

# An Automated, High Throughput Method for the Quantification of Endogenous Cellular Kinase Activity

Brad Larson<sup>1</sup>, Nicolas Pierre<sup>2</sup>, Suzanne Graham<sup>2</sup>, Laurence Jacquemart<sup>2</sup>, Francois Degorce<sup>2</sup>, and Peter Banks<sup>1</sup>

<sup>1</sup>BioTek Instruments, Inc., Winooski, Vermont, USA • <sup>2</sup>Cisbio US, Inc., Bedford, Massachusetts, USA



## Introduction

Cellular kinases play an important role in relaying signals from activated receptors residing at the cell membrane to the interior of the cell, through signal transduction. The cellular processes in which they are associated include angiogenesis, cell growth, cell migration, and apoptosis. Overexpression of the kinase, or constitutive activity, has also been linked to a number of disease states, including vascular disease, bone disorders, and multiple forms of cancer. Therefore, cellular kinases continue to be an important target for small molecule and large molecule drug development. The ability to screen compounds or antibodies for their antagonistic effects on these signaling molecules, in a rapid and robust way, is critical to current efforts in this area.

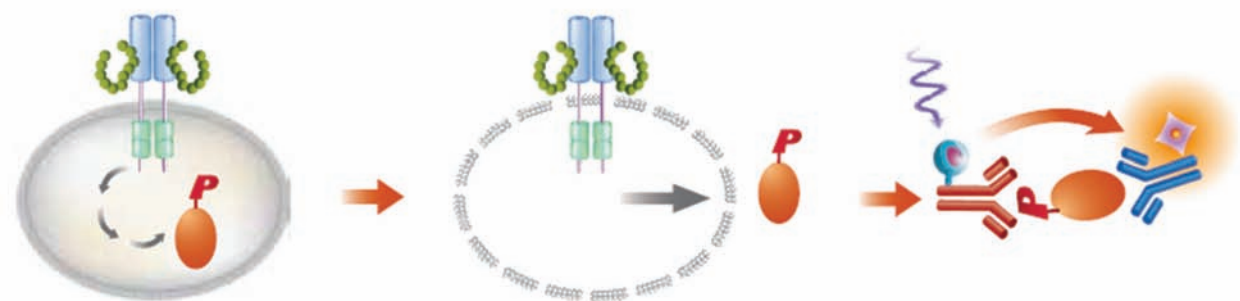
Here we describe a method using HTRF<sup>®</sup> technology to quantify cellular kinase activity without overexpression of kinase or its substrate. Two specific kinases were included in the study:

1. Vascular Endothelial Growth Factor Receptor 2 (VEGFR2), one of multiple subtypes of the receptor tyrosine kinase VEGFR, is involved in mediation of almost all cellular responses to VEGF.
2. Signal Transducer and Activator of Transcription 3 (STAT3), acts as a transcription activator, and is phosphorylated by receptor-associated kinases in response to cytokines and growth factors.

Detection of phosphorylated kinase involves a sandwich immunoassay which includes an anti-phospho specific antibody coupled with an anti-total antibody labeled with either donor or acceptor fluor. Activation of the appropriate signaling cascade causes an increase in HTRF signal, whereas inhibition will demonstrate the opposite effect.

Automation of all assay component dispense steps were performed in a high throughput format using a non-contact dispenser. Simultaneous, rapid detection of the fluorescent signal from the donor and acceptor molecules, in LV384-well format, was accomplished through the combination of a high performance Xenon flash lamp and dual matched PMTs incorporated into a novel HTS microplate reader. Optimization, validation, and pharmacology data demonstrate how the combination of assay chemistry, automation, and detection technologies create a robust method for high throughput screening of potential modulators of these important cell signaling pathways.

## HTRF Cellular Kinase Assay



**Figure 1** – The HTRF phospho-STAT3 (Tyr705) and phospho-VEGFR2 (Tyr1175) assays are designed for detecting and studying activated kinases directly in whole cells. Upon receptor activation, the kinases are activated, leading to kinase phosphorylation. After the lysis of the cell membrane, phosphorylated kinase can be detected upon the addition of two monoclonal antibodies: an anti-total antibody labeled with d2 and an anti-phospho-kinase antibody labeled with d1. In the presence of phosphorylated kinase, upon excitation of the Eu<sup>3+</sup>-cryptate, energy is transferred to the d2 molecule, and emission at 665 nm is seen. In the absence of the phosphorylated kinase, no energy is transferred, and no 665 nm signal is detected.

