu\text{L}$  of the A375 lysate to the corresponding wells labeled '1.1  $\mu\text{g}$  Lysate'
  - Add 5  $\mu\text{L}$  of 2.5 nM GSK3A/B (5X) to the wells labeled '0.5 nM GSK3beta'.
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.

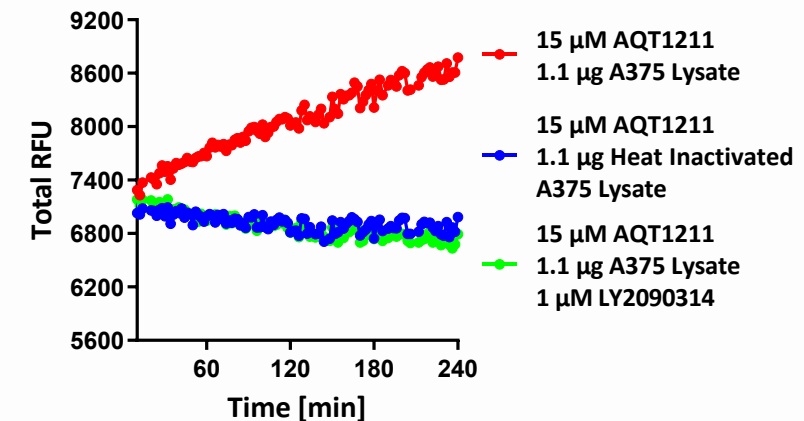
# Plate Additions and Plate Map For a Simple Lysate Activity Assay

Component	Volume to add to wells
Tool Compound or 100% DMSO	0.5 µL
Master mix	19.5 µL
Lysate, GSK3beta, BLANK, or 1X EDB	5 µL

*Total volume 25 µL*

	Plate map for Lysate activity assay							
	A375 Cell Lysate			With recombinant enzyme	EMPTY			
	No tool compound	With tool compound (LY2090314)	Heat inactivated (HI) lysate	No tool compound	With tool compound (LY2090314)	No tool compound	With tool compound (LY2090314)	Phosphopeptide control
	1	2	3	4	5	6	7	8
A	BLANK	BLANK	BLANK	1X EDB	1X EDB	EMPTY	EMPTY	AQT1213
B	BLANK	BLANK	BLANK	1X EDB	1X EDB	EMPTY	EMPTY	AQT1213
C	1.1 ug Lysate	1.1 ug Lysate	1.1 ug HI lysate	0.5 nM GSK3beta	0.5 nM GSK3beta	EMPTY	EMPTY	AQT1213
D	1.1 ug Lysate	1.1 ug Lysate	1.1 ug HI lysate	0.5 nM GSK3beta	0.5 nM GSK3beta	EMPTY	EMPTY	AQT1213

## Progress Curves for Total Fluorescence with AQT1211



It is unclear why there is a slight downward slope with the Heat-inactivated and LY2090314 compound, but these rates are close to zero.

AQT1213 data can be found on slide 10

**This test uses 32 wells (28 wells for the AQT1211 sensor peptide substrate and 4 wells for the AQT1213 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.



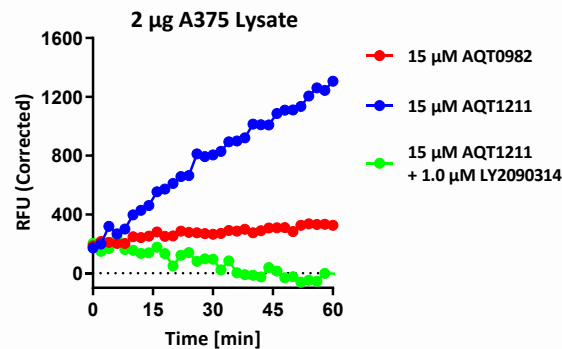
# Data for a Simple Activity Assay Validation

A375 Cell Lysate or Recombinant GSK3beta with AQT1211 Sensor Peptide Substrate or AQT1213 Control

