

AQT1076 – MAPK3/1 (ERK1/2) Assay Validation

PhosphoSens[®]-Lysate Assay Format

Outline for this Study

PhosphoSens-Lysate Assay Validation

Lysate Source:

NIH-3T3 Cells (ATCC, CRL-1658.2) +/- 25 ng/mL PDGF-bb (ThermoFisher, 100-14B-10UG)

See slides 11-13 for Preparation of Crude Cell Lysates from NIH3T3 Cells Treated +/- PDGF to Show Activation of ERK1/2

Reference Compound Information:

SCH772984 (MedChemExpress, HY-50846)

Vx-11e (MedChemExpress, HY-14178)

Experiments to be run:

NIH-3T3 cell lysate (activated with 25 ng/mL PDGF) titration

AQT1076 substrate Km determination

DMSO Tolerance Test

Reference Compound IC₅₀ Determinations

Lysate 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

0.54 mM EGTA

10 mM MgCl₂

15 μM AQT1076

0, 0.020, 0.039, 0.078, 0.16, 0.31, 0.63, 1.3, 2.5, 5.0, 10, and 20 μg/well
NIH-3T3 (activated with 25 ng/mL PDGF)

Reaction Set Up:

20 μL Reaction Mix with AQT1076, ATP, & DTT

5 μL Enzyme dilution buffer (EDB) with ERK1/2 Lysate Buffer (1x) or ERK1/2 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 round 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.

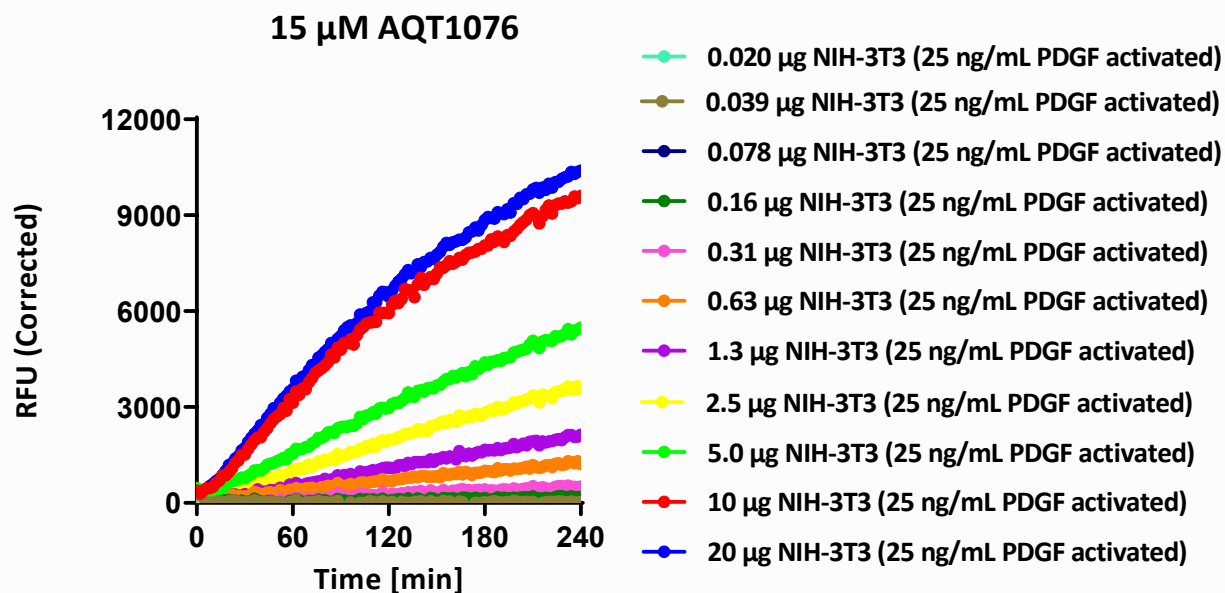
Notes:

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.

