

HTRF from well to clinic: development of a multi-purpose phos-VASP assay

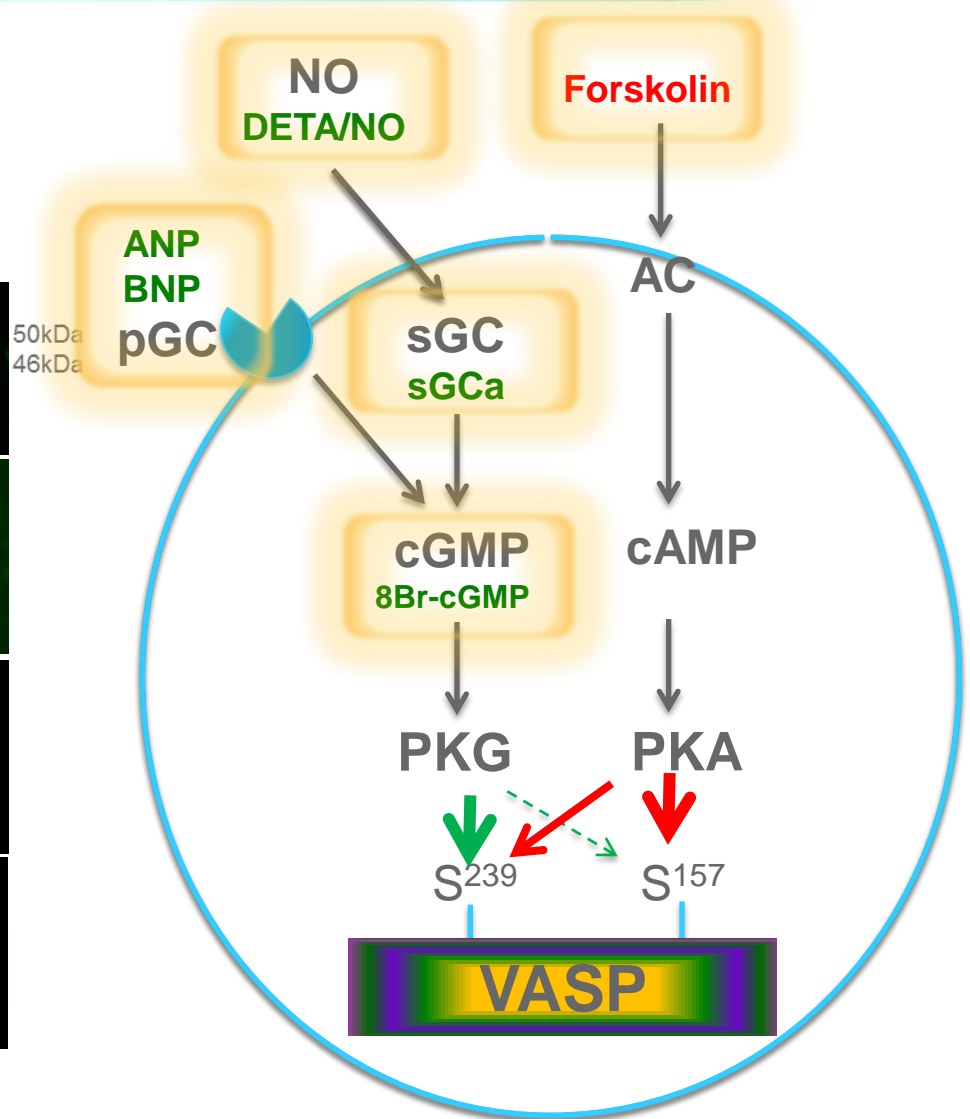
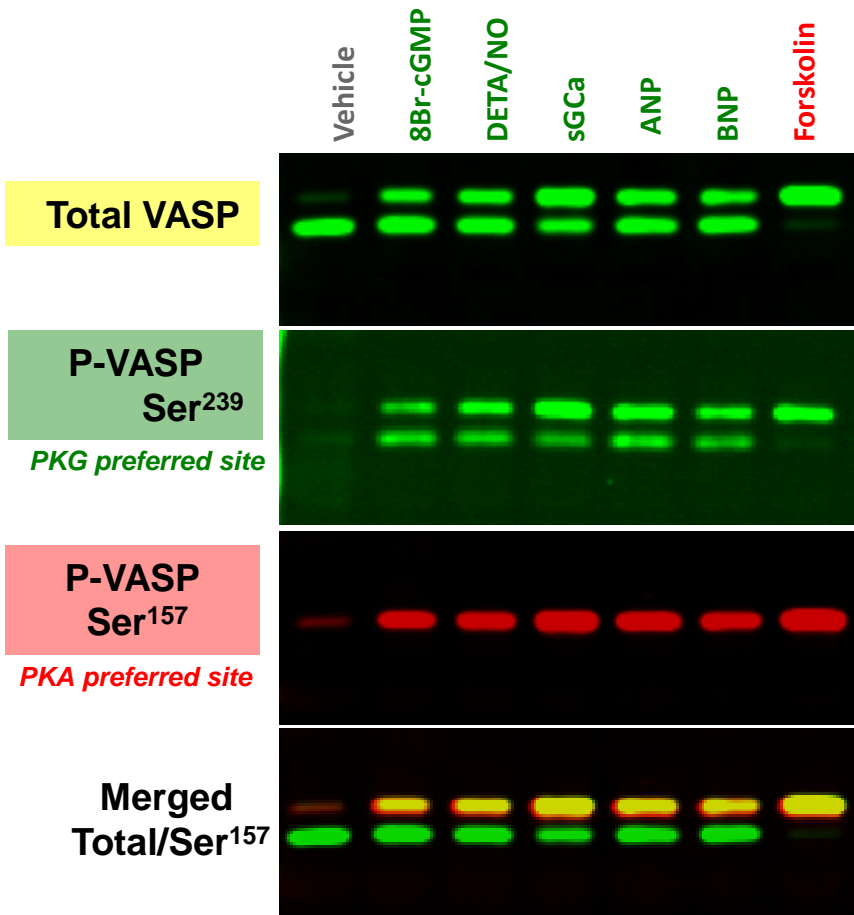
Andrea Basso¹, David Bauman², Beata Zamlynny¹, Debra McGuinness¹, Richard Visconti¹, Aida Rivera-Gines¹ Steve Fried¹, Frederick Monsma¹

