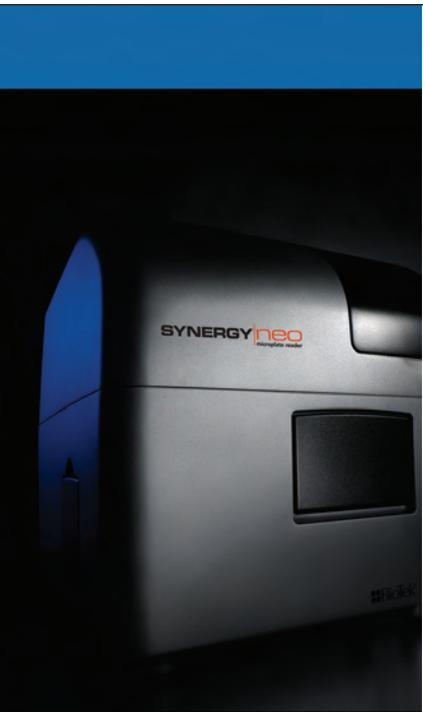


High Throughput Screening Applications



Introduction

High throughput screening (HTS), or the process by which libraries of small molecule compounds are individually assessed for binding, activating or inactivating biological activity in drug target molecules, has been part of the drug discovery process for more than two decades. Its primary purpose is to rapidly select which compounds possess the desired activity and thereby undergo further testing and optimization. Since its inception, HTS methods have undergone significant change: the early dependence on “home-brewed” assay chemistries and basic research instruments evolved into industrialized methods incorporating millions of compounds that could be screened against purified drug targets using sophisticated liquid handling, detection and robotics.

The HTS market remains strong, accounting for 35% of the global drug discovery technologies market, and is projected to reach \$19.9 billion by 2017¹. However, the industry is pressured to reduce the cost of the drug discovery process, and improve the number of true lead compounds coming out of screening campaigns; eventually becoming a marketed drug. This pressure has caused a shift in how HTS is conducted across the industry. Today, HTS methods have moved away from the “industrialized” emphasis, and adopted a “smarter screening” precept. Screening campaigns typically no longer involve millions of compounds, but use focused libraries that offer not only diversity but also known pharmacophores of the gene family of interest. Biochemical assays using purified drug targets are being supplanted by functional and phenotypic assays involving cells. Some of the phenotypic assays used today involve the kinetic monitoring of cellular response in real time. In addition, immortalized cell lines with over expressed drug targets are in turn being supplanted by the use of human primary cells with endogenously-expressed drug targets. Sourcing issues of sufficient quantities of human primary cells for HTS campaigns are beginning to be addressed by the use of induced pluripotent stem cells².

Smarter screening also extends to the instrumentation used in HTS. The liquid handling and detection instrumentation have to be able to perform a number of high level tasks from biochemical and cell-based assay procedures, to deliver accurate, efficient, and repeatable results using a diverse repertoire of detection modes. BioTek Instruments understands the evolution of HTS and has designed instrumentation that facilitates the smarter screens run in today’s HTS labs. In the following sections we describe key BioTek liquid handling and detection instrumentation for HTS methods and provide four HTS case studies. The case studies use commercially available compound libraries and assay chemistries suitable for nuclear receptor, GPCR, cellular kinase, as well as epigenetic drug targets.

BioTek Instrumentation

MultiFlo™ Microplate Dispenser

The MultiFlo™ Microplate Dispenser combines “multiple” dispensers in one compact unit, thus reducing overall instrument costs, robotic footprint, scheduling complexity and processing time. Its unique parallel dispense technology combines up to two eight-tip, non-contact peristaltic dispensers and two variable-tube syringe pump dispense manifolds, to maximize assay flexibility in 96-, 384-, and 1536-well microplates. For use with cell-based assays, the MultiFlo’s dispense lines and manifolds may be sterilized via chemical means or autoclaving. The instrument may also be placed in a laminar flow hood when used for sterile dispensing in standalone mode. In the assays detailed here, MultiFlo was used to dispense cells and medium, other assay components, and detection reagents to assay wells in 384- and low volume 384-well (LV384) format.

EL406™ Microplate Washer Dispenser

The EL406™ Microplate Washer Dispenser combines fast, full microplate washing along with a single eight-tip, non-contact peristaltic dispenser and two variable-tube syringe pump dispense manifolds in one compact unit. Therefore, valuable bench space is saved and, by reducing the expense and maintenance of multiple dedicated instruments, plate scheduling routines are simplified and processes are run faster. The instrument is also ideal for use with cell-based assays, affording the same sterilization capabilities and small footprint as the MultiFlo. In addition, the aspiration capabilities of the wash manifold, combined with any of the three dispensers, provide for fast, accurate medium exchanges and buffer washes. In the assays detailed here, the EL406 was used to dispense cells and medium, aspirate medium and perform buffer washes, and dispense detection reagents to assay wells in 96- and 384-well formats.

Precision™ Microplate Pipetting System

The Precision™ Microplate Pipetting System combines an 8-channel pipetting head and an 8-channel bulk reagent dispenser in one instrument. The use of sterile, disposable tips, combined with a small footprint, allows Precision to be inserted into biological safety cabinets and used with cell-based assays to perform routine procedures such as dilutions or serial titrations. For the projects described here, the instrument was used to dilute library compounds, create compound dose response curves for validation and secondary screening, and transfer all compounds to the 96- or 384-well assay plates.

Synergy™ Multi-Mode Microplate Readers

BioTek's Synergy™ Multi-Mode Microplate Reader family is ideally suited to the demands of today's screening laboratories. For applications such as high throughput screening of small molecules, the Synergy Neo is the optimal choice. For applications such as smaller small molecule library screens or biotherapeutics drug discovery, Synergy H1 is a more cost-effective choice. Both microplate readers use filter cubes and minimal optical components (no fiber optics) to ensure high light transmission and high sensitivity. They also incorporate BioTek's patented Hybrid Technology™, offering the flexible choice of detection with a monochromator system or a filter/dichroic mirror system.

