

# AQT0732 – MTOR Assay Validation

---

*PhosphoSens<sup>®</sup>*-Kinetic Assay Format

# Outline for this Study

## PhosphoSens–Kinetic Assay Validation

### Enzyme Source, Construct, and Lot Information:

Carna mTOR/MLST8 (Cat#/Lot#: 11-431/21CBS-0121B) amino acids 1362-2549(end); N-terminal DYKDDDDK tag

### Reference Compound Information:

Wortmannin MedChemExpress (Cat#: HY-10197)

### Experiments to be run:

Enzyme Titration

Sensor Peptide  $K_m$  Determination

ATP  $K_m$  Determination

DMSO Tolerance Test

Reference Compound  $IC_{50}$  Determination at ATP  $K_m$

# Enzyme Titration

## 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

0.55 mM EGTA

10 mM MgCl<sub>2</sub>

20 μM AQT0732

0, 0.040, 0.080, 0.16, 0.31, 0.63, 1.3, 2.5, 5.0, 10, 20, and 40 nM mTOR/MLST8

### Reaction Set Up:

2 or 2.5 μL

10x Sensor Peptide

14 or 17.5 μL

Reaction Mix with ATP & DTT

4 or 5 μL

1x EDB or Kinase dilutions (5x in EDB)

20 or 25 μL

Final reaction volume

Reactions were run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat round 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.

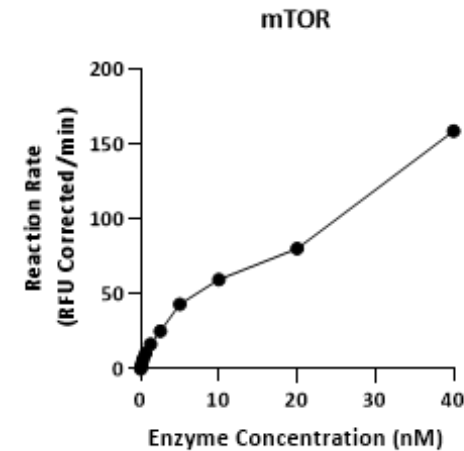
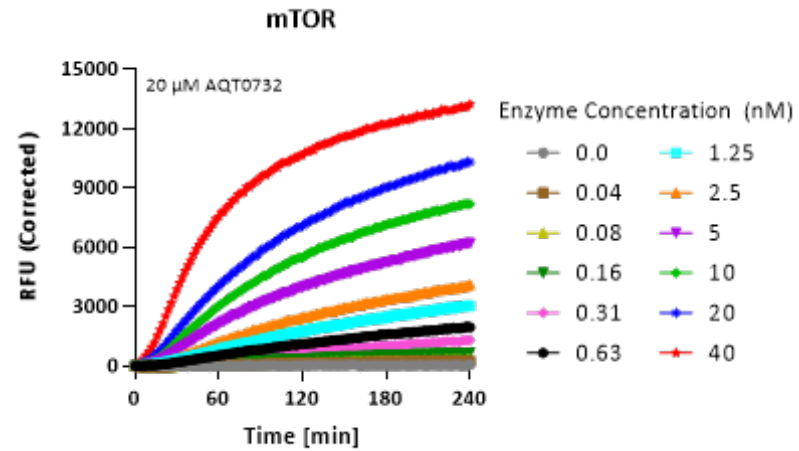
### Notes:

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.

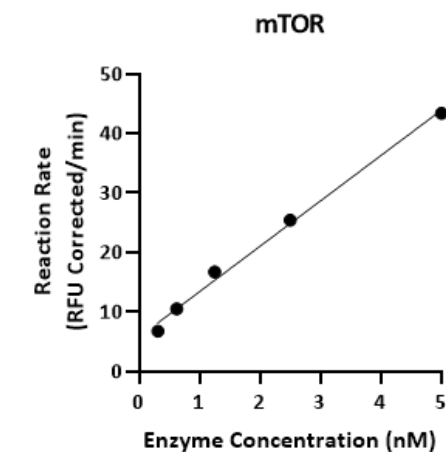
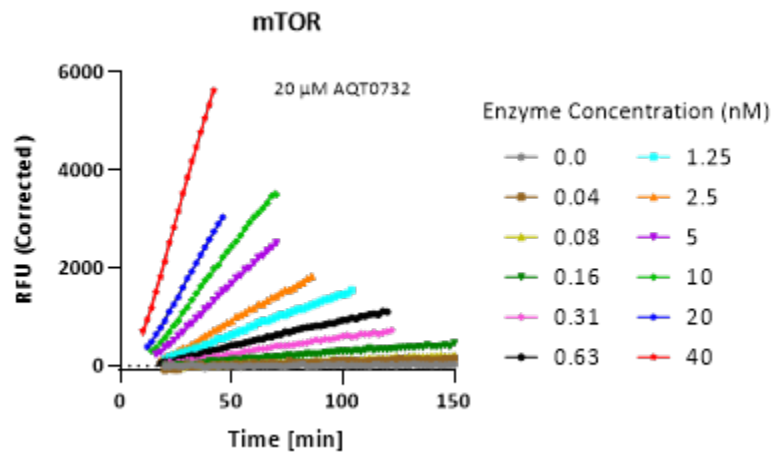
# Enzyme Titration