## Materials and Methods

### Materials

A431 cells (Catalog No. 85090402) were purchased from Sigma-Aldrich Corporation (St. Louis, MO). Dulbecco's Modified Eagle Medium (DMEM, Catalog No. 11995-065), Penicillin-Streptomycin, liquid (Catalog No. 15140-148), and Fetal Bovine Serum (FBS, Catalog No. 10437-028) were purchased from Life Technologies (Carlsbad, CA). Cryopreserved Human Umbilical Vein Endothelial Cells (HUVEC, Catalog No. CC-2517, Lot 152470), EBM<sup>®</sup> (Endothelial Base Medium, Catalog No. CC-3121), and EGM<sup>®</sup> SingleQuots<sup>®</sup> (Catalog No. CC-4133) were donated by Lonza (Walkersville, MD). Human Serum Albumin (HSA), 25% Solution (Catalog No. 1500233) was purchased from Baxter (Deerfield, IL). HTRF Phospho-STAT3 (Tyr705, Catalog No. 62AT3PEG) and Phospho-VEGFR2 (Tyr1175) assay kits were provided by Cisbio Bioassays (Codolet, FR). Hank's Balanced Salt Solution (HBSS) was purchased from Mediatech, Inc. (Manassas, VA). Recombinant Human Epidermal Growth Factor (EGF, Catalog No. cyt-217) and Recombinant Human Vascular Endothelial Growth Factor (VEGF, Catalog No. cyt-241) were purchased from ProSpec-Tany Technogene Ltd. (Ness Ziona, Israel). SD 1008 (Catalog No. 3035), Stattic (Catalog No. 2798), Cryptotanshinone (Catalog No. 3713), Ki 8751 (Catalog No. 2542), DMH4 (Catalog No. 4471), and Human VEGFR2 Antibody (Catalog No. AF357) were purchased from R&D Systems (Minneapolis, MN). 384 Well, Low Volume, Non-treated plates (Catalog No. 3674) were donated by Corning Life Sciences (Kennebunk, ME).

### Methods

**Cell Propagation:** A431 cells were propagated in complete medium, as described above. The cells were serum-starved 18 hours prior to running the assay by trypsinizing the cells from the flask, removing the complete medium, and replacing with serum-free medium (SFM) (DMEM plus 0.1% HSA). HUVECs were propagated in EBM plus EGM SingleQuots.

**Cell Plating:** Following the serum starvation period, A431 cells were removed from the flask and resuspended to a concentration of 2.5x10<sup>6</sup> cells/mL in HBSS buffer and plated in a 4 µL volume. HUVECs were trypsinized from the flask and resuspended to a concentration of 1.25x10<sup>6</sup> cells/mL in EBM. The cells were then plated in a 200 µL volume.

**Ligand/Compound Preparation:** EGF and VEGF were resuspended in Milli-Q water to a concentration of 100 µg/mL and then diluted appropriately in HBSS buffer (STAT3 assay) or EBM (VEGFR2 assay) respectively. Compounds were resuspended in 100% DMSO at a concentration of 10 mM, and then titrated in the same solvent. VEGFR2 Ab was resuspended in sterile PBS and titrated in EBM. The compounds and antibody were further diluted in HBSS buffer or medium as previously described to create the final assay concentrations.