## A. Crude Lysate Samples (2.0 µg/well)

### PROGRESS CURVES

#### 1) Linear Time Course (0–60 min)

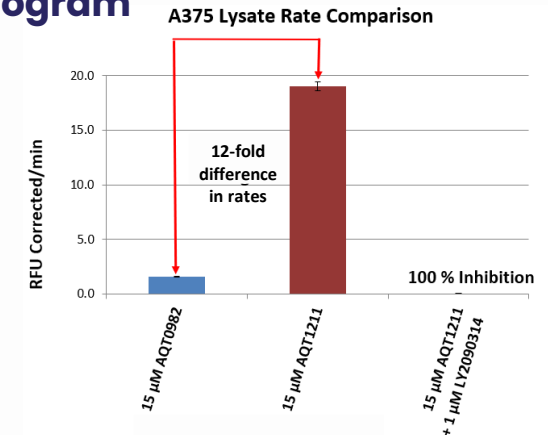


### QUANTITATIVE ASSESSMENT

#### 2) Reaction Rates to Assess Change

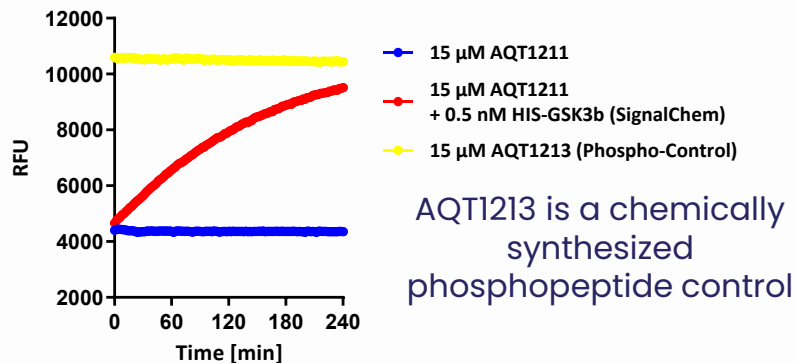
	Reaction Rate (RFU/min)	Change
2.0 µg A375, 15 µM AQT0982 (AKT1/2/3 Substrate)	$1.6 \pm 0.045$	
2.0 µg A375, 15 µM AQT1211 (GSK3A/B Substrate)	$19 \pm 0.41$	12-fold difference in rates
2.0 µg A375, 15 µM AQT1211 (GSK3A/B Substrate) + 1 µM LY2090314 (GSK3A/B Inhibitor)	$0 \pm 0.074$	100% inhibition of GSK3A/B Activity

#### 3) Histogram



## B. Recombinant GSK3beta & AQT1213 Control

#### 1) Full Time Course (0–240 min.)



**A. A375 Crude Lysate Sample:** The AQT1211 and AQT0982 sensor peptides were used to generate **1)** A linear progress curve time course (0–60 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 12-fold difference in rates comparing the activity for AKT1/2/3 with the AQT0982 substrate to the activity for GSK3A/B with the AQT1211 substrate since phosphorylation of GSK3 by AKT is inhibitory. The GSK3A/B signal was eliminated by adding the selective GSK3A/B inhibitor LY2090314.

**Note:** It is unclear why there is a slight downward slope with the LY2090314 compound (this also occurs with heat-inactivated lysate), but these rates are close to zero. The amount of activation depends on several factors, including cell type, the serum deprivation pretreatment used to make cells quiescent, as well as the nature, concentration, and duration of the activating stimulus, if applicable. These conditions can be varied to determine the effect on GSK3A/B activity. The total amount of GSK3beta 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 GSK3beta enzyme & AQT1213 Control:** The GSK3beta protein (0.5 nM) fully phosphorylated the AQT1211 sensor peptide substrate by 240 min., as shown by convergence with the signal obtained with the AQT1213 phosphopeptide positive control (a flat horizontal line defining the maximum RFU with this sensor peptide). The signal with the GSK3beta enzyme was eliminated by adding 1 µM LY2090314 inhibitor. The signal with AQT1213 is used to convert RFU (Corrected) values to nmoles of phosphopeptide.