## Progress Curves

Complete  
Progress  
Curves



Linear  
Region of  
Progress  
Curves



Linear  
Range

# Enzyme Titration

## Reaction Rate Table

Enzyme Conc. (nM)	Reaction Rate (RFU/min)	Normalized Reaction Rate (RFU/pmole/min)
0.04	1.4	1788
0.08	1.9	1240
0.16	3.4	1089
0.31	6.8	1082
0.63	11	842
1.3	17	669
2.5	25	508
5.0	43	434
10	60	299
20	81	201
40	159	199

The reaction is linear from 0.63–5.0 nM

# 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

0.55 mM EGTA

10 mM  $MgCl_2$

0, 0.20, 0.39, 0.78, 1.6, 3.1, 6.3, 13, 25, 50, 100, and 200  $\mu M$  AQT0732

8 nM mTOR/MLST8

### Reaction Set Up:

2 or 2.5  $\mu L$

10x Sensor Peptide

14 or 17.5  $\mu L$

Reaction Mix with ATP & DTT

4 or 5  $\mu L$

1x EDB or Kinase dilutions (5x in EDB)

20 or 25  $\mu L$

Final reaction volume

Reactions were run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat round bottom polystyrene NBS microplates (Cat. #3824) at 25  $\mu L$  final well volume or in PerkinElmer, ProxiPlate-384 Plus, white shallow well microplates (Cat. #6008280) at 20  $\mu L$  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.

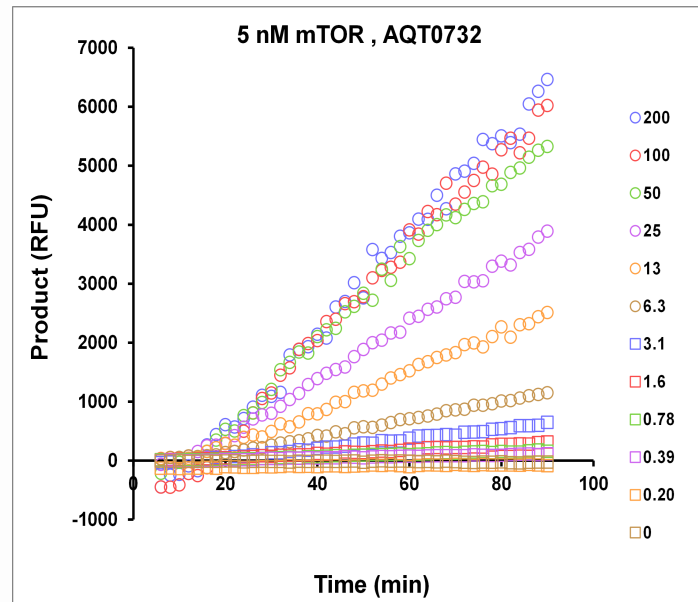
### Notes:

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.