Synergy™ Neo

Synergy Neo was designed specifically for ultra high performance detection from microplates required by today's HTS labs. For dual emission assays, such as time-resolved fluorescence resonance energy transfer (TR-FRET), the unit uses a filter cube design that allows simultaneous measurement of donor and acceptor fluorescence using dual photomultiplier tubes (PMTs), one which is a red-shifted PMT with high sensitivity for far red spectral emissions (Figure 1A). For luminescence, one side of the lower excitation cube is empty to allow transmission of the luminescent signal. The top emission cube contains a dichroic mirror to channel the signal to the low noise PMT (Figure 1B). Configurations for the TR-FRET and luminescent assays can be accomplished using the same excitation and emission filter cubes. The optical path is devoid of components that limit light transmission, such as fiber optic cables, therefore enabling high sensitivity.

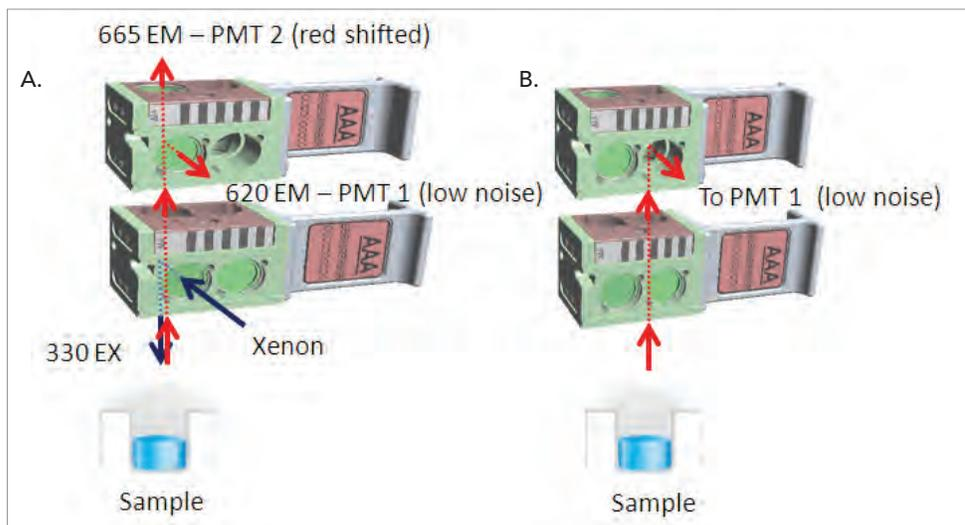


Figure 1. Schematic of the dual filter cubes in Synergy Neo used for the simultaneous detection of donor and acceptor fluorescence from (A) TR-FRET or (B) luminescent assays.

Synergy™ H1

Synergy H1 was designed to provide high sensitivity using a single filter block and PMT. Sequential detection of donor and acceptor fluorescence (Figure 2A) is therefore incorporated for TR-FRET assays, while luminescent signal detection is as previously described. Similar to the Synergy Neo design, the optical path (Figure 2B) is devoid of components that limit light transmission, such as fiber optic cables, therefore enabling high sensitivity.

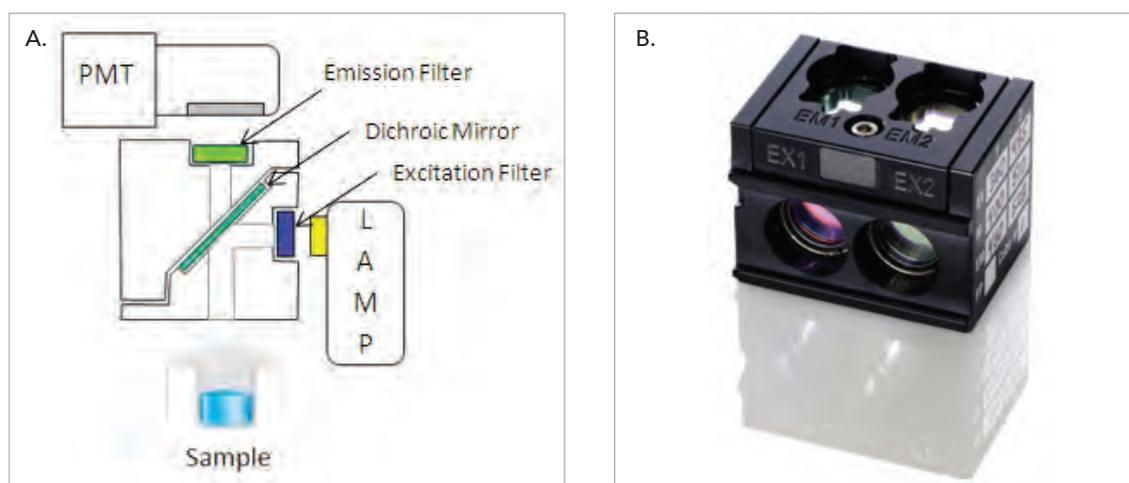


Figure 2. (A) Exploded schematic of the Synergy H1 filter cube used for sequential detection of donor and acceptor fluorescence. (B) Synergy H1 filter cube. Filters and dichroic mirror removed from position 2 for luminescence detection.

Application Overview

The following biochemical or cell-based screening applications analyze the effect of small molecules on a variety of current popular drug discovery targets. The 76 compound Screen-Well® Nuclear Receptor Ligand Library (Cat. No. BML-2802), 502 compound Screen-Well Natural Product Library (Cat. No. BML-2865), or 43 compound Screen-Well Epigenetics Library (Cat. No. BML-2836), donated by Enzo Life Sciences, Inc. (Farmingdale, NY), were used in each application. The assay chemistries incorporated either HTRF® (Cisbio Bioassays, Codolet, France) or luminescent detection technologies. Each screen was automated using the liquid handling instruments previously referenced. Quantification of the emitted signals was accomplished using the appropriate Synergy microplate reader.