¹In Vitro Pharmacology ²Cardiometabolic Diseases,
Merck Research Laboratories, Kenilworth NJ

VASP

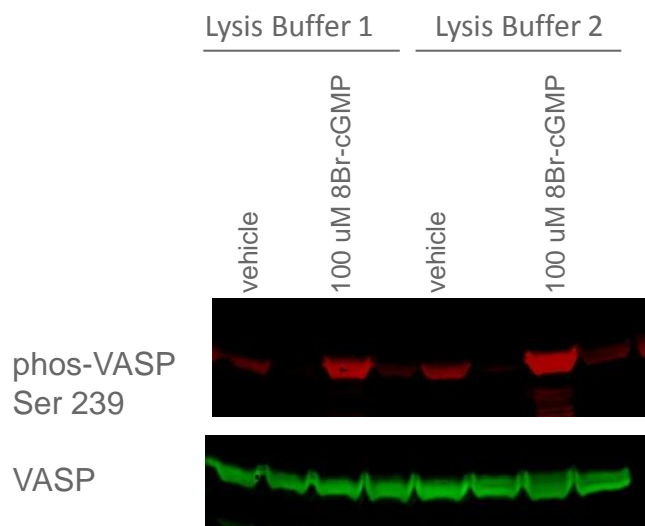
- ▶ VASP= vasodilator-stimulated phosphoprotein
- ▶ Regulates actin cytoskeleton dynamics and plays a role in smooth muscle contraction.
- ▶ Modulated by cAMP and cGMP pathways
- ▶ We sought to develop a phos-VASP HTRF assay for both Ser157 and Ser239 phosphorylation in order to evaluate signaling downstream of cAMP and cGMP pathways.
- ▶ Goals:
 - Develop an assay to be utilized for routine screening
 - Miniaturize
 - Minimize plate handling
 - Robust signal
 - High Z-prime
 - Assay platform translatable to physiologically relevant cells
 - Assay platform translatable to in vivo studies
 - Potential for human clinical biomarker studies

Differentiating cGMP and cAMP Pathways



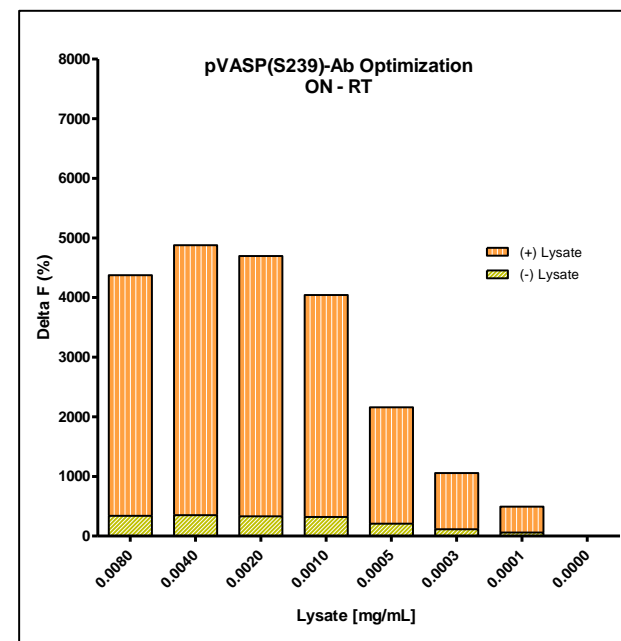
Custom Assay Development for VASP Phos-Ser239

Hek293 Cells VASP Transient Transfection



- ▶ Tested 2 total VASP and 3 phos-VASP antibodies; each labeled with donor and acceptor using 3 concentrations of lysate samples
- ▶ Optimized concentration of each antibody using 3 concentrations each antibody in matrix fashion on 7 concentrations of lysate samples
- ▶ Final selection tested on titration of lysate samples (graphed)
 - Monoclonal phospho-VASP (Ser239) labeled d2
 - Monoclonal VASP labeled K
- ▶ Significant signal-to-noise between treated and untreated samples

ANTIBODIES		
ID#	Name	Clonality
1	VASP Antibody-K	Poly
2	VASP Antibody-d2	Poly
3	VASP-K	Mono
4	VASP-d2	Mono
5	Anti-pVASP (Ser239)-K	Mono
6	Anti-pVASP (Ser239)-d2	Mono
7	VASP Antibody (Ab-239)-K	Poly
8	VASP Antibody (Ab-239)-d2	Poly
9	VASP (phospho-Ser239)-K	Mono
10	VASP (phospho-Ser239)-d2	Mono