## BioTek Instrumentation

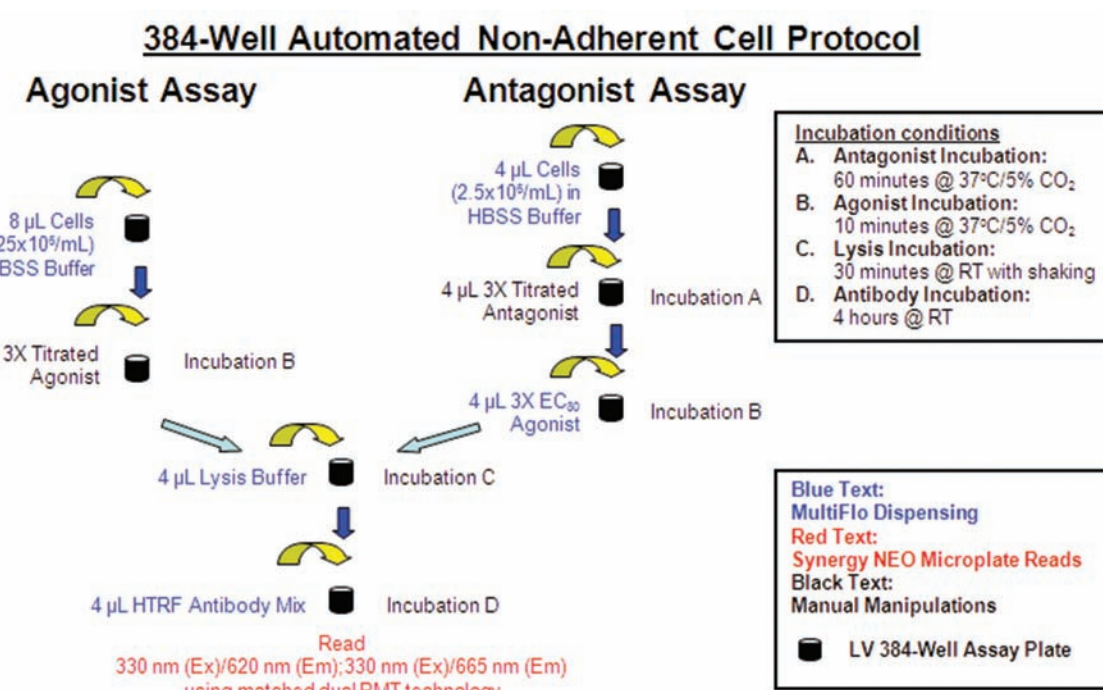
**MultiFlo<sup>™</sup> Microplate Dispenser.** The dispenser offers fast, accurate plate dispensing capabilities through its two peristaltic and two syringe pumps, with volumes ranging from 500 nL – 3000 µL. The instrument was used to dispense cells, agonist, lysis buffer, and antibody mixes in volumes as low as 4 µL.

**EL406<sup>™</sup> Combination Washer Dispenser.** The EL406 offers fast, accurate media removal and plate washing capabilities through its Dual-Action<sup>™</sup> Manifold. It also offers reagent dispensing capabilities through the use of its peristaltic or syringe pumps, with volumes ranging from 500 nL – 3000 µL/well. The instrument was used for cell dispensing, media removal and plate washing, as well as to dispense EC<sub>50</sub> agonist and HTRF assay components.

**Precision<sup>™</sup> Microplate Pipetting System.** The Precision combines an 8-channel pipetting head and an 8-channel bulk reagent dispenser in one instrument. The instrument was used to titrate and dilute inhibitors, as well as transfer diluted inhibitors to the 96-well cell plate, and a portion of the lysate to the LV384-well plate in the VEGFR2 assay procedure.

**Synergy<sup>™</sup> NEO Multi-Mode Microplate Reader.** The reader combines a filter-based and monochromator-based detection system in one unit. The HTRF certified reader uses a filter-based system, high performance xenon flash lamp, as well as dual matched PMTs to simultaneously detect the 665 nm and 620 nm fluorescent emissions from the assay chemistry.