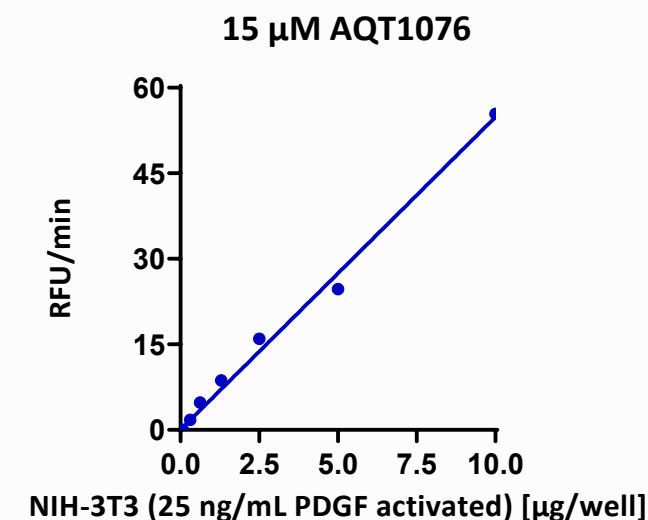
Lysate Titration

Progress Curves

Complete Progress Curves



Linear Region of Progress Curves



The ERK1/2 Lysate Assay is linear from 0.63 – 10 μ g/well lysate

ERK1/2 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

10 mM MgCl₂

0, 0.20, 0.39, 0.78, 1.6, 3.1, 6.3, 13, 25, 50, and 100 μ M AQT1076

1.0 μ g/well NIH-3T3 (activated with 25 ng/mL PDGF)

Reaction Set Up:

5 μ L 5X AQT1076 Substrate dilutions

15 μ L Reaction Mix with ATP & DTT

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

25 μ L Final reaction volume

Reaction was 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 μ L final well volume or in PerkinElmer, ProxiPlate-384 Plus, white shallow well microplates (Cat. #6008280) at 20 μ 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.

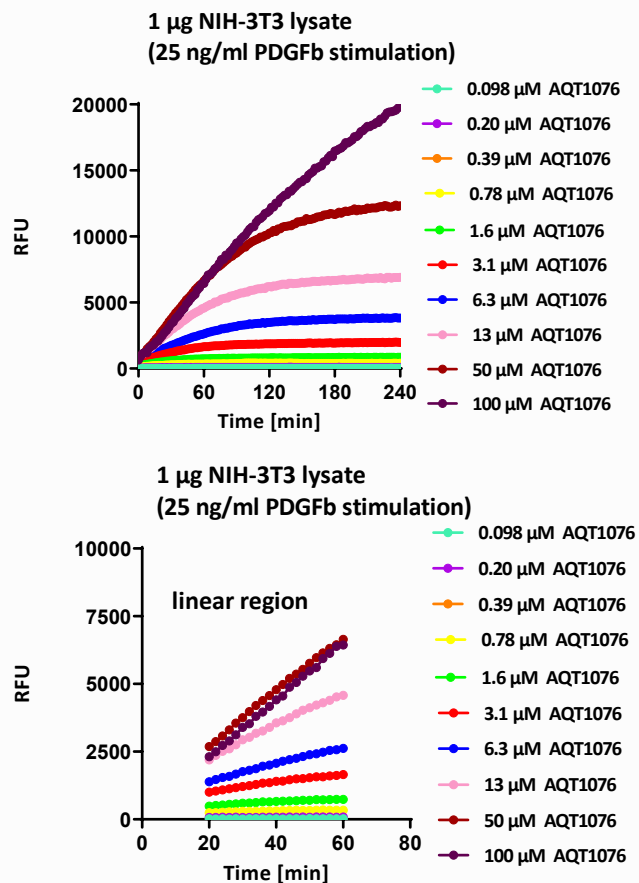
Notes:

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.