Nuclear Receptors

Nuclear receptors make up a super family of signal mediating, intra-cellular receptors that are found in multiple species including insects and vertebrates such as humans. Being transcriptional regulators, they function as metabolic ligand activated transcription factors, regulating a number of functions, including homeostasis, development, and metabolism, and also play a role in diseases such as cancer and diabetes. Additionally, they are shown to bind small molecules modified by drug design³. This combination makes nuclear receptors popular pharmacological targets. Approximately 13% of drugs approved for sale in the United States are nuclear receptors with 15 of these drugs in the top 200 prescribed medicines. These top drugs represented \$27.5 billion of sales revenue in 2009⁴.

Androgen Receptor Agonist/Antagonist Studies using a Cell-based Nuclear Receptor Assay

The nuclear receptor assay system from INDIGO Biosciences (State College, PA) utilizes proprietary, non-human mammalian cells engineered to provide constitutive, high-level expression of a full length, unmodified nuclear receptor of choice. INDIGO's reporter cells include the luciferase reporter gene functionally linked to a bonafide nuclear receptor-responsive promoter. The readout from INDIGO's reporter cells demands the same orchestration of all intracellular molecular interactions and events that can be expected to occur in vivo. Thus, quantifying changes in luciferase expression in the treated reporter cells provides a sensitive surrogate measure of the changes in nuclear receptor activity.

The principle assay application is screening test samples to quantify the functional activities, either agonist or antagonist, that they may exert against the nuclear receptor. In the project described here, luciferase gene expression occurs after ligand-bound Androgen Receptor (AR) undergoes nuclear translocation, DNA binding, recruitment and assembly of the co-activators and accessory factors required to form a functional transcription complex, culminating in expression of the target gene.

A Z'-factor⁵ test was performed to validate the automated 384-well AR assay (Figure 3A). 6 α -Fluoro-testosterone (6 α -FIT) was used as the control compound. Twenty-four replicates of 10 μ M or 0 μ M compound were used as the positive and negative control, respectively. Dose response curves were also created using the control agonist, 6 α -FIT, and the control antagonist, Bicalutamide (BIC), to validate the ability of the automated assay to generate correct compound pharmacology (Figure 3B).

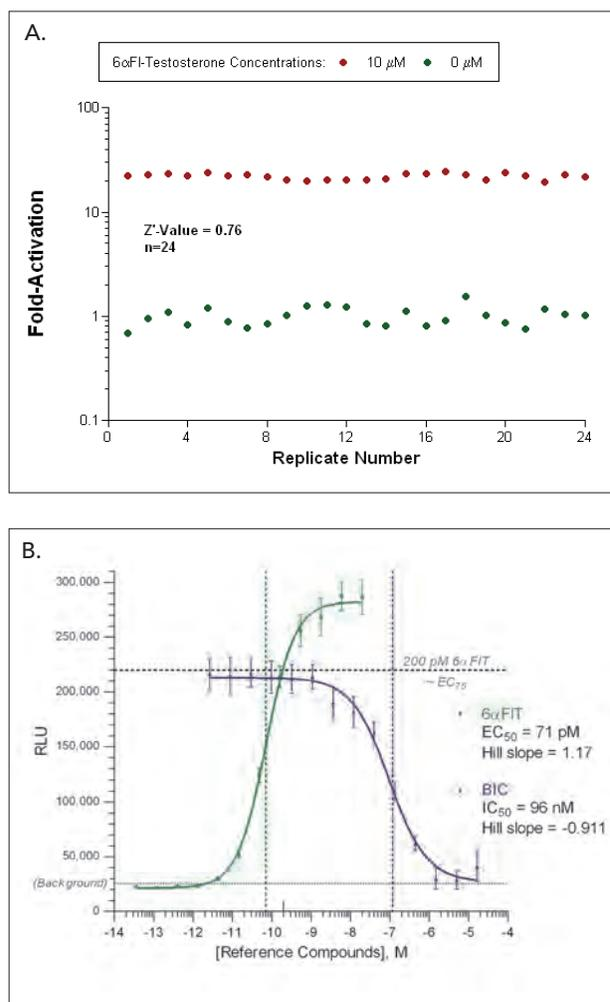
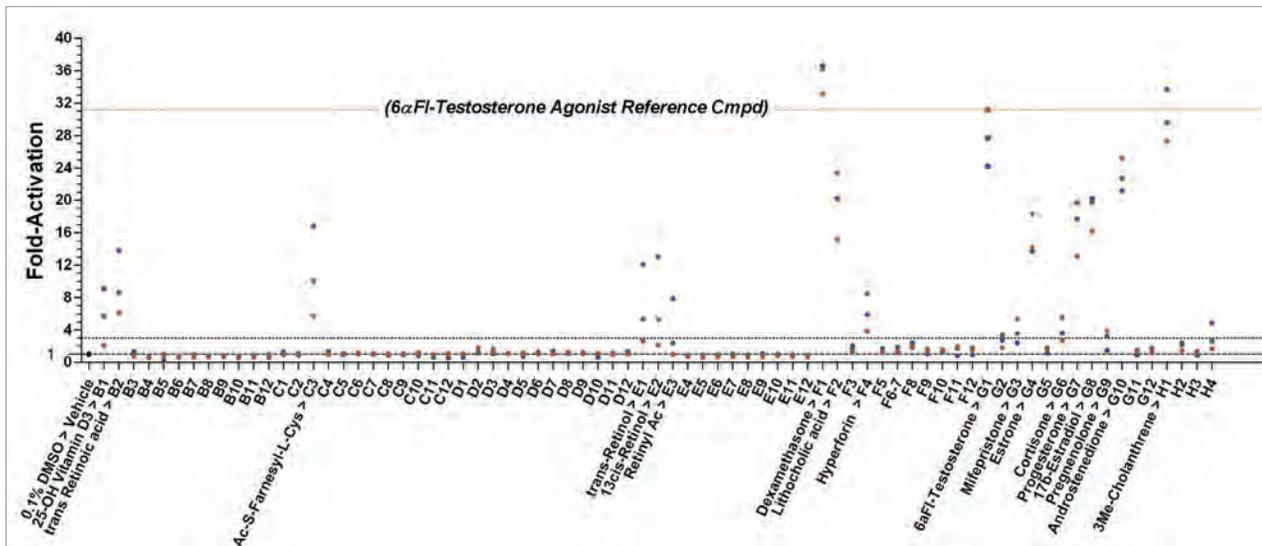