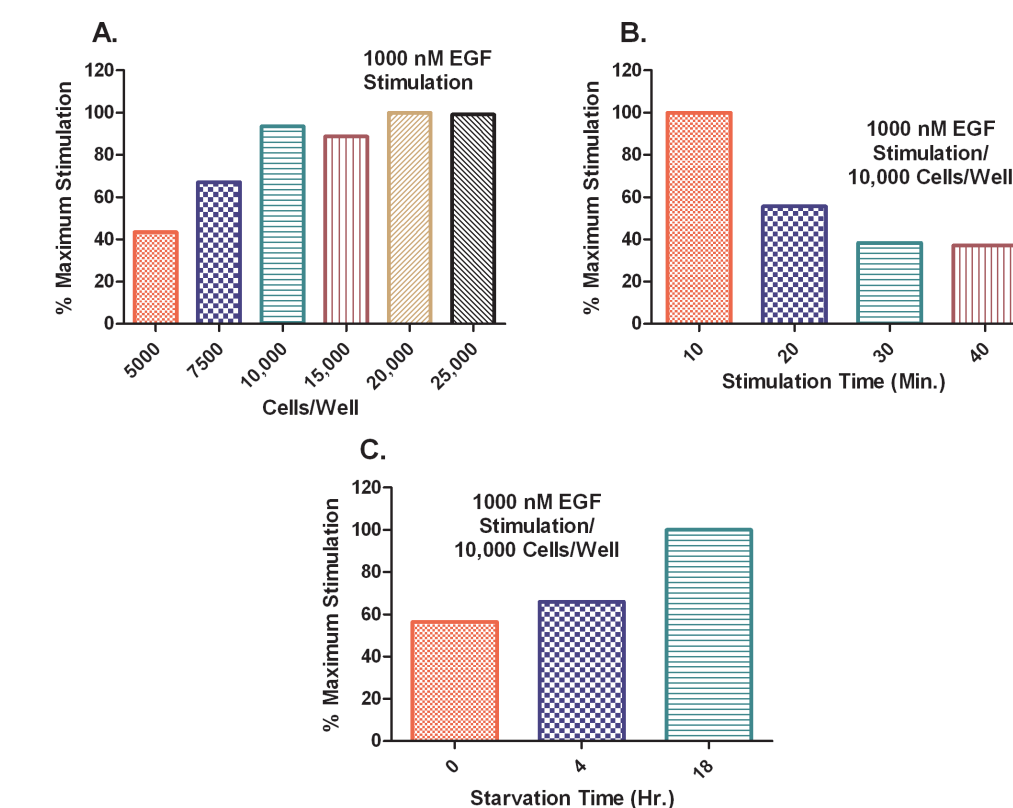
## Phospho-STAT3 (Tyr705) Assay



**Figure 2** – Automated phospho-STAT3 assay workflow. HTRF signal detection was carried out using the Xenon Flash Lamp of the Synergy NEO with the following settings: Delay after plate movement: 0 msec; Delay after excitation: 150 µsec; Integration time: 500 µsec; Read height: 8.0 mm.

### Assay Optimization

Optimization of assay parameters is essential to ensure the generation of robust, reliable experimental data. When monitoring the modulation of kinase signaling pathways, typical parameters include cell concentration, agonist stimulation time, and serum starvation length. Each condition tested was analyzed for its effect on the difference in signal between stimulated and unstimulated cells (assay window).

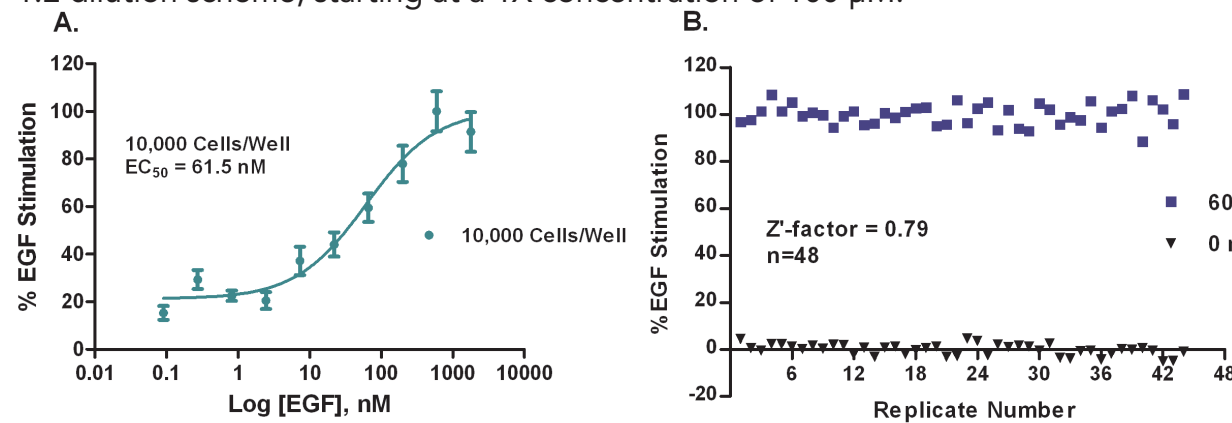


**Figure 3** – phospho-STAT3 Assay optimization results. A. Cell concentration analysis (5,000-25,000 cells/well tested). B. Signaling pathway stimulation (10-40 minute stimulations). C. Serum starvation (0, 4 hour, or overnight starvations).

The results shown in Figure 3A demonstrate that cell concentrations above 10,000 cells/well yield no appreciable increase in assay window. Therefore, the 10,000 cells/well concentration was chosen to lower cell propagation time and cost. Furthermore, Figure 3B illustrates that rapid stimulation of the signaling pathway is seen at 37°C, following agonist addition. A 10 minute post agonist incubation time was then used for subsequent experiments. Finally, it is apparent from the data shown in Figure 3C that serum starving the A431 cells prior to assay performance serves to lower the basal level of STAT3 signaling pathway activity, thereby creating a larger assay window between stimulated and unstimulated cells.