# Reagent Preparation for a 384-well Format (25 $\mu$ L 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  $\mu$ L per well

**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  $\mu$ L of 10 mM ATP by adding 12  $\mu$ L of 100 mM ATP to 108  $\mu$ L of ultrapure deionized water.
  - 10 mM DTT**: Make 300  $\mu$ L of 10 mM DTT by adding 3  $\mu$ L of 1M (1000 mM) DTT to 297  $\mu$ L of ultrapure deionized water.
  - 5.5 mM EGTA**: Make 300  $\mu$ L of 5.5 mM EGTA by adding 3  $\mu$ L of 550 mM EGTA to 297  $\mu$ L of ultrapure deionized water.
  - 150  $\mu$ M Sensor Peptide**: Make 105  $\mu$ L of 150  $\mu$ M [AQT1211](#) by adding 15.75  $\mu$ L of 1 mM AQT0982 to 89.25  $\mu$ L of ultrapure deionized water.
- Prepare final 1X *PhoshoPreserve* Cell Extraction Buffer (CEB) by adding 1  $\mu$ L of 1M DTT, 16.7  $\mu$ L of the provided AQT Inhibitor Cocktail for Proteases and 16.7  $\mu$ L of the provided AQT Inhibitor Cocktail for Phosphatases to 1000  $\mu$ L 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  $\mu$ L (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 AQT1211**

Components for 1.25X Master Mix:	For 1 Well:	For 40 Wells:
Enzyme Reaction Buffer (10X)	2.5 $\mu$ L	100 $\mu$ L
ATP (10 mM)	2.5 $\mu$ L	100 $\mu$ L
DTT solution (10 mM)	2.5 $\mu$ L	100 $\mu$ L
EGTA Solution (5.5 mM)	2.5 $\mu$ L	100 $\mu$ L
Selective Sensor Peptide Substrate <a href="#">AQT1211</a> (150 $\mu$ M)	2.5 $\mu$ L	100 $\mu$ L
Ultrapure deionized water	7.5 $\mu$ L	300 $\mu$ L
<b>Total volume</b>	<b>20 <math>\mu</math>L</b>	<b>800 <math>\mu</math>L</b>

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

# Preparation of Crude Cell Lysates

## Example: Preparing A375 Cell Lysates

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

### 1. Cell Culture and Stimulation:

- a.  $4.9 \times 10^6$  A375 cells were plated in 40 mL in a T-175 tissue culture-treated flask and incubated at 37 °C in EMEM Medium with 10% FBS and 1% PenStrep in an atmosphere of 5% CO<sub>2</sub>.
- b. Cells at ~80% confluency (~  $18 \times 10^6$  cells) were washed with PBS, the residual liquid 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, adding DNAase, 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 600 µL per well of 3.6 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 until the stability is demonstrated.**

# 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  $\mu$ 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  $\mu$ 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  $\mu$ L of 1X EDB, add 1000  $\mu$ L 5X EDB Base along with 5  $\mu$ L of 1M DTT, 5  $\mu$ L of 550 mM EGTA, and 3990  $\mu$ L of ultrapure deionized water to create the final composition shown on slide 5. **Keep on ice.**
4. Prepare 60  $\mu$ L of 0.50 mg/mL of your stimulated lysate (for example, combine 30  $\mu$ L of a 1 mg/mL lysate stock with 30  $\mu$ 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% CEB) for diluting the lysate prepared above, by mixing 200  $\mu$ L of CEB with 200  $\mu$ 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  $\mu$ L of 50% EDB/50% CEB per well to wells 1-11 in a single row. Add 30  $\mu$ 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  $\mu$ L to well 10. Mix the contents of well 10 and transfer 30  $\mu$ 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% CEB.
8. Transfer 5  $\mu$ 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 $\mu$ L
Lysate	5.0 $\mu$ L