Figure 3. A. Z'-factor validation data. Z'-factor values ≥ 0.5 are indicative of an excellent assay⁵. B. EC₅₀ value of 71 pM for 6 α -FIT, and IC₅₀ value of 96 nM for BIC are equivalent to values previously generated using manual methods.

The 76 member Screen-Well Nuclear Receptor Ligand Library, containing compounds with known bioactivity to nuclear receptors, was used to validate the automated 384-well Human Androgen Receptor Assay setup for agonist and antagonist screening purposes (Figure 4).

A.



B.

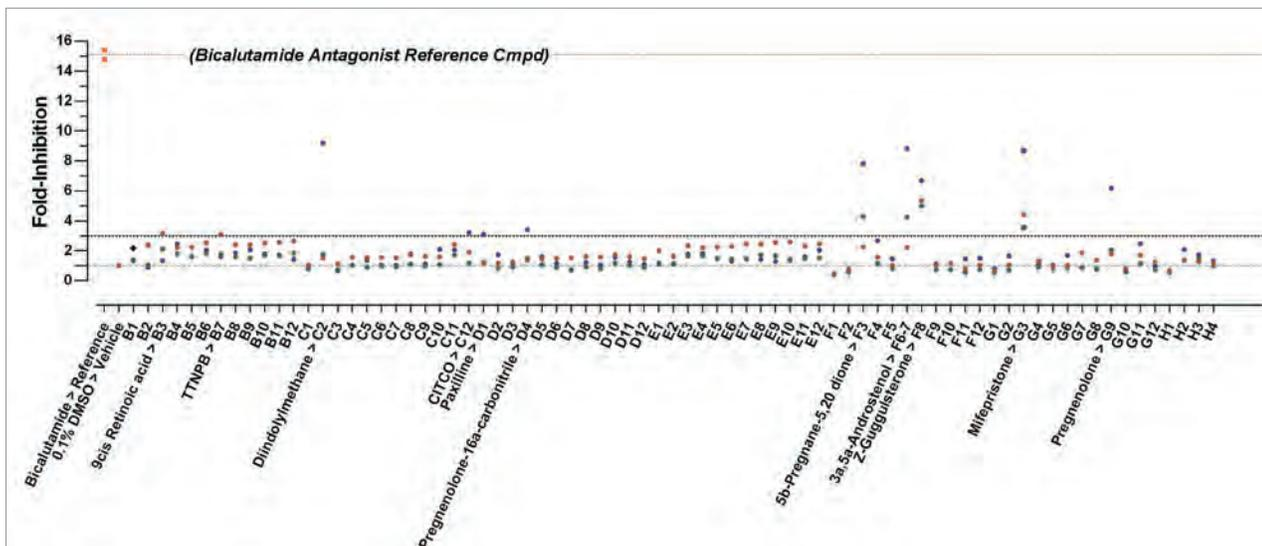


Figure 4. Human Androgen Receptor Assay Validation Screen. Test compounds were prepared at three intermediate concentrations, then screened ($n=6$) at the following concentrations: • 10 μM , • 2.0 μM , and • 0.4 μM for either (A) agonist activity or (B) antagonist activity. 0.1% DMSO was included as a vehicle control. For the antagonist assay, cells were pre-treated with 200 pM ($\sim\text{EC}_{75}$) 6α -Fluoro-Testosterone as the challenge agonist. Fold activation is defined as $[\text{RLU}_{\text{Test Cmpd}} / \text{RLU}_{\text{Vehicle Control}}]$. Fold inhibition is defined as $[\text{RLU}_{\text{Vehicle Control}} / \text{RLU}_{\text{Test Cmpd}}]$.

The luminescent signal from library compounds was compared to those generated from vehicle control wells containing 0.1% DMSO only, to generate fold activation and fold-inhibition values. Hits were defined as compounds having fold values ≥ 3 for two of the three concentrations tested. Results from known AR agonists and antagonists, such as 6α -Fluoro-Testosterone and mifepristone, demonstrate the utility of the automated 384-well Human Androgen Receptor Assay to accurately detect AR modulators.

The 506 compound Screen-Well Natural Product Library was then screened using the automated 384-well Human Androgen Receptor Assay (Figure 5).