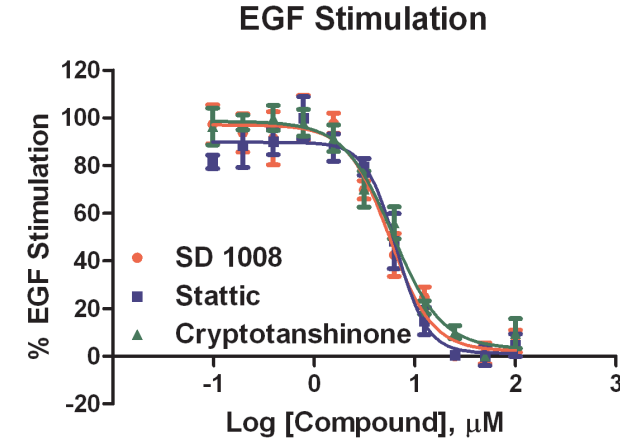
### Assay Validation

Z'-factor<sup>1</sup>, agonist, and antagonist titration experiments were performed to validate the optimized phospho-STAT3 automated assay procedure. Serial 1:3 EGF titrations were created in HBSS buffer. Concentrations tested ranged from 5000-0 nM. The EGF concentration yielding the largest EGF stimulation, 600 nM, was then used as the positive control, along with 0 nM as the negative control, in a Z'-factor experiment to measure assay robustness. The Z'-factor value takes into account the difference in signal between a positive and negative control, as well as the variation in the signal amongst replicates. A scale of 0-1 is incorporated, with values ≥0.5 being indicative of an excellent assay. Finally, a pharmacology validation was performed using the three known STAT3 pathway inhibitors, SD 1008, Stattic, and Cryptotanshinone. 11-point titrations were created using a 1:2 dilution scheme, starting at a 1X concentration of 100 µM.



**Figure 4** – EGF Titration and Z'-factor results. A. Data shown represents the average and standard deviation of four replicates tested at each EGF concentration. B. Assay run in agonist mode as previously described. Forty-four replicates of 600 nM and 0 nM EGF used as positive and negative control, respectively.

### Inhibition of EGF Stimulation



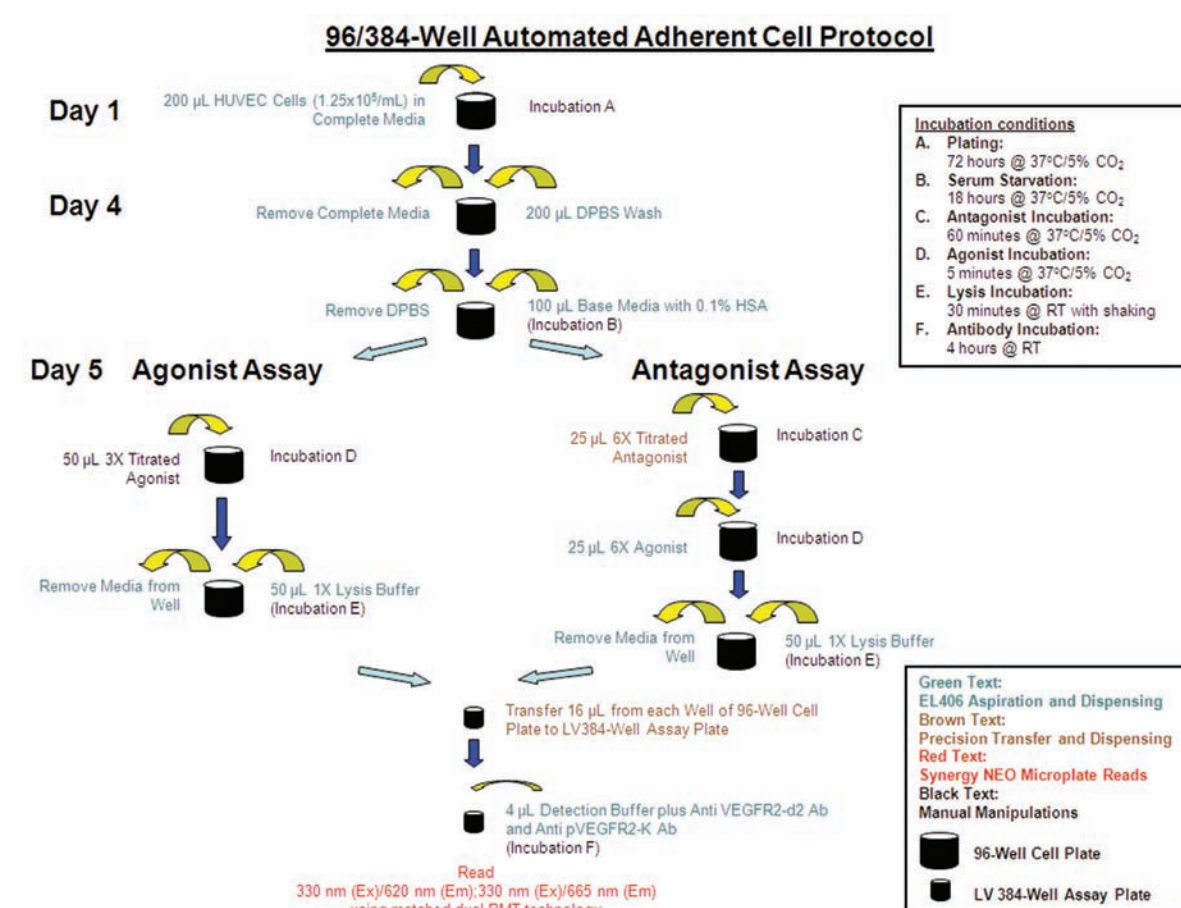
**Figure 5** – Compound inhibition curves. Percent EGF stimulation plotted for each concentration of inhibitor tested. Data shown represents average and standard deviation for n=4 replicates.