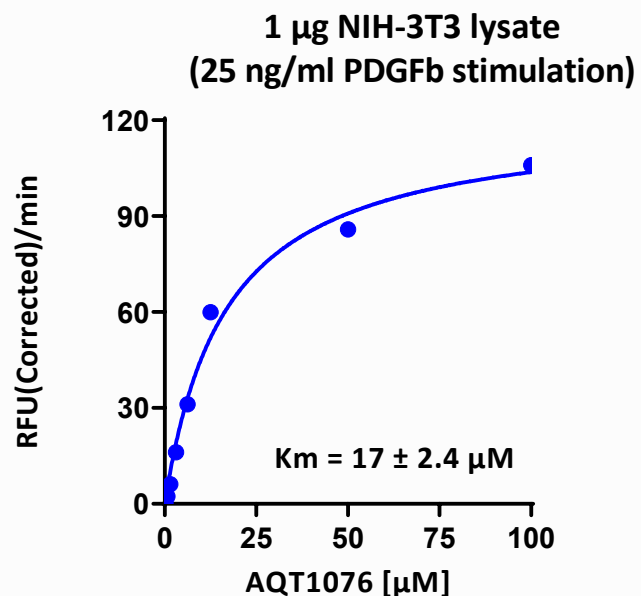
Sensor Peptide K_m Determination

Titration Curves and K_m Plot and Table

Sensor Peptide Titration Curves



Sensor Peptide K_m Plot



Sensor Peptide K_m Table

Michaelis-Menten
Best-fit values

Vmax 121.0
Km 16.65

Std. Error
Vmax 5.875
Km 2.430

95% CI (asymptotic)
Vmax 107.5 to 134.6
11.04 to 22.25

Km
Goodness of Fit 0.9898
Degrees of Freedom 8
R squared 0.9898
Sum of Squares 142.2
Sy.x 4.216

	AQT1076 K_m (μ M)
NIH-3T3 (25 ng/mL stimulated) Lysate	17
Recombinant FL tagless ERK1 (Sino, M29-10U)	9.3
Recombinant FL GST-ERK1 (Sino, M29-10G)	9.3
Recombinant FL GST-ERK2 (Sino, M28-10G)	14

The K_m value for AQT1076 is 17 μ M, similar to the K_m values observed for recombinant ERK1 and ERK2.

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

0.54 mM EGTA

10 mM MgCl₂

15 μM AQT1076

1.0 μg/well NIH-3T3 (activated with 25 ng/mL PDGF)

0-10% DMSO

Reaction Set Up:

2.5 μL 10X DMSO Titration

17.5 μL Reaction Mix with CSx Substrate, ATP & DTT

15 minutes incubation at 30°C

5 μL Enzyme dilution buffer (EDB) (1x) or Kinase (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 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.

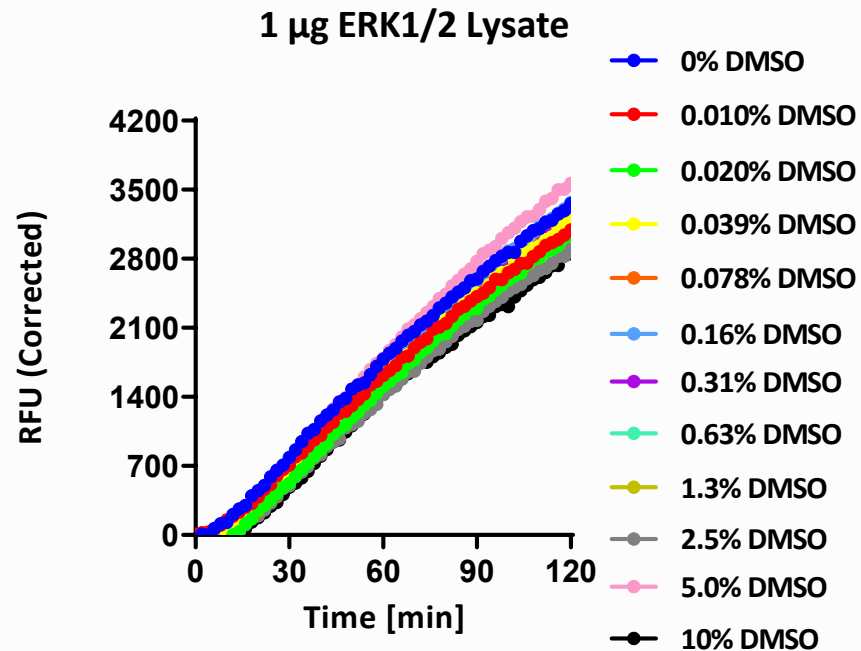
Notes:

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.

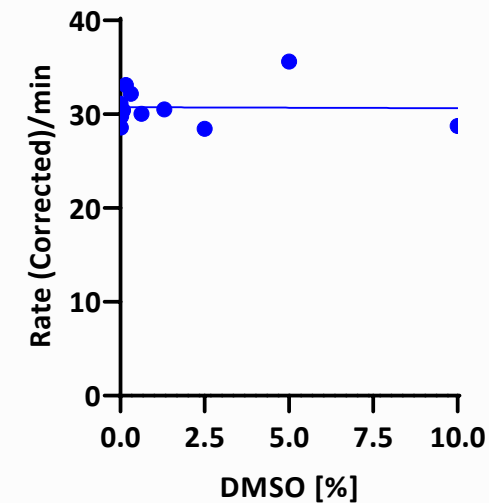
DMSO Tolerance Test

Titration Curves and Inhibition Plot

Complete Progress Curves



Reaction Rate vs [DMSO] Plot



No change in enzyme activity out to 10% DMSO

IC₅₀ Determination

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
10 mM MgCl₂
15 μM AQT1076
2% DMSO
0-1.0 μM SCH772984 (MedChemExpress, HY-50846)
0-1.0 μM Vx-11e (MedChemExpress, HY-14178)
1.0 μg/well NIH-3T3 (activated with 25 ng/mL PDGF)