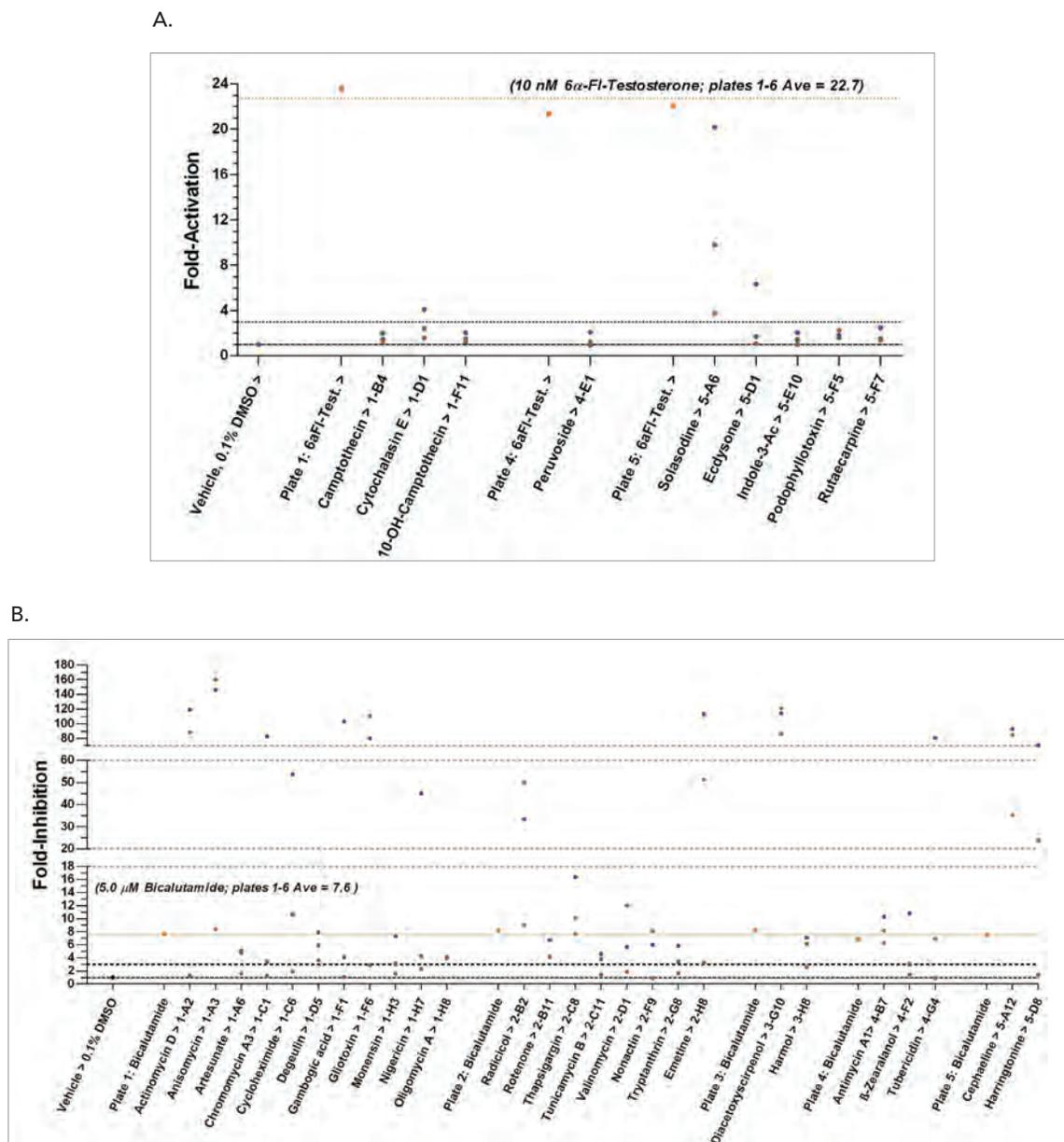


Figure 5. Summary of (A) agonist positive and (B) antagonist positive compounds from Screen-Well Natural Product Library. Test compounds were prepared at three intermediate concentrations, then screened ($n=1$) at the following concentrations: • 1:1,000, • 1:5,000, and • 1:25,000. 0.1% DMSO was included as a vehicle control. Per the antagonist assay, cells were pre-treated with 200 pM ($\sim EC_{70}$) 6α -Fluoro-Testosterone as the challenge agonist.

Following tests to remove false agonists and antagonists (data not shown), full dose-response curves were generated using the confirmed positive AR agonists and antagonists to determine EC₅₀ or IC₅₀ values (Figure 6).