Compound	Generated IC <sub>50</sub> Value	Literature IC <sub>50</sub> Value
SD 1008	5.6	Inhibition seen at 10 µM in as little as 30 minutes <sup>2</sup>
Stattic	6.6	5.1 <sup>3</sup>
Cryptotanshinone	6.1	4.6 <sup>4</sup>

A Z' value of 0.79 was generated, which as explained previously, is indicative of an excellent, robust assay. The EC<sub>50</sub> value of 61.5 nM for EGF stimulation of the STAT3 signaling pathway in A431 cells agrees with the value previously generated by the assay manufacturer of 180 nM. An EC<sub>80</sub> value of approximately 200 nM was then used for subsequent antagonist testing.

Results from the antagonist test were in close agreement to published literature values (Table 1), demonstrating the ability of the automated assay to produce accurate results when testing other unknown STAT3 pathway inhibitors.

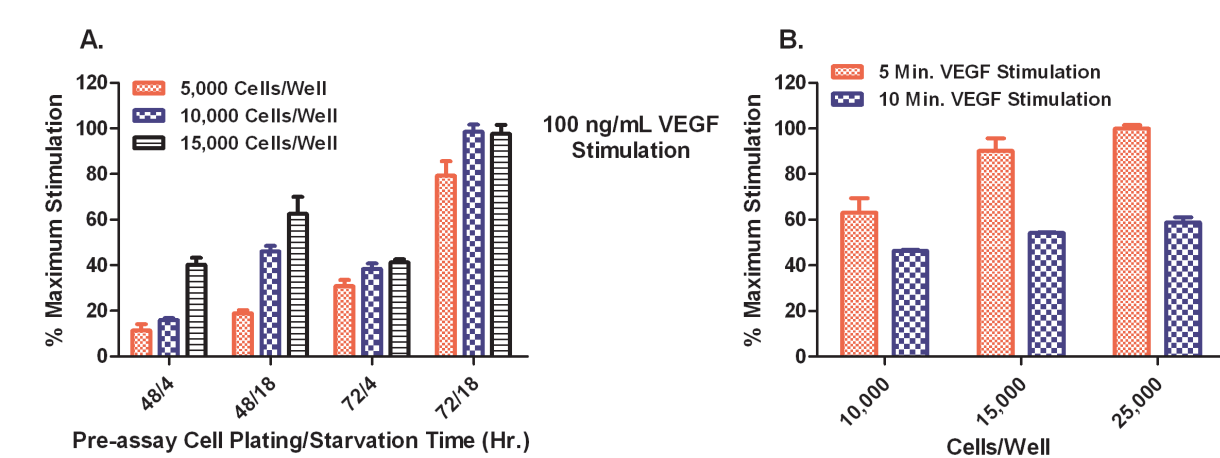
## Phospho-VEGFR2 (Tyr1175) Assay



**Figure 6** – Automated phospho-VEGFR2 assay workflow. HTRF signal detection settings on the Synergy NEO were as previously described.