Reaction Set Up:

0.5 μL 50X Staurosporine dilutions in 100% DMSO
19.5 μL Reaction Mix with CSx Substrate, ATP & DTT
15 minutes incubation at 30°C
5 μL Enzyme dilution buffer (EDB) (1x) or Kinase (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 round 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.

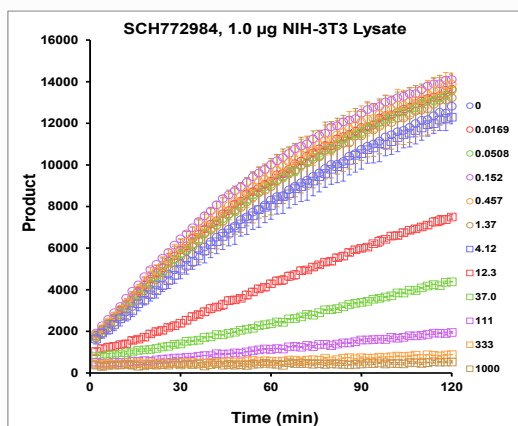
Notes:

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.

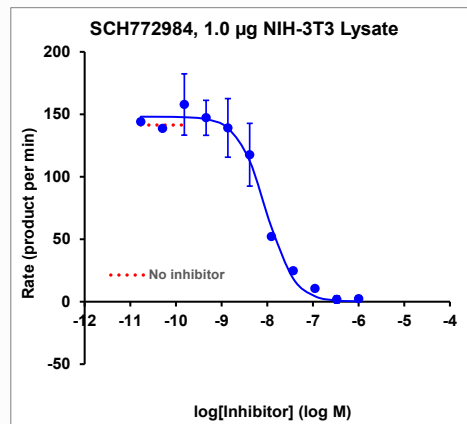
IC₅₀ Determination

Progress Curves

SCH772984



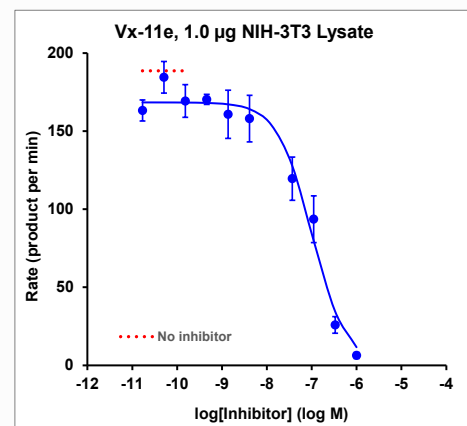
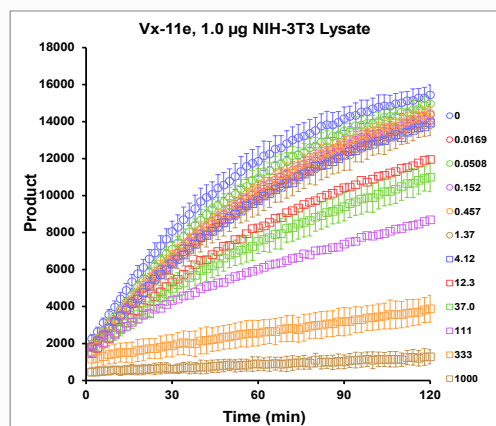
IC₅₀ Curve



IC₅₀ Table

Compound ID	Highest conc. (mM)	Data points	IC ₅₀ (nM)	% Inh @ curve min	% inh at highest conc.	Min rate (product / min)	Max rate (product / min)	R squared
SCH772984	1.0	22	9.3	2.4	98	0	148	0.993
Vx-11e	1.0	20	101	6	96	0	169	0.980