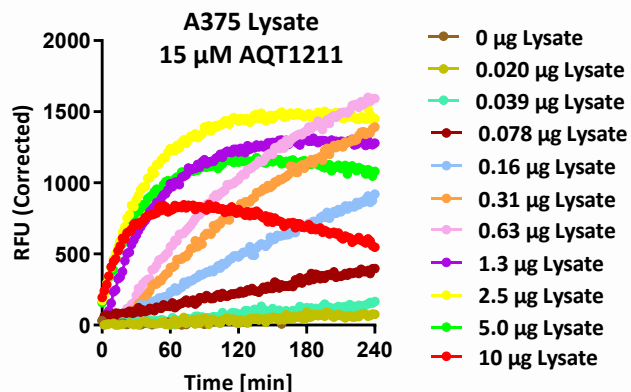
*Total volume 25  $\mu$ L*

Plate Map for Lysate Titration to Assess Dose Dependence													
		1	2	3	4	5	6	7	8	9	10	11	12
A375 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

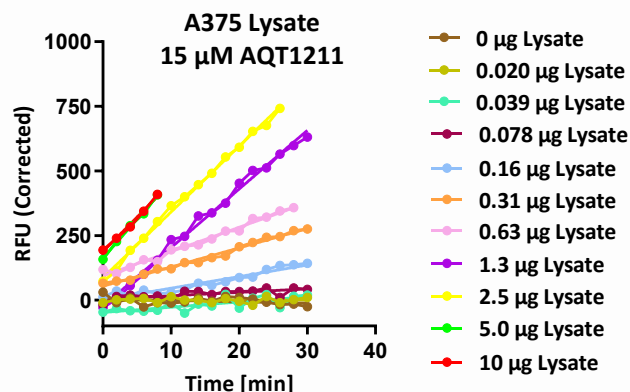
**This test uses 36 wells and only the AQT1211 sensor peptide substrate and the lysate from A375 cells. Column 1 is the Blank. All conditions are tested in triplicate.**

# Lysate Titration for A375 Cells and GSK3A/B Activity Measured with AQT1211

## 1. Full Time Course Progress Curves

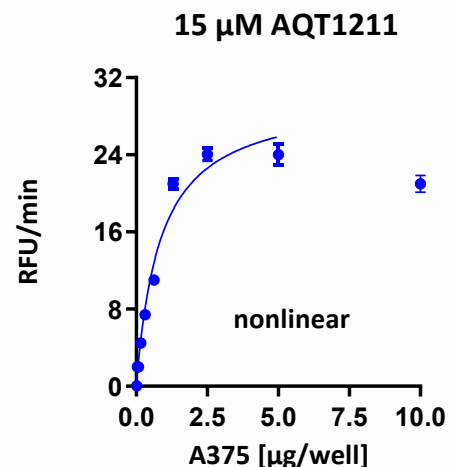


## 2. Linear Range of Progress Curves

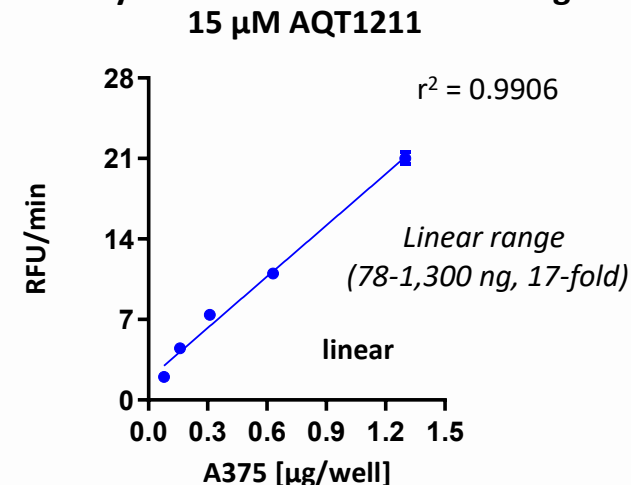


## 3. Reaction Rate vs Amount of Lysate per Well

### A. All Lysate Amounts



### B. Lysate Amounts in Linear Range



The AQT1211 sensor peptide was used at 15 µM with an increasing amount of lysate from A375 cells treated. 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-B 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 for GSK3A/B provides a selective, highly quantitative, and accurate measure of kinase activity in a complex sample.**