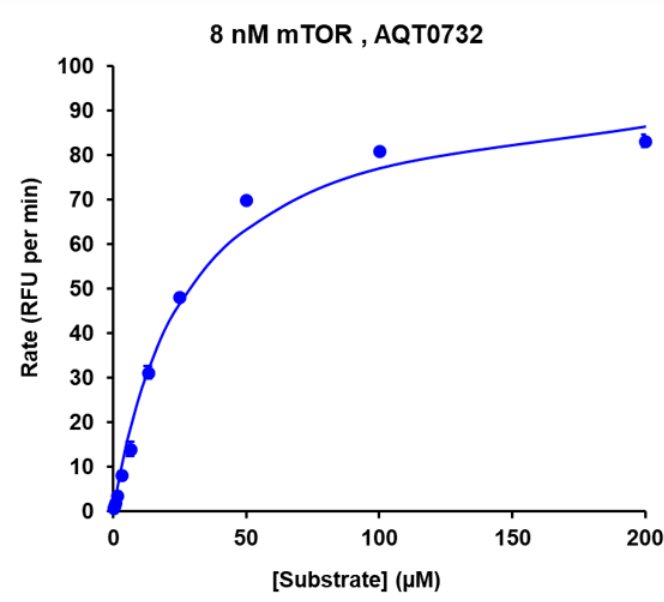
# Sensor Peptide $K_m$ Determination

Titration Curves and  $K_m$  Plot and Table

## Sensor Peptide Titration Curves



## Sensor Peptide $K_m$ Plot



## Sensor Peptide $K_m$ Table

Parameter	Value	Approx SE
Vmax (RFU per min)	98.5	3.3
Vmax (RFU per pmol per min)	616	21
Km ( $\mu\text{M}$ )	27.8	0.9
R squared	0.992	

Sensor Peptide  $K_m$  is 28  $\mu\text{M}$

# ATP $K_m$ Determination

## Reaction Conditions and Set Up

### Reaction Conditions:

54 mM HEPES, pH 7.5  
0, 2.0, 3.9, 7.8, 16, 31, 63, 125, 250, 500, 1000, and 2000  $\mu$ M ATP  
1.2 mM DTT  
0.012% Brij-35  
1% glycerol  
0.2 mg/ml BSA  
0.55 mM EGTA  
10 mM  $MgCl_2$   
20  $\mu$ M AQT0732  
8 nM mTOR/MLST8

### Reaction Set Up:

2 or 2.5 $\mu$ L	10x Sensor Peptide
14 or 17.5 $\mu$ L	Reaction Mix with ATP & DTT
<u>4 or 5 <math>\mu</math>L</u>	1x EDB or Kinase dilutions (5x in EDB)
20 or 25 $\mu$ L	Final reaction volume

Reactions were run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat round bottom polystyrene NBS microplates (Cat. #3824) at 25  $\mu$ L final well volume or in PerkinElmer, ProxiPlate-384 Plus, white shallow well microplates (Cat. #6008280) at 20  $\mu$ L 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:

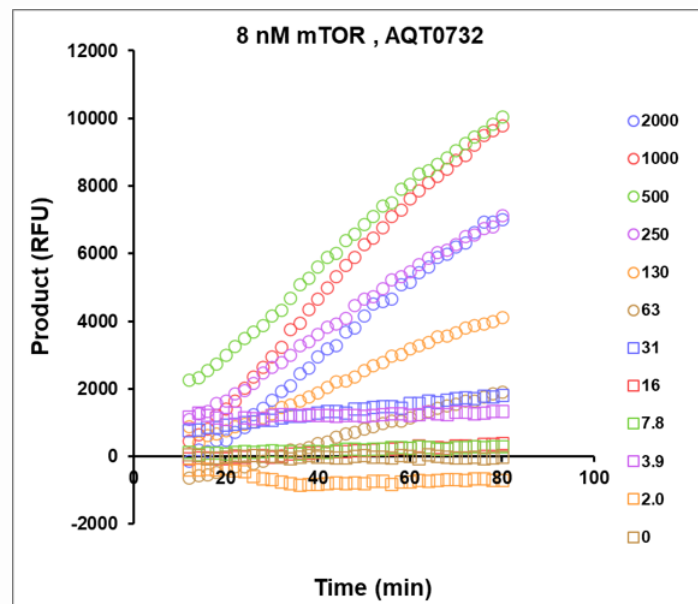
Enzyme Dilution Buffer (EDB): 20 mM HEPES, pH 7.5, 0.01% Brij-35, 5% Glycerol, 0.5 mM EGTA, 1 mM DTT, 1 mg/ml Bovine Serum Albumin.