Custom Assay Development for VASP Phos-Ser157

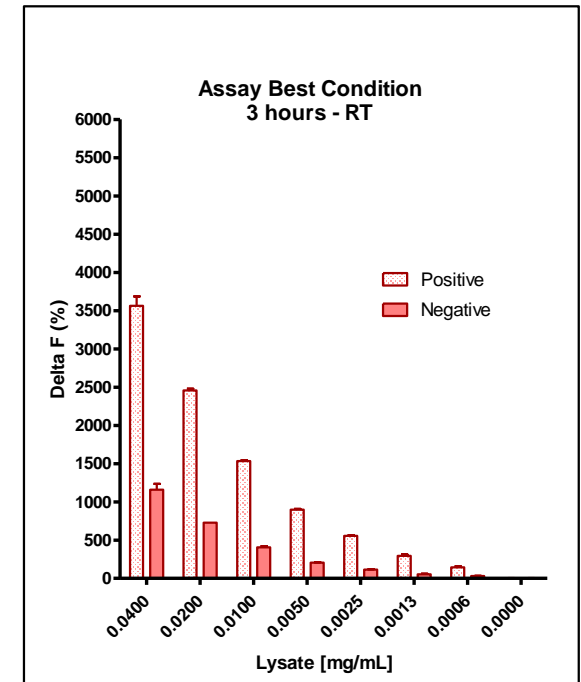
Hek293 Cells VASP Transient Transfection



- ▶ Tested 2 total VASP and 3 phos-VASP antibodies; each labeled with donor and acceptor using 3 concentrations of lysate samples
- ▶ Optimized concentration of each antibody using 3 concentrations each antibody in matrix fashion on 7 concentrations of lysate samples
- ▶ Final selection tested on titration of lysate samples (graphed)
 - Monoclonal phosho-VASP labeled K
 - Monoclonal VASP labeled d2
- ▶ Significant signal-to-noise between treated and untreated samples

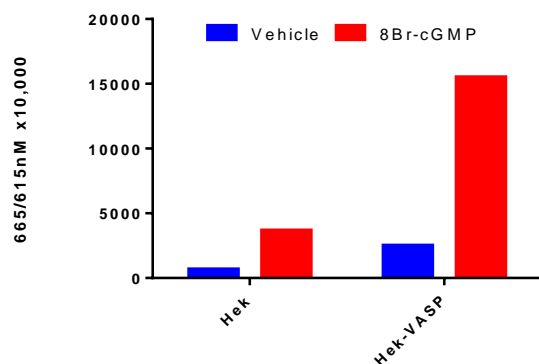
HTRF Assay Development (Cisbio)

ANTIBODIES		
ID#	Name	Clonality
1	VASP Antibody-K	Poly
2	VASP Antibody-d2	Poly
3	VASP-K	Mono
4	VASP-d2	Mono
5	Phospho-VASP (Ser157)-K	Poly
6	Phospho-VASP (Ser157)-d2	Poly
7	Anti-VASP (phospho S157)-K	Poly
8	Anti-VASP (phospho S157)-d2	Poly
9	Anti-VASP (phospho S157)-K	Mono
10	Anti-VASP (phospho S157)-d2	Mono

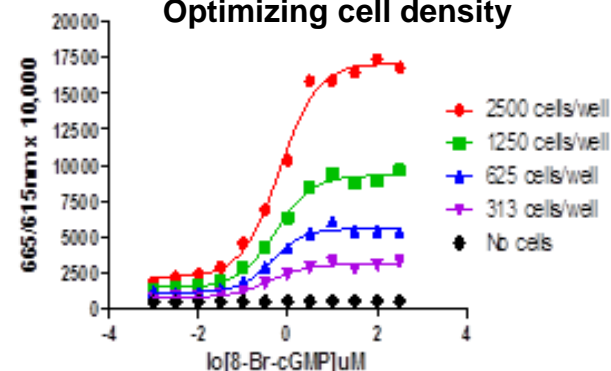


Optimizing Cell Type, Cell Density and Time Course

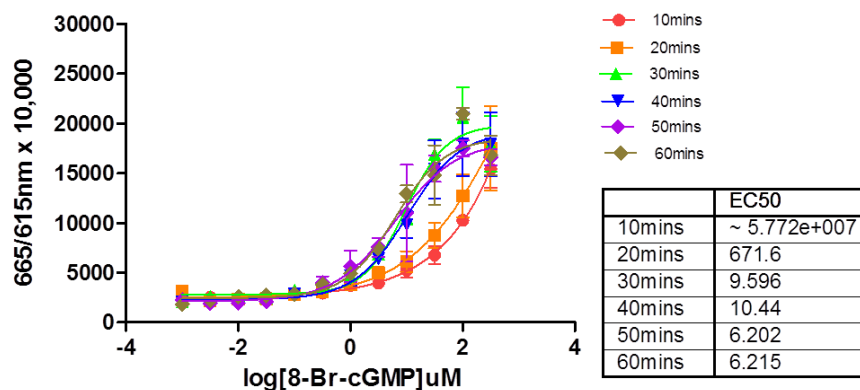
Phos-VASP Ser239
Hek vs. Hek VASP Stable cell line



Phos-VASP Ser239
Optimizing cell density



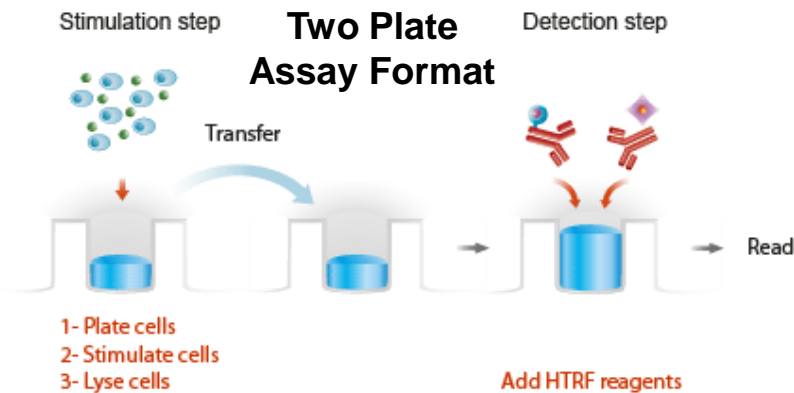
Phos-VASP Ser239
Optimizing Cell Treatment Time