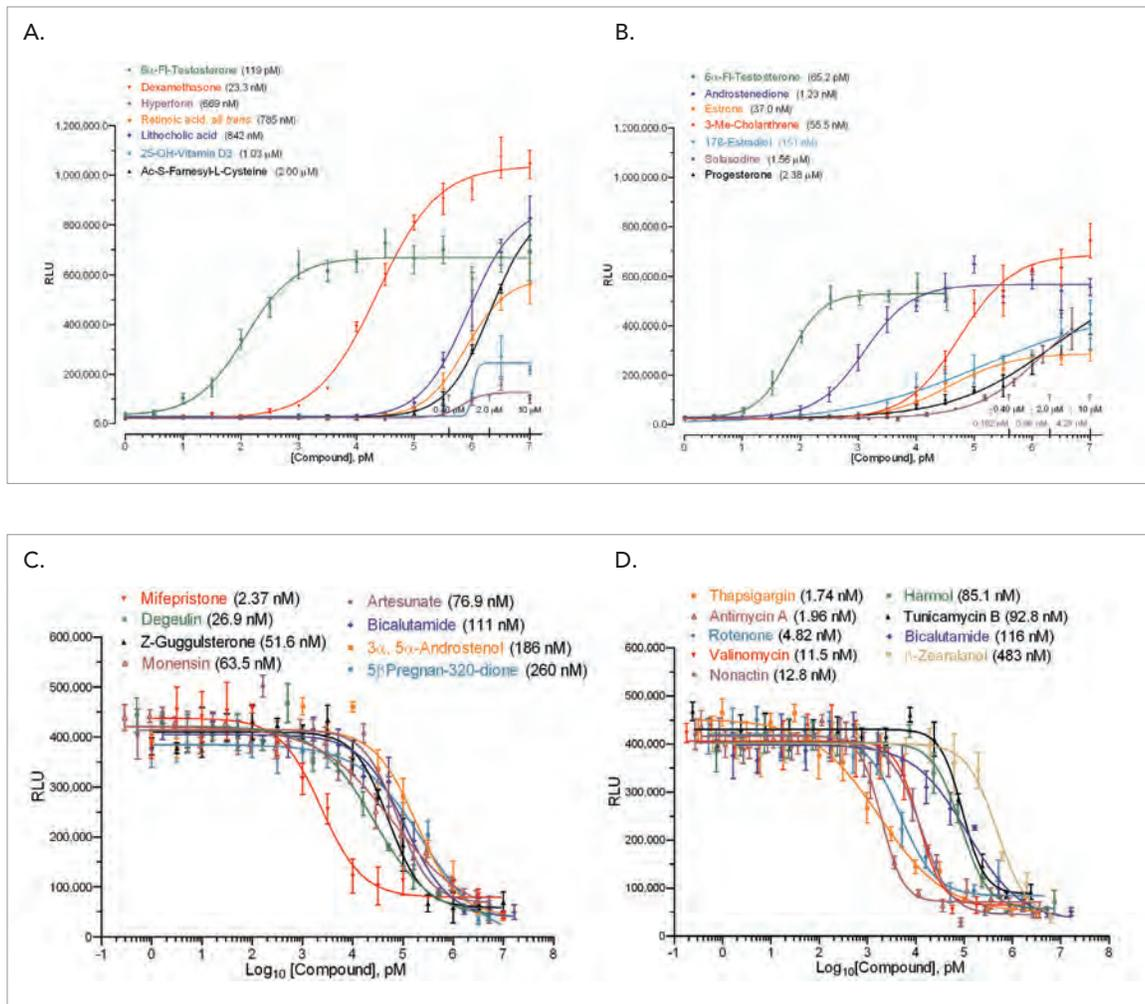


Figure 6. Positive (A, B) agonist and (C, D) antagonist compound dose response test from the Screen-Well Nuclear Receptor Ligand and Natural Product Libraries. Compound titrations were begun at 10 μ M and a 1:1000 dilution of the original stock concentration for the agonist and antagonist assays, respectively, with the exceptions of 6 α -Fluoro-Testosterone (Panel B, 33.3 nM) and Solasodine (4.8 μ M). EC₅₀ or IC₅₀ values are reported for each dose response curve.

GPCRs

G-protein-coupled receptors (GPCRs) are membrane proteins characterized by a seven-transmembrane α -helix structure. They constitute one of the largest and most diverse protein families, whose primary function is to transduce extracellular stimuli into intracellular signals⁶. GPCRs are involved in numerous physiological processes, such as behavior and mood regulation, immune response, autonomic nervous system transmission and tumor growth and metastasis. Abnormal signal transduction is related to various serious conditions, such as allergy, heart trouble, cancer, high blood pressure, and inflammation⁷. Therefore they are one of the most popular drug targets, accounting for approximately one-third of all approved drugs⁸. Continued growth in this market is expected, reaching approximately \$122 billion by 2018⁹.

GLP-1 Ligand Binding Analysis using a Cell-Based HTRF Tag-lite® Assay

The Glucagon GLP-1 Receptor Ligand Binding Assay from Cisbio Bioassays uses HEK293 cells transfected with the pSNAP-GLP-1 plasmid (PSNAPGLP1) for twenty-four hours, then subsequently labeled with the small fusion tag, SNAP-Lumi4 Tb (SSNPTBC) (Figure 7). The labeled cells are then frozen in liquid nitrogen and 10% DMSO. In the application described here, the cells were dispensed into assay plates along with sequential additions of compounds to be tested and Exendin 4-red (L0030RED), a fluorescent derivative of Exendin 4. When the fluorescent ligand was bound to the GLP-1 receptor, a time-resolved fluorescence resonance energy transfer occurred between the Lumi4 Tb donor bound to the GPCR and the red emitting labeled ligand. Competition between non-labeled compounds and the Exendin 4-red (at 4 nM) diminished energy transfer, thus leading to a decrease in HTRF signal.

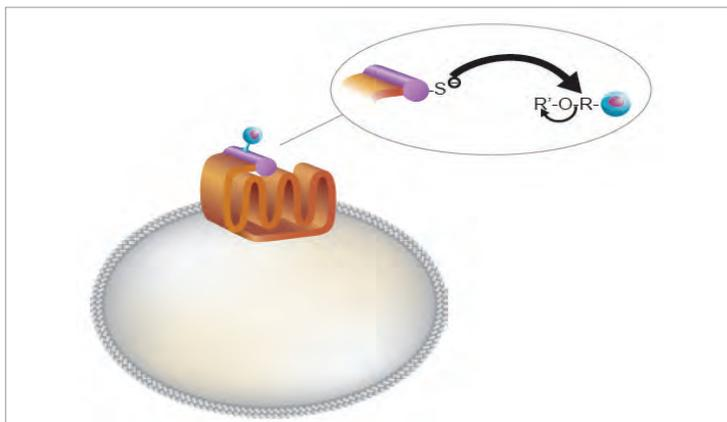


Figure 7. Tag-lite system comprised of HTRF donor fluorophore labeled GPCR. GPCR with SNAP-tag was cloned into the cell line, which can then be specifically labeled with donor fluorophore.

A total of 384 compounds from the Screen-Well Natural Product Library plates 1-4 were screened in duplicate. Compounds were pre-diluted 1:250 from the original 100% DMSO stocks and then screened at a final 1:1000 dilution. A no compound (0% inhibition) control was included as well as Exendin-4, the unlabeled GLP-1 receptor agonist, and the known antagonists GLP-1 (9-36) and Exendin-3 (9-39). When plotting the compound percent inhibition data, a distribution profile similar to that seen in larger compound library primary screens was observed (Figure 8).