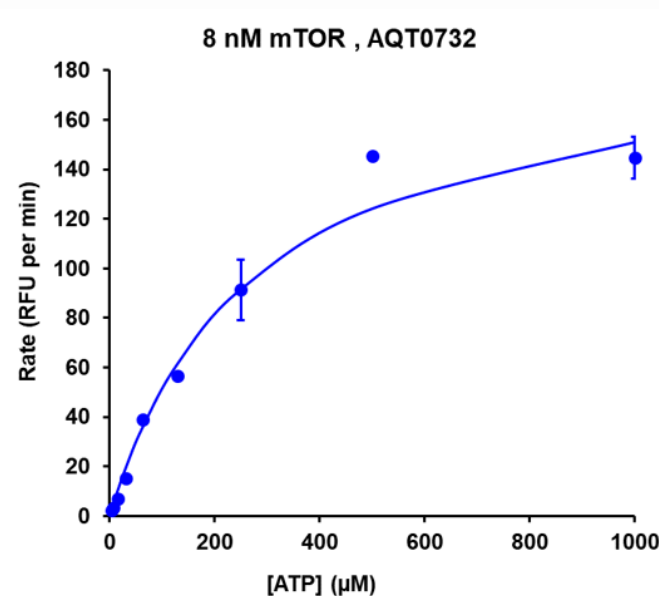
# ATP $K_m$ Determination

Titration Curves and  $K_m$  Plot and Table

## ATP Titration Curves



## ATP $K_m$ Plot



## ATP $K_m$ Table

Parameter	Value	Approx SE
$V_{\text{max}}$ (RFU per min)	192.8	3.7
$V_{\text{max}}$ (RFU per pmol per min)	1205	23
$K_m$ ( $\mu\text{M}$ )	276.6	23.4
R squared	0.982	

ATP  $K_m$  is 277  $\mu\text{M}$

# DMSO Tolerance Test

## 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

0.55 mM EGTA

10 mM MgCl<sub>2</sub>

0, .01, .02, .04, .08, .16, .31, .63, 1.3, 2.5, 5.0, and 10% DMSO

40 μM AQT0732

5 nM mTOR/MLST8

### Reaction Set Up:

2 or 2.5 μL

10x DMSO dilutions

14 or 17.5 μL

Reaction Mix with Sensor Peptide, ATP & DTT

4 or 5 μL

1x EDB or Kinase dilutions (5x in EDB)

20 or 25 μL

Final reaction volume

Reactions were run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat round bottom polystyrene NBS microplates (Cat. #3824) at 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.

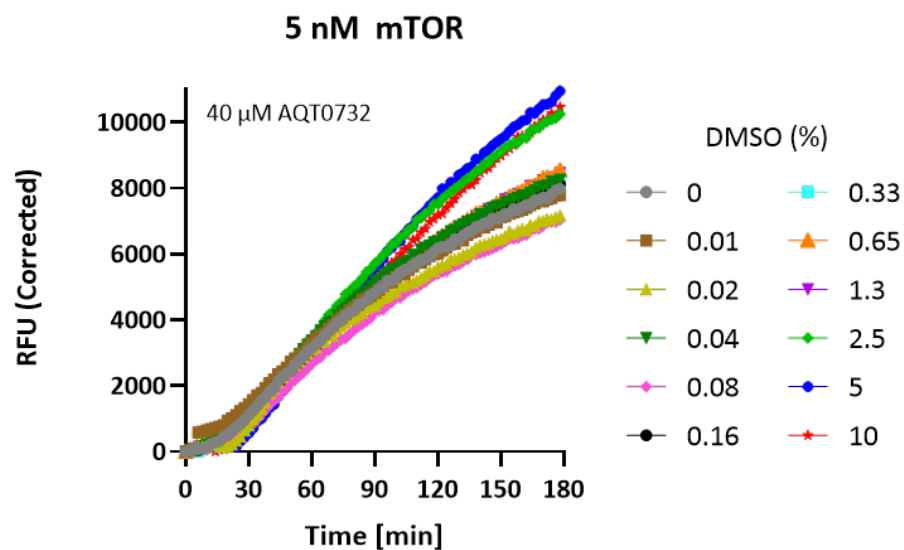
### Notes:

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.