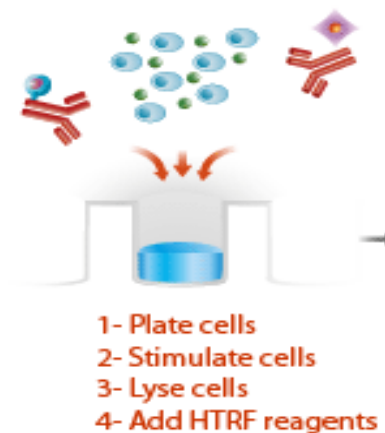
	EC50
10mins	~ 5.772e+007
20mins	671.6
30mins	9.596
40mins	10.44
50mins	6.202
60mins	6.215

Preferred Method: Hek-VASP cell line, 1250 cells/well, 30 min treatment

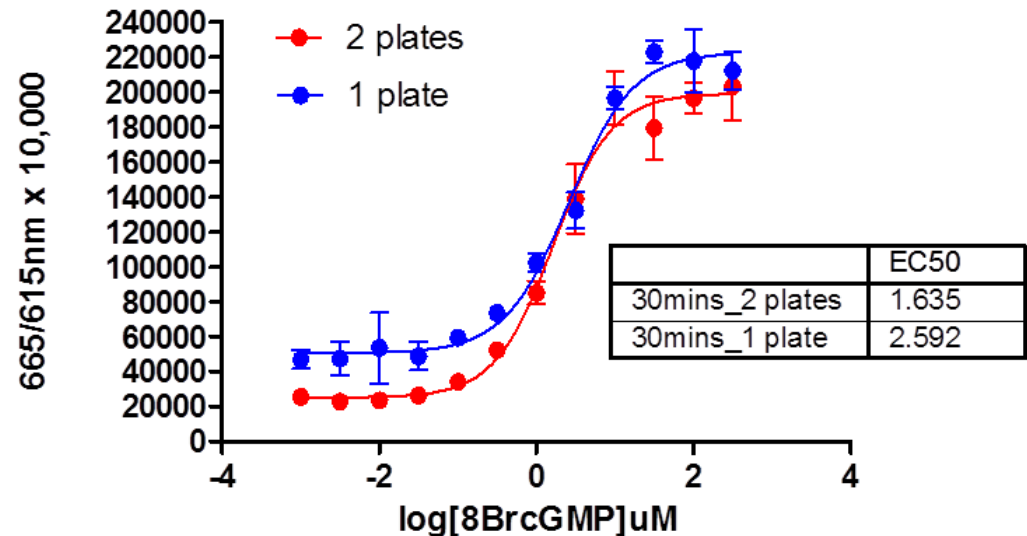
Two vs. One Plate Assay Format



One Plate Assay Format



Phos-Ser239 VASP: 2 vs. 1 Plate



- ▶ Cells were treated with 8Br-cGMP for 30 minutes
- ▶ Media removed and cells lysed at RT and then transferred to new second plate or
- ▶ Cells lysed directly in media and assay continued in same plate