Vx-11e

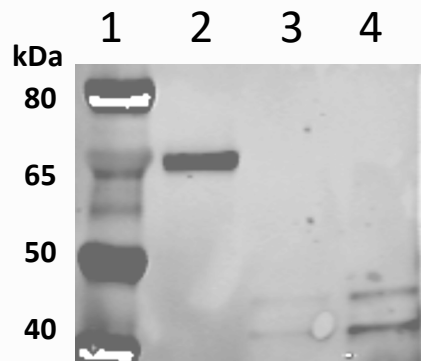


Preparation of Crude Cell Lysates from NIH3T3 Cells Treated +/- PDGF to Show Activation of ERK1/2

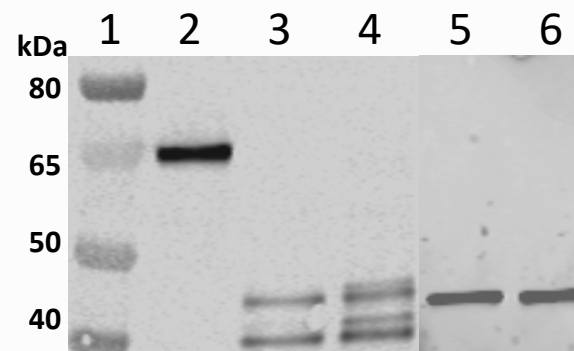
- 1) NIH-3T3 Cells (ATCC, CRL-1658.2) were plated in 6-well tissue culture-treated plates and incubated for 48 hours at 37°C in DMEM Medium with 10% FBS and PenStrep in an atmosphere of 5% CO₂. Cells were then serum-starved in culture medium with 0.1% FBS (ThermoFisher, A56708-01) for 24 hours and incubated for 15 minutes with or without 25 ng/mL PDGF-bb (ThermoFisher, 100-14B-10UG). Cells were then washed with PBS, and lysed with lysis buffer containing:
 - 50 mM HEPES, pH 7.4
 - 150 mM NaCl
 - 2 mM EGTA
 - 1 mM DTT
 - 1% Triton X-100
 - 30 mM NaF
 - 10 mM Na₄P₂O₇
 - 100 μM Na₃VO₄
 - 50 mM β-glycerophosphate
 - Protease Inhibitor Cocktail diluted 60-fold into lysis buffer (Sigma-Aldrich product no. #P8340)
 - Phosphatase Inhibitor Cocktail diluted 60-fold into lysis buffer (Sigma-Aldrich product no. #P2850)
- 2) The DNA strands were broken by briefly sonicating on ice for 2 seconds on low power. Lysates were used immediately or aliquoted and frozen at -80°C. 1 μg of each Lysate was run on a gel, transferred to a nitrocellulose membrane, and the signal developed by Western blotting using antibodies to total ERK1/2 (Cell Signaling, #4696) and phospho-ERK1/2 (T202/Y204) (Cell Signaling, #4370).