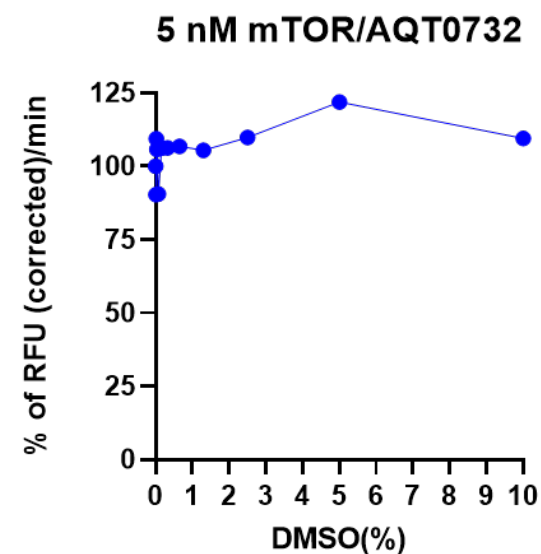
# DMSO Tolerance Test

Titration Curves and Inhibition Plot

## Complete Progress Curves



## Reaction Rate vs [DMSO] Plot



No change in enzyme activity out to 3% DMSO

# IC<sub>50</sub> Determination

## Reaction Conditions and Set Up

### Reaction Conditions:

54 mM HEPES, pH 7.5

ATP at K<sub>m</sub>

1.2 mM DTT

0.012% Brij-35

1% glycerol

0.2 mg/ml BSA

0.55 mM EGTA

10 mM MgCl<sub>2</sub>

2% DMSO

40 μM AQT0732

5 nM mTOR/MLST8

0, 0.17, 0.51, 1.5, 4.6, 14, 41, 123, 370, 1,110, 3,330, and 10,000 nM Wortmannin

### Reaction Set Up:

16 μL            Reaction Mix with Sensor Peptide and Inhibitor

4 μL            1x EDB or Kinase dilutions (5x in EDB)

20 μL            Final reaction volume

Reactions were run at 30°C for 240 minutes in either Corning, low volume 384-well, white flat round bottom polystyrene NBS microplates (Cat. #3824) at 20 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.

Inhibitors are added via direct (0.4 μL of 50X stock in 100% DMSO) or intermediate dilutions (2.0 μL of 10X stock in 10% DMSO).

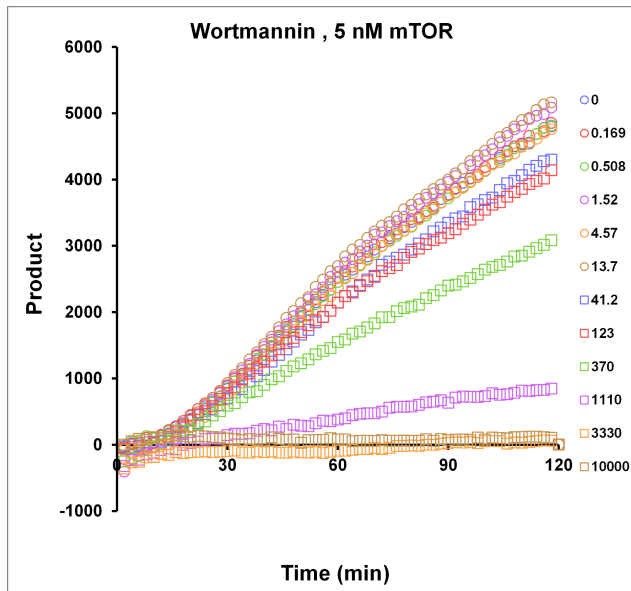
### Notes:

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.