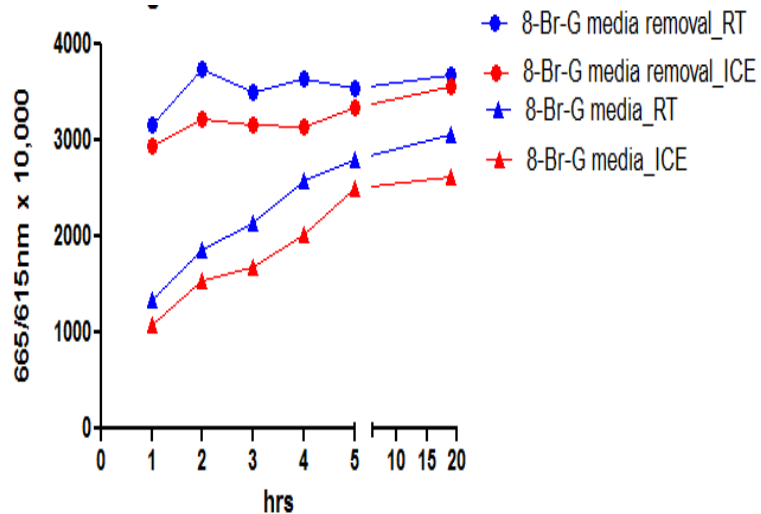
Preferred Method: One plate format

Optimizing Lysis Conditions

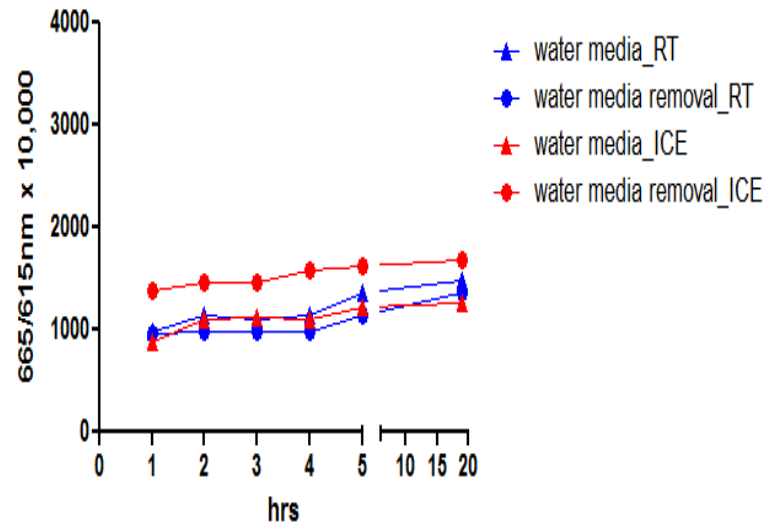
Exploring Signal Stability Over Time

Phos-Ser239 VASP

Cells treated with 8Br-cGMP



Cells treated with vehicle (water)



- ▶ Cells treated with 100 μ M 8Br-cGMP
- ▶ Cells lysed by adding lysis buffer directly into the media, or the media was removed and then added to the cells
 - ▶ Later switched to treat cells in HBSs instead of media and were able to get maximum signal by 2 hours
- ▶ Cells lysed either on ice or RT for 30 minutes
- ▶ Signal read over time between 1 and 20 hours

Preferred Method: Lyse cells at RT and read plate after 2 hours

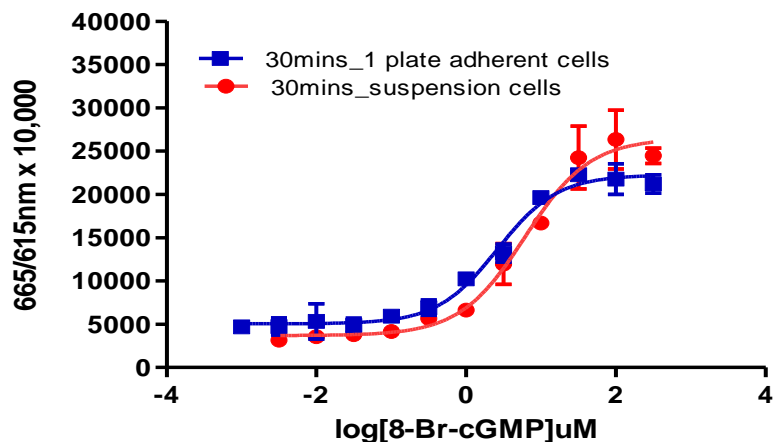
Assay Ready Frozen Cells in Suspension

DMSO Tolerability

Phos-Ser239 VASP

Cultured cells:

Plated day before or day of assay

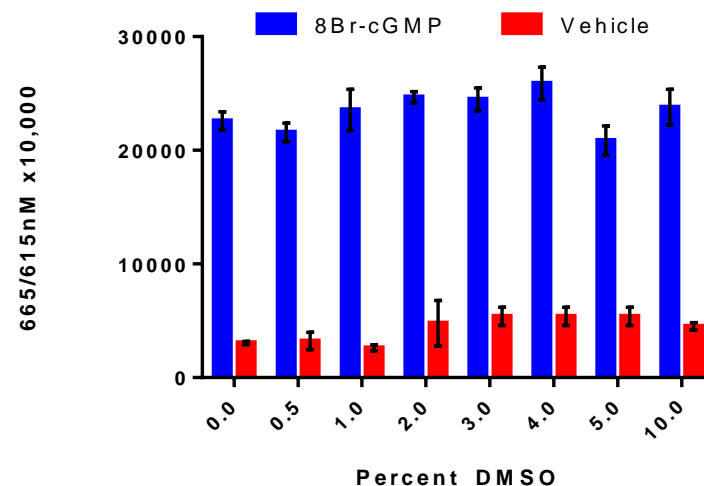


	EC50
30mins_suspension cells	6.099
30mins_1 plate adherent cells	2.612

- ▶ Cells were plated day before or day of assay
- ▶ Cells were treated with 8Br-cGMP for 30 minutes

Frozen cells in suspension:

DMSO tolerability

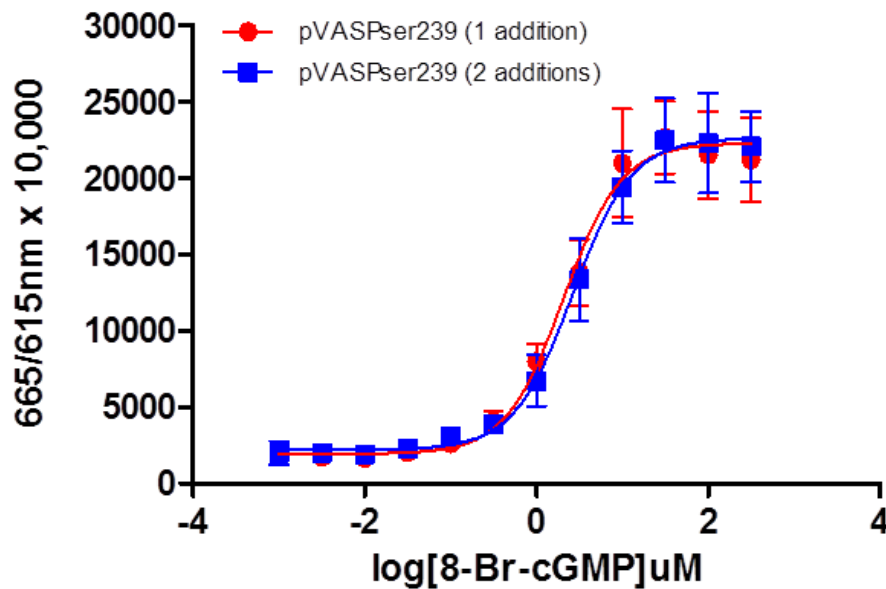


- ▶ Assay ready frozen cells in suspension plated day of assay
- ▶ Cells were treated with 100 uM 8Br-cGMP for 30 minutes