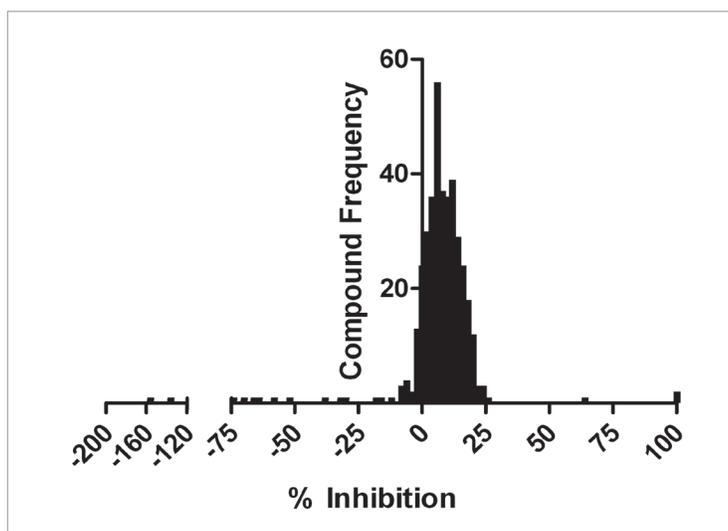


Figure 8. Natural product compound library screen percent inhibition distribution.

Most compounds exhibited little to no inhibition, while a small percentage demonstrated high positive or negative inhibition. The donor molecule fluorescent signal assessment may be used to ensure that compounds exhibited true positive or negative red-labeled agonist binding inhibition (Figure 9). This control, inherent in HTRF technologies, provided a rapid method to remove false hits. An example of a false result is Chelerythrine.

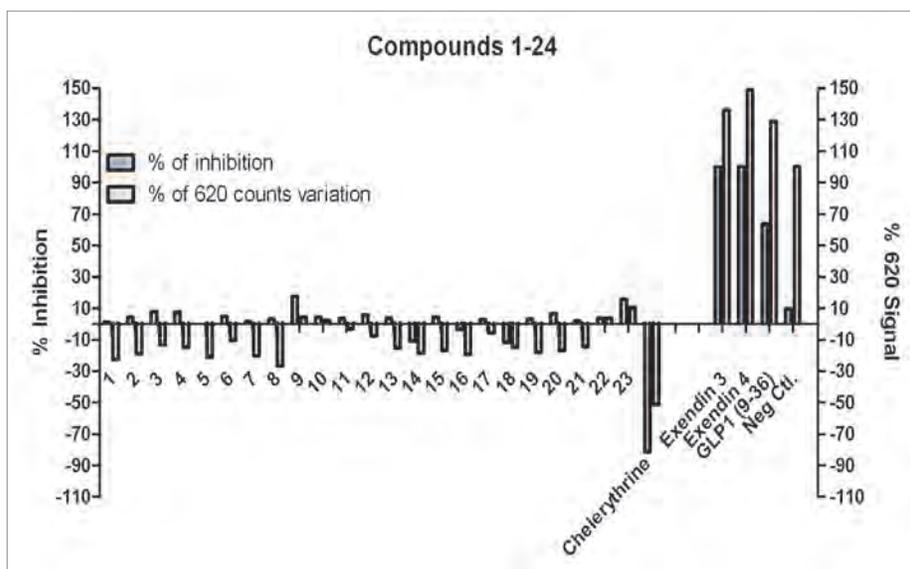


Figure 9. Representative GLP-1 receptor and assay inhibition data. Percent inhibition of red labeled Exendin-4 agonist and percent of donor molecule fluorescent values from 0% inhibition wells shown for natural product library compounds 1-24.

This compound exhibited high negative inhibition, but also decreased the donor molecule signal by approximately 50%. In contrast, Exendin-3, Exendin-4 and GLP-1 all showed high positive inhibition with no negative effect on the donor molecule's fluorescent signal.

Dose response curves were generated for those compounds exhibiting positive red agonist binding inhibition with no effect on the donor molecule fluorescent signal in the primary compound library screen (Figure 10). Eight-point 1:10 titration curves were created for each compound, starting with a 10 μ M concentration. K_i data were calculated from IC_{50} results using the Cheng-Prusoff equation¹⁰ (Table 1).

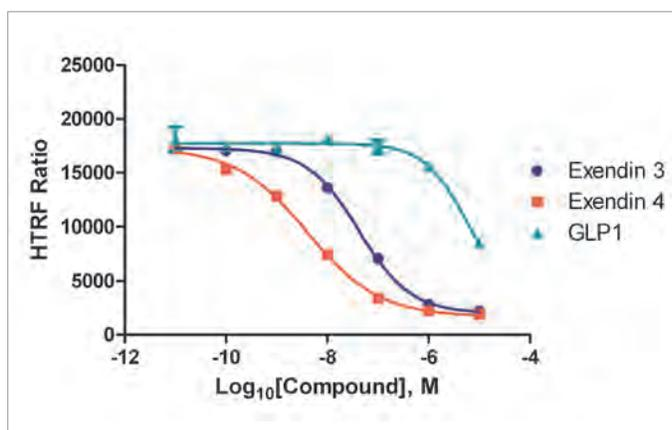


Figure 10. Dose response curves generated for positive inhibitor compounds.

GLP1 Ligand K_i Values (nM)		
Compound	Generated K_i Value	Literature K_i Value (using Radioactivity)
Exendin-3 (9-39)	11.0	7.3 ¹¹
Exendin-4	1.0	1.0 ¹¹
GLP1 (9-36)	1444	Low Binding Affinity ¹²

Table 1. K_i values for positive inhibitor compounds.

Kinases

Cellular kinases transfer phosphate groups from donors, such as ATP, to specific substrates through phosphorylation. As such, they play an important role in relaying signals from activated receptors residing at the cell membrane to the cell's interior 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, is also 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 drug development. Current global revenues for kinase inhibitors were nearly \$29.1 billion in 2011 and are projected to reach nearly \$40.3 billion in 2015¹³.