### Assay Optimization

Optimization of assay parameters was once again performed for the phospho-VEGFR2 assay. Cell concentration, plating time, serum starvation, and stimulation time were again analyzed to determine their effect on assay window. Signaling pathway stimulation was carried out using 100 ng/mL VEGF.

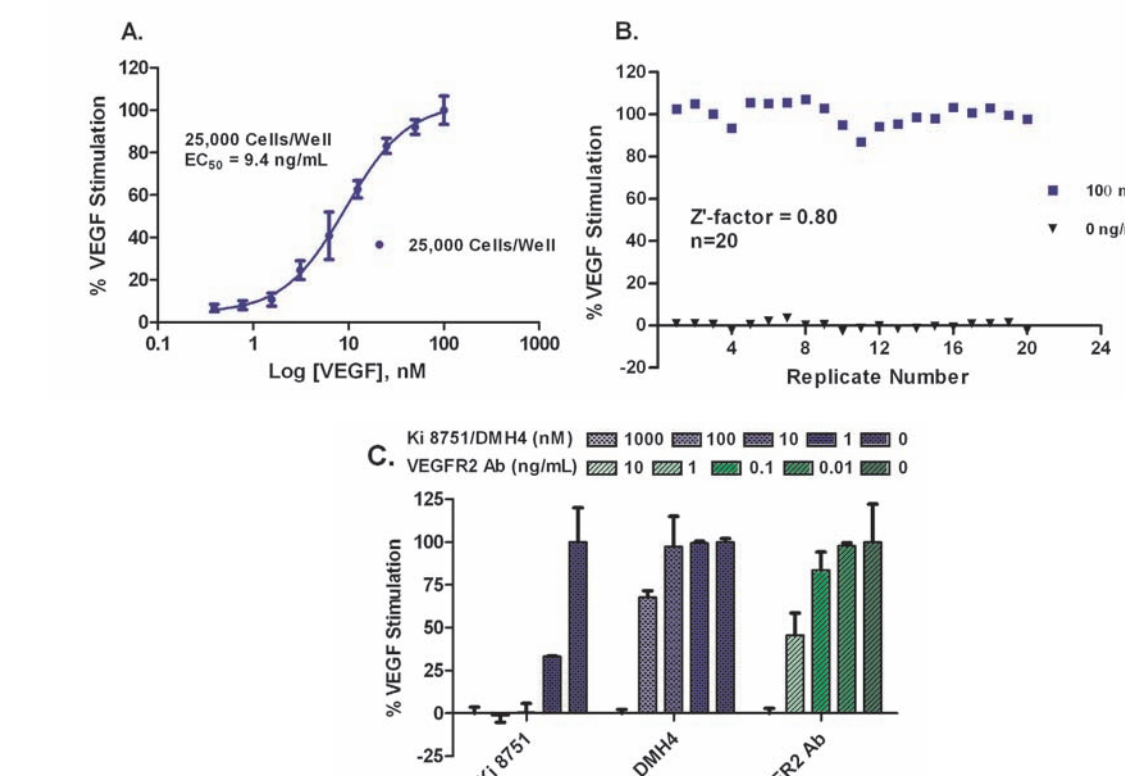


**Figure 7** – phospho-VEGFR2 Assay optimization results. A. Cell plating and serum starvation time analysis (2 and 3 day plating/4 hour and overnight starvations). B. Cell concentration and stimulation time test (5 and 10 minute stimulations/10,000-25,000 cells/well).

The data in Figure 7A demonstrates that allowing the cells to recover from plating and propagating for a period of three days, followed by an overnight serum starvation provides the largest assay window for each of the three cell concentrations tested (5,000-15,000 cells/well). Further optimization verified that higher cell concentrations, in addition to a minimized stimulation period, provide larger differences between stimulated and unstimulated cells. The final conditions chosen included a 3 day plating and overnight (18 hour) serum starvation of 25,000 cells/well, followed by a 5 minute VEGF stimulation time.

### Assay Validation

Z'-factor, agonist, and antagonist titration validation experiments were performed as previously described with the automated phospho-VEGFR2 assay procedure. Serial 1:2 VEGF titrations were created in SFM. Concentrations tested ranged from 100-0 ng/mL. 100 and 0 ng/mL VEGF were also used as the positive and negative controls, respectively, in the Z'-factor experiment. Finally, pharmacology validations were once again performed using two known VEGFR2 small molecule inhibitors, Ki 8751 and DMH4, as well as a VEGFR2 neutralizing antibody. 5-point titrations were created using a 1:10 dilution scheme, starting at a 1X concentration of 1000 nM (small molecule) or 10 µg/mL (antibody).