Preferred Method: Assay ready frozen cells in suspension with 1% DMSO

One vs. Two Additions for Antibody Reagents

Phos-Ser239 VASP: 1 vs. 2 step Antibody Addition

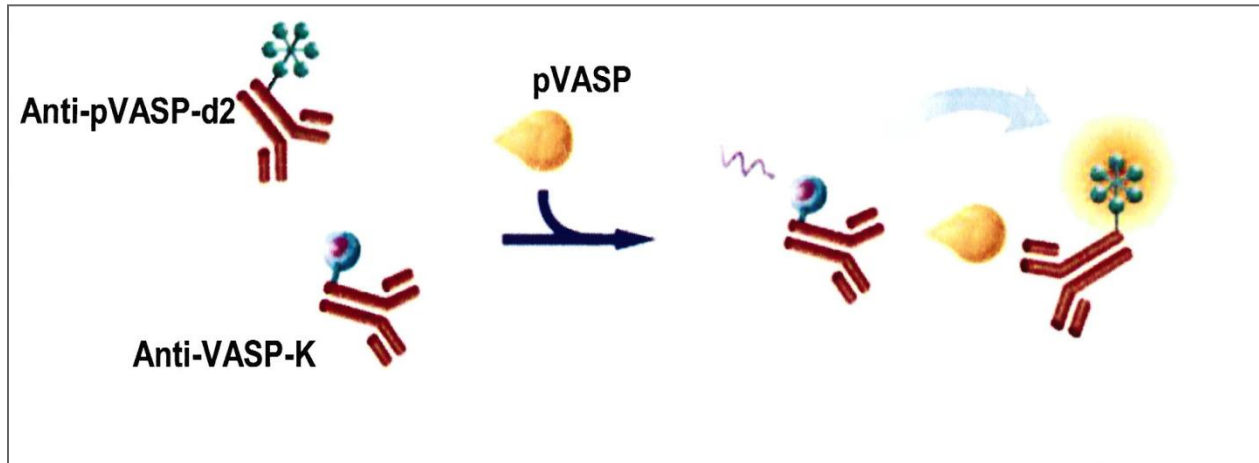


- ▶ Cells treated with 8Br-cGMP for 30 minutes;
- ▶ Antibodies added one at a time sequentially or pre-mixed and then added together

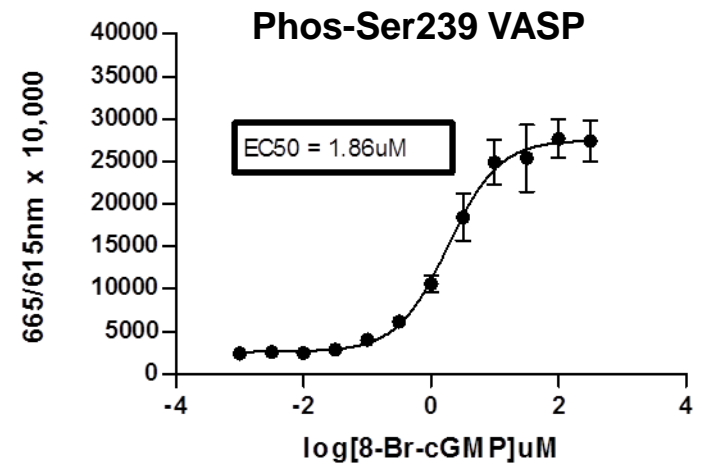
	EC50
pVASPser239 (1 addition)	2.071
pVASPser239 (2 additions)	2.657

Preferred Method: Pre-mix antibodies and one step addition

Phos-VASP Ser239 and Ser157 HTRF Assay

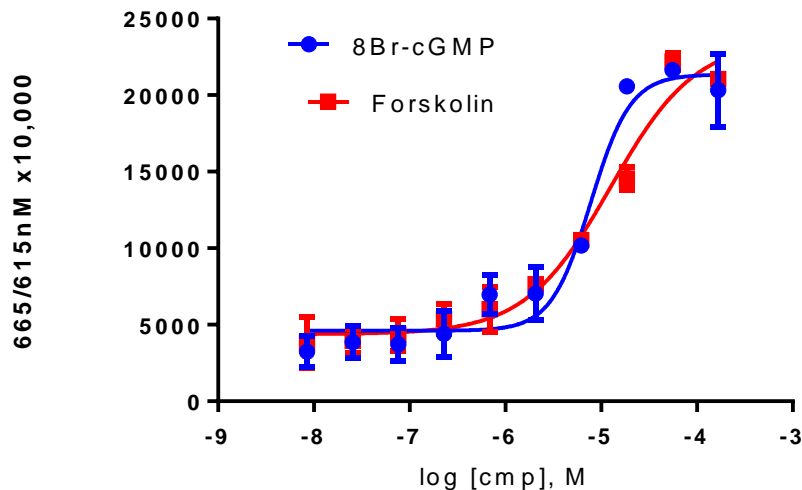


- Jump-In Hek293 or U2OS overexpressing VASP cell lines
- Assay ready frozen cells used in suspension day of assay
- 384 well low volume proxiplate
- 1250 cells per well in 6 ul HBSS
- Compound treatment in 1 % DMSO
- Lyse cells (2 ul) directly into HBSS at room temperature for 30 minutes
- Mix antibodies and add to plate (2 ul)
- Read two hours later, up to 20+ hours later
- ~0.7 Z-prime