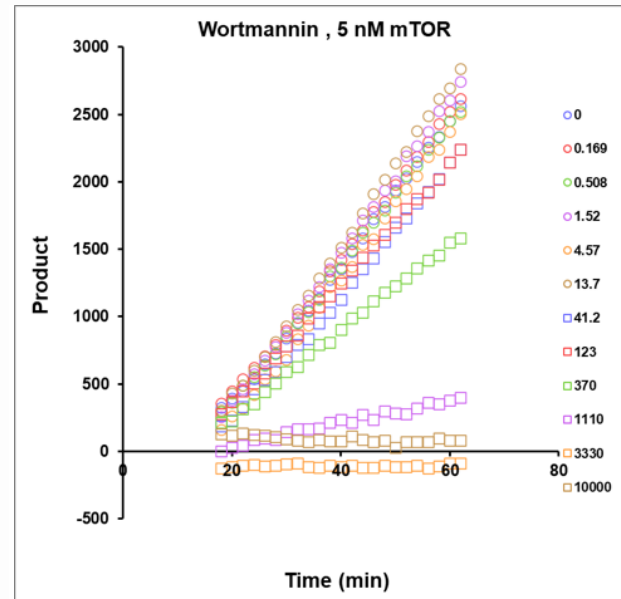
# IC<sub>50</sub> Determination

## Progress Curves and IC<sub>50</sub> Curves and Table

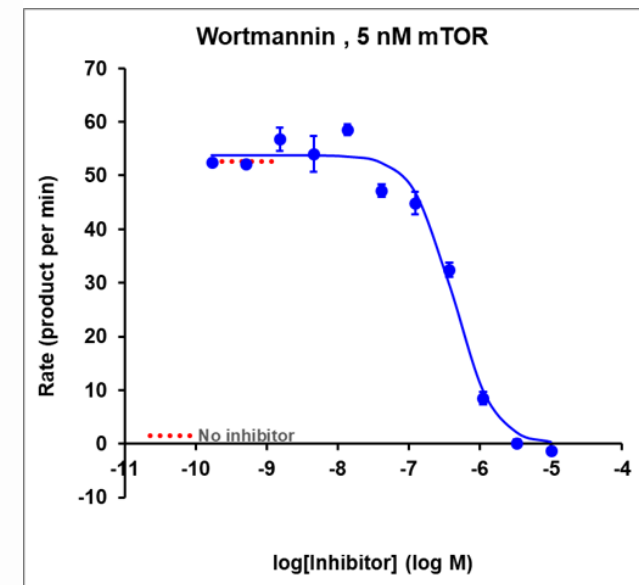
### Inhibitor Titration Progress Curves



### Linear Region of Progress Curves



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



The Y-axis label is RFU/min.

### IC<sub>50</sub> Table

Parameter	Value
Bottom	0.0
Top	53.7
log IC50	-6.37
IC50 (nM)	425.93
Ki (nM)	212.97
Slope	-1.544
R squared	0.985
IC50 approx SE (nM)	17.80
50% inhibition (nM)	425.93

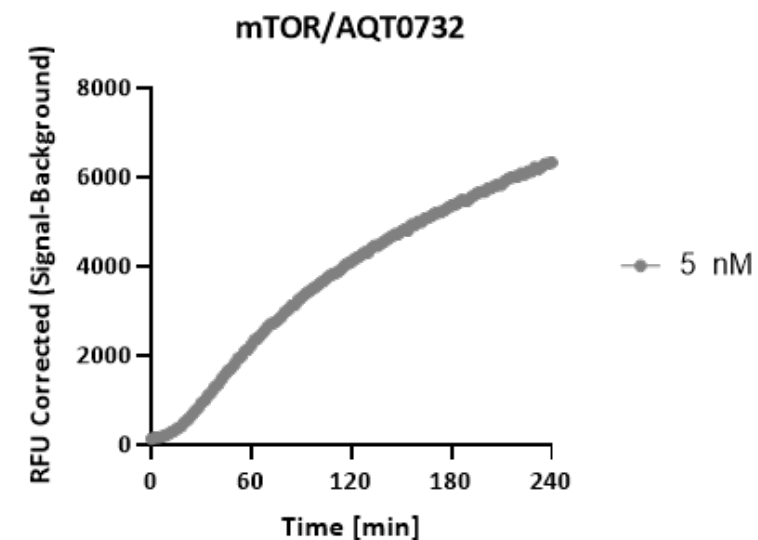
Wortmannin IC<sub>50</sub> at ATP K<sub>m</sub> is 426 nM

# Summary

## Assay Validation Results and Progress Curve and Assay Strength at 1 mM ATP

Experiment	Result
Enzyme Titration Linear Range	0.63-5.0 nM
Sensor Peptide $K_m$ Value	28 $\mu$ M
ATP $K_m$ Value	277 $\mu$ M
DMSO Tolerance (highest % recommended)	1
Wortmannin $IC_{50}$ Determination at ATP $K_m$	426 nM

## Progress Curve



Kinase Name	Conc. (nM)	Sox-based Substrate Name	Normalized Reaction Rate (RFU/pmole/min)	Normalized Rate Std Error (RFU/pmole/min)
mTOR	5.0	AQT0732	434	3

Assay Strength Key	
Very Strong	> 1,000 (RFU/pmole/min)
Strong	300 to 999 (RFU/pmole/min)
Moderate	100 to 299 (RFU/pmole/min)
Weak	30 to 99 (RFU/pmole/min)

Under the conditions utilized for this experiment, the assay is Strong