ERK1/2 Lysate Western Blots

Phospho-ERK1/2



Total ERK1/2 and Actin Loading Control

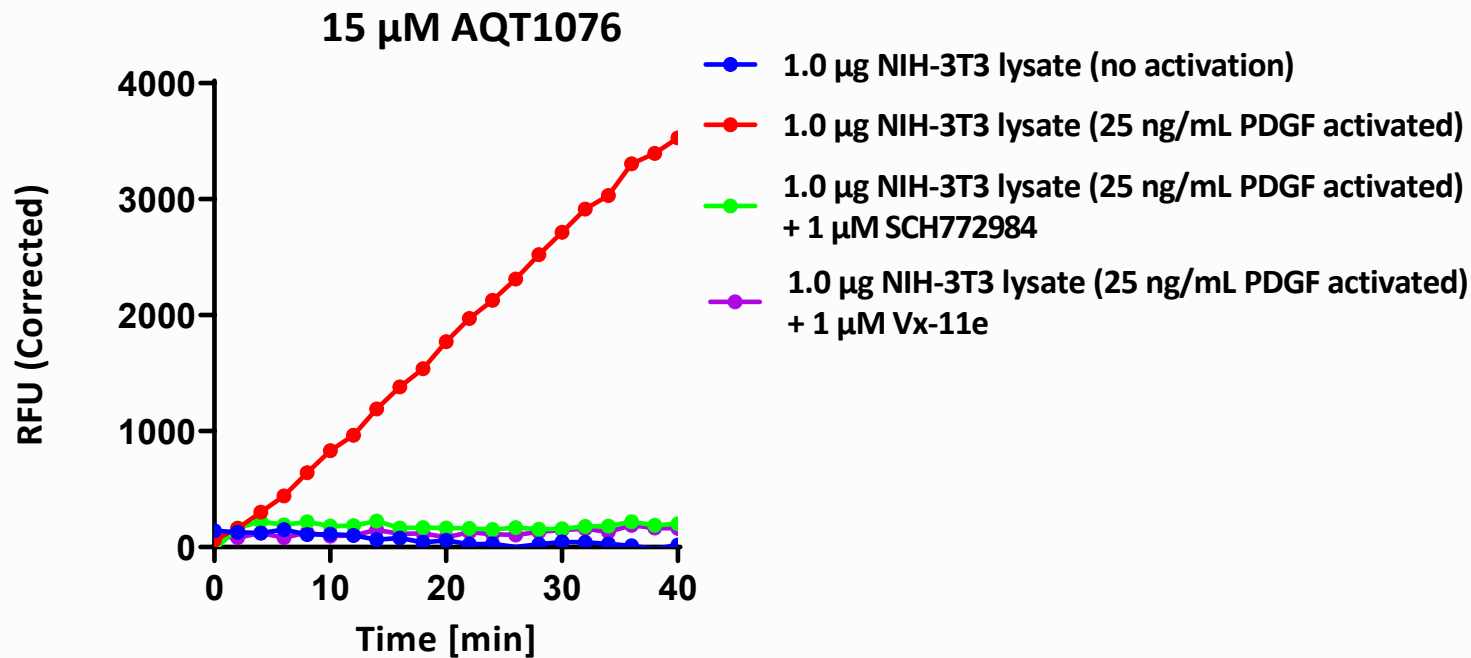


Lane Description:

1. MW Markers
2. Recombinant ERK2
3. NIH-3T3 (no activation)
4. NIH-3T3 (+ 25 ng/mL PDGF for 15 minutes)
5. NIH-3T3 (no activation) – Loading Control
6. NIH-3T3 (+ 25 ng/mL PDGF for 15 minutes) – Loading Control

Membranes were blocked in TBST with 5% BSA for 2 hours and then incubated for 2 hours with anti-ERK1/2 (Cell Signaling Technology, 4696) and anti-phospho-ERK1/2 (Cell Signaling Technology, 4370) diluted 2000-fold in TBST with 2% BSA (Lanes 2-4) or anti-beta actin (Cell Signaling Technology, 4970) diluted 5000-fold in TBST with 2% BSA (Lanes 5-6). Membranes were washed 3 times for 10 minutes each in TBST and then incubated for 2 hours with goat anti-mouse antibody (LI-COR, IR Dye 800CW 926-32210) and/or donkey anti-rabbit antibody (LI-COR, IR Dye 680RD 926-68073) diluted 20000-fold in TBST with 2% BSA. Membranes were washed 3 times for 10 minutes each in TBST and then read in a LI-COR Odyssey.

ERK1/2 Lysate Activity Assay



NIH-3T3 PDGF activation significantly increases ERK1/2 lysate activity which is inhibited by 1.0 μ M of both SCH772984 and Vx-11e.

	Reaction Rate (RFU/min)	Change
1.0 μ g NIH-3T3 lysate (no activation)	0	-
1.0 μ g NIH-3T3 lysate (+25 ng/mL PDGF activated)	92	92-fold increase above control
1.0 μ g NIH-3T3 lysate (+25 ng/mL PDGF activated), + 1 μ M SCH772984	1.1	99% inhibition of PDGF activation
1.0 μ g NIH-3T3 lysate (+25 ng/mL PDGF activated), + 1 μ M Vx-11e	1.8	98% inhibition of PDGF activation

Summary

NIH-3T3 cell activation with 25 ng/mL PDGF-bb increases ERK1/2 phosphorylation and subsequent activity which is inhibited fully by 1 μ M of both SCH772984 and Vx-11e

The ERK1/2 Lysate Assay with the AQT1076 selective sensor peptide demonstrates a robust and more physiologically relevant assay using the PhosphoSens Platform:

- ❖ The ERK1/2 Lysate titration was linear from 0.31 to 10 μ g/well.
- ❖ Sensor peptide substrate AQT1076 has a K_m of 17 μ M.
- ❖ The IC_{50} values for SCH772984 and Vx-11e were 9.3 nM and 101 nM, respectively.