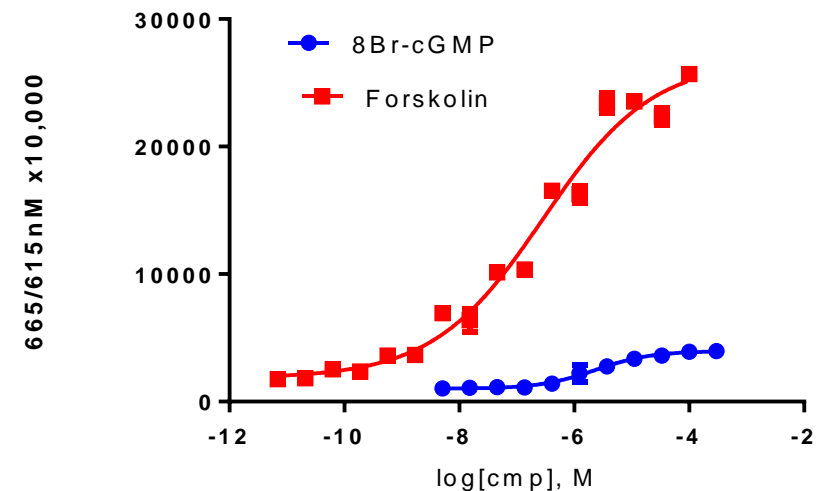


Recombinant Cell Line for Screening: Hek293 Cells Overexpressing VASP

Phos-Ser239 VASP



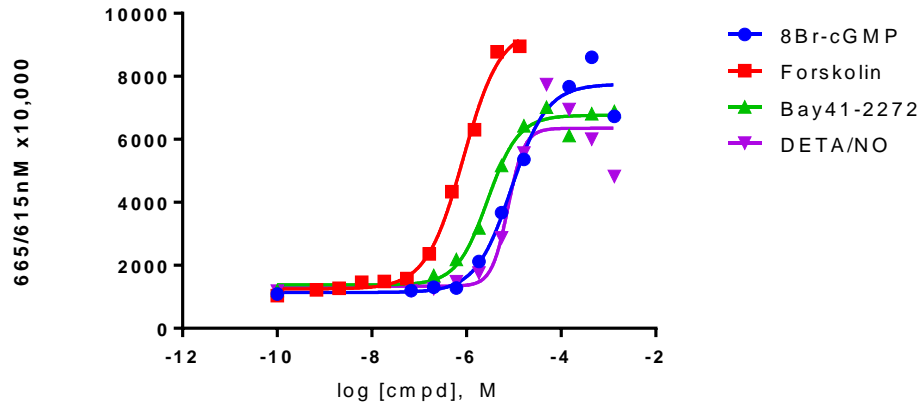
Phos-Ser157 VASP



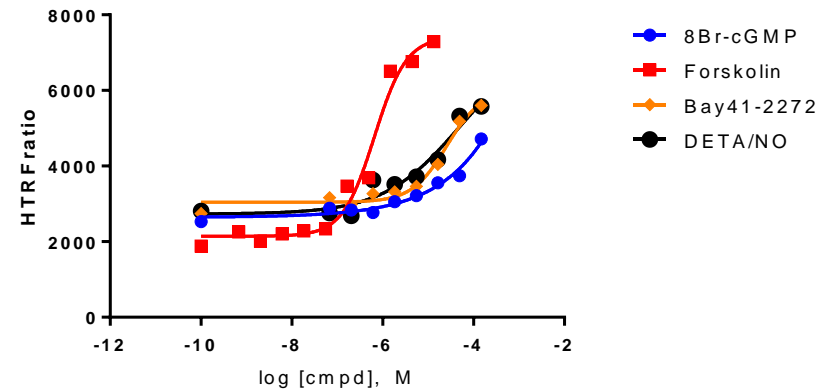
cAMP pathway leads to Ser239 and Ser157 VASP phosphorylation
cGMP pathway leads to Ser239 VASP phosphorylation