Ligand/Inhibitor	Observed Value	Literature Value
Human VEGF (EC <sub>50</sub> )	9.4 ng/mL	1-6 ng/mL <sup>3</sup>
Ki 8751 (IC <sub>50</sub> )	< 1 nM	0.9 nM <sup>6</sup>
DMH4 (IC <sub>50</sub> )	100<IC <sub>50</sub> <1000 nM	161 nM <sup>7</sup>
VEGFR2 Antibody (IC <sub>50</sub> )	0.1<IC <sub>50</sub> <1 µg/mL	0.05-0.25 µg/mL <sup>8</sup>

**Figure 8** – Assay validation results. A. Data shown represents the average and standard deviation of two replicates tested at each VEGF concentration. B. Assay run in agonist mode as previously described. Twenty replicates of 100 and 0 ng/mL VEGF used as positive and negative control, respectively. C. Percent VEGF stimulation plotted for each concentration of inhibitor tested. Data shown represents average and standard deviation for n=2 replicates.

The Z' value of 0.80 once again demonstrates the robustness of the automated process. VEGFR2 ligand and inhibitor pharmacology (Table 2) also illustrate the ability to generate appropriate EC<sub>50</sub> or IC<sub>50</sub> values using the optimized procedure.

## Conclusions

- The HTRF phospho-STAT3 and phospho-VEGFR2 assays afford sensitive, functional cell-based formats for the assessment of kinase signaling pathway modulation
- Each assay procedure can be easily automated in 384-well or 96-well format using the aspiration and non-contact dispensing capabilities of the EL406 or MultiFlo
- The xenon-based excitation and filter-based detection system of the Synergy NEO microplate reader is able to simultaneously quantify the emitted signals from the donor and acceptor fluor
- The combination of assay chemistry, as well as liquid handling and detection instrumentation, create an ideal solution for analysis of STAT3 and VEGFR2 pathway activators, in addition to small and large molecule inhibitors

The authors would like to thank Mark Rothenberg and Corning Life Sciences for the generous donation of 384-well low volume microplates, as well as Stephanie Nickles and Lonza Walkersville for their donation of Human Umbilical Vein Endothelial Cells and Medium. Both contributed greatly to the success of the project.

<sup>1</sup>Zhang, et al.: A Simple Statistical Parameter for Use in Evaluation and Validation of High Throughput Screening Assays. *J Biomol Screen* 1999 4(2): 67-73. | <sup>2</sup>Duan, et al.: 8-Benyl-4-oxo-8-azabicyclo[3.2.1]oct-2-ene-6,7-dicarboxylic Acid (SD-1008), a Novel Janus Kinase 2 Inhibitor, Increases Chemotherapy Sensitivity in Human Ovarian Cancer Cells. *Mol Pharmacol* 2007 72(5): 1137-45. | <sup>3</sup>Schust, et al.: Stattic: A Small-Molecule Inhibitor of STAT3 Activation and Dimerization. *Chem Biol* 2006 13(11): 1235-42. | <sup>4</sup>Shin, et al.: Cryptotanshinone Inhibits Constitutive Signal Transducer and Activator of Transcription 3 Function through Blocking the Dimerization in DU145 Prostate Cancer Cells. *Cancer Res* 2009 69(11): 193-202. | <sup>5</sup>Conn, et al.: Purification of a glycoprotein vascular endothelial cell mitogen from a rat glioma-derived cell line. *Proc Natl Acad Sci U S A*. 1990 87(4): 1323-1327. | <sup>6</sup>Kubo, et al.: Novel potent orally active selective VEGFR-2 tyrosine kinase inhibitors: synthesis, structure-activity relationships, and antitumor activities of N-phenyl-N'-[4-(4-quinolyl)oxy]phenylureas. *J Med Chem* 2005 48(5): 1359-1366. | <sup>7</sup>Hao, et al.: In vivo structure-activity relationship study of dorsomorphin analogues identifies selective VEGF and BMP inhibitors. *ACS Chem Biol* 2010 5(2): 245-253. | <sup>8</sup>Ferrara, et al.: The Biology of Vascular Endothelial Growth Factor. *Endocr Rev* 1997 18(1): 4-25. HTRF is a registered trademark of Cisbio Bioassays.