Physiological Relevant Cells: Rat Aortic Smooth Muscle Cells

Rat AoSMC- VASP Phos-Ser239

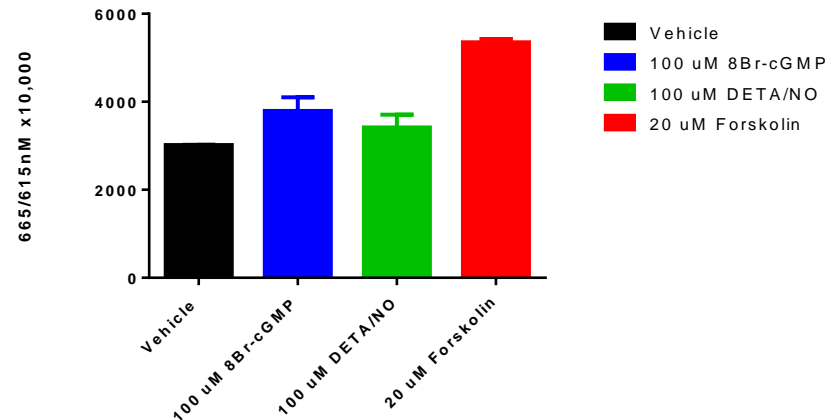


Rat AoSMC- VASP phos-Ser157



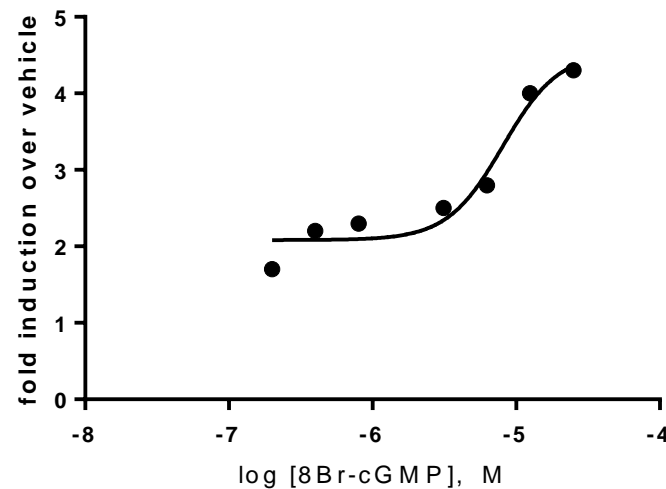
- cAMP pathway leads to Ser239 and Ser157 VASP phosphorylation and ~2-fold induction phos-CREB
- cGMP pathway leads to Ser239 VASP phosphorylation

Rat AoSMC- phos-CREB



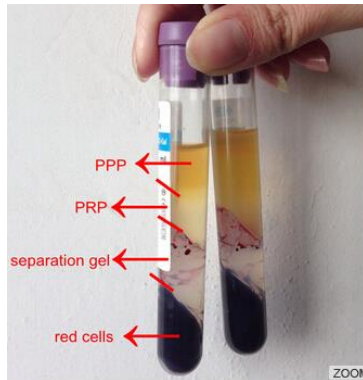
Mouse Whole Blood Ex Vivo Assay

Phos-Ser239 VASP: Ex Vivo Treatment



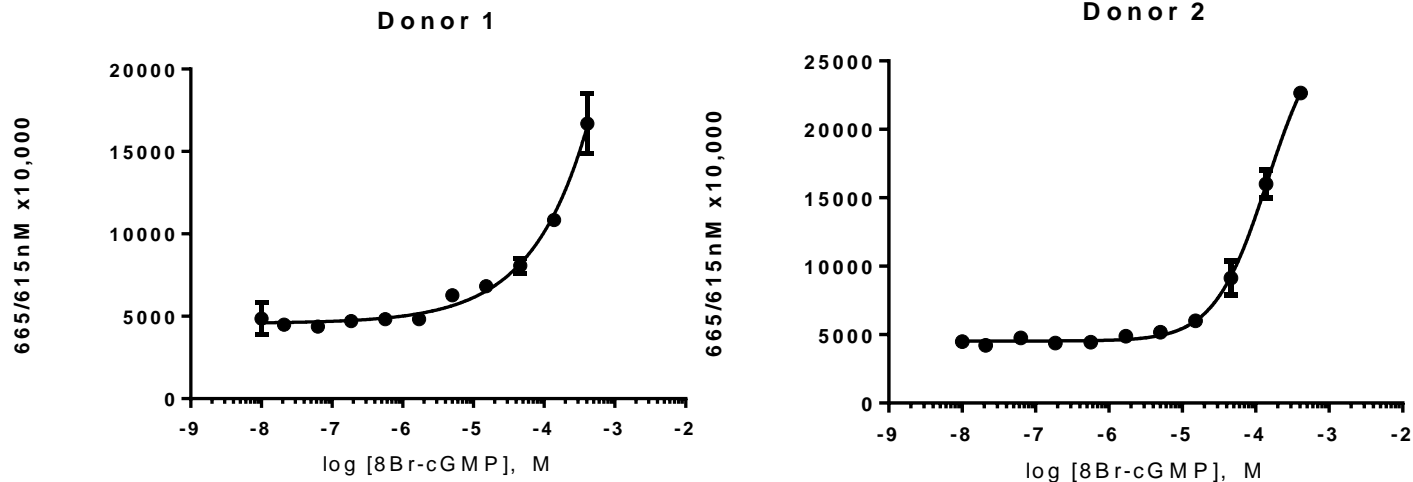
VASP HTRF can be used as target engagement assay
to help establish in vitro in vivo correlations

Human Platelet Ex Vivo Assay



- Human donor blood collected in citrate tubes
- Blood centrifuged to separate RBCs and plasma
- Plasma centrifuged in a second step
- Remove plasma and retain platelet pellet
- Platelets treated with Prostaglandin E1
- Platelets counted and plated in 384-well plate
- Platelets treated ex-vivo
- HTRF assay performed

Phos-Ser239 VASP: Ex Vivo Treatment



VASP HTRF can be used on human platelet samples for potential biomarker
Potential to look at PBMC and whole blood

Conclusion

- ▶ We developed VASP HTRF assays to measure phos-Ser239 and phos-Ser157 in cells.
 - 384 well format, 10 ul final volume, robust signal
- ▶ Assay can be used to differentiate compounds affecting cGMP and cAMP pathways
- ▶ Advantages of HTRF for Target Engagement
 - Assay can be run in high throughput in a streamlined fashion with few liquid additions, no liquid exchanges and with minimal steps
 - Robust assay in recombinant cells for routine screening
 - Assay translates to physiologically relevant cells (ex. Rat AoSMCs)
 - Assay can be performed on ex vivo and in vivo animal model samples and utilized to help establish in vitro in vivo correlations
 - Assay has potential as biomarker assay for human clinical samples