Quantification of Endogenous Cellular Kinase Activity using a Cell-based HTRF® Protein Kinase Assay

The HTRF cellular kinase assays from Cisbio Bioassays are designed to detect and study activated kinase directly in whole cells. Upon receptor activation, the kinases are activated, leading to kinase phosphorylation. After cell membrane lysis, 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 Eu³⁺-cryptate. The assay is based on a sandwich immunoassay principle. In the presence of phosphorylated kinase, and upon excitation of the Eu³⁺-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 (Figure 11).

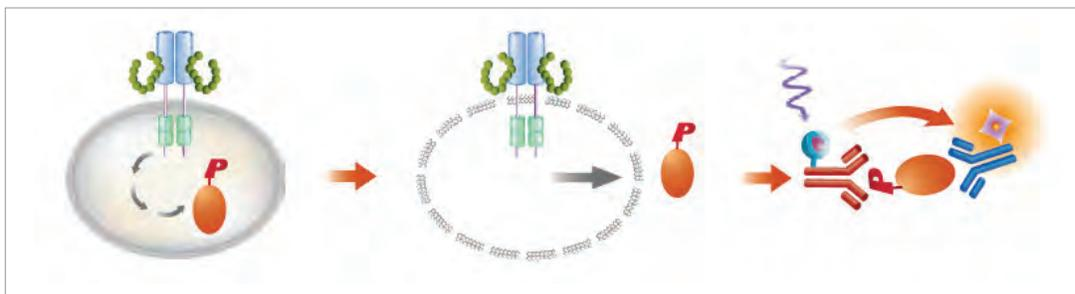


Figure 11. The HTRF Cellular Kinase assay principle. Upon receptor activation, kinases are activated, leading to kinase phosphorylation. After cell membrane lysis, phosphorylated kinase can be detected upon the addition of two monoclonal antibodies: an anti-total antibody labeled with d2 and an anti-phospho-kinase antibody labeled with Eu³⁺-cryptate.

In the work described here, the phospho-STAT3 (Tyr705) assay was used to assess potential inhibitors of the EGFR signaling pathway leading to phosphorylation of STAT3 kinase at the Tyr705 site using the human epithelial carcinoma cell line, A431. Initial testing demonstrated that the automated assay procedure was accurate and robust, as demonstrated by the EC₅₀ value for EGF stimulation of 61.5 nM (compared to 180 nM previously generated by the assay manufacturer), and a Z'-factor score of 0.79 (Figure 12).

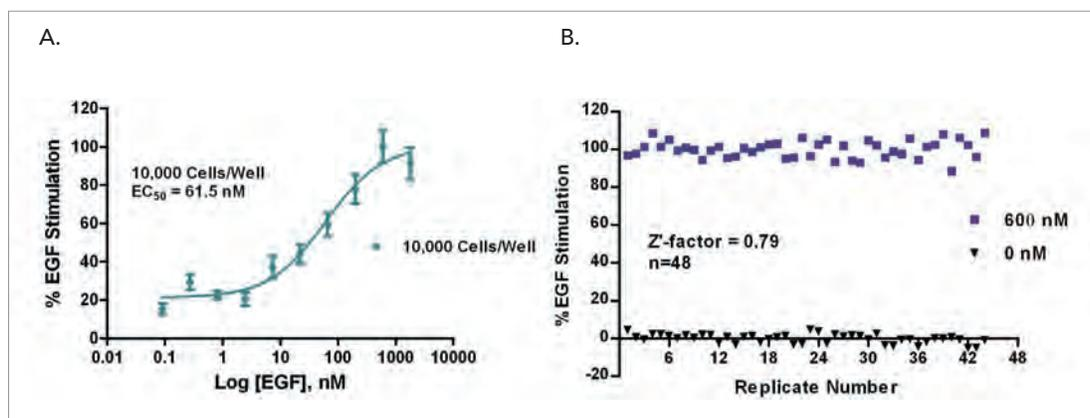


Figure 12. (A) EGF stimulation of STAT3(Tyr705) phosphorylation demonstrated as a full dose response; (B) Z'-factor score from 40 replicate measurements each at [EGF] of 600 nM and vehicle.

It is clear that many compounds demonstrate an effect on the 620 signal at higher concentrations, which may lead to false inhibition data. Further testing is necessary to determine any cytotoxic effects that potential library inhibitory compounds might have on the A431 cells, and to determine each compound's full inhibitory characteristics.

Initial dose response studies were conducted on the three control compounds included in the screen. 11-point titrations were created using a 1:2 dilution scheme starting at a 1X concentration of 100 μM (Figure 15). IC_{50} values (Table 2) were calculated from each inhibition curve.

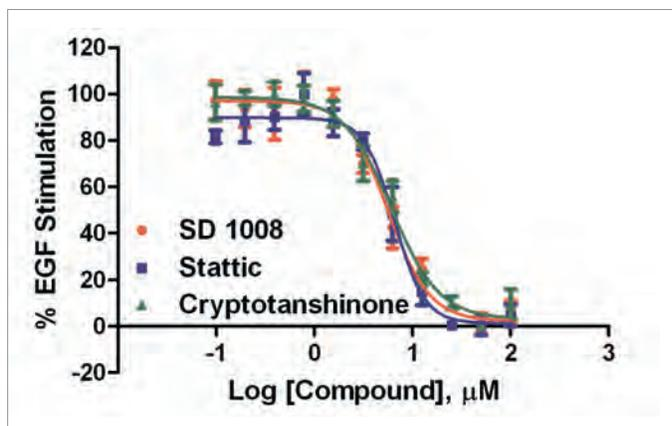


Figure 15. Inhibition of EGF-induced STAT3(Tyr705) phosphorylation using the small molecule inhibitors, SD 1008, stattic and cryptotanshinone.

STAT3 Inhibitor IC_{50} Values (μM)		
Compound	Generated IC_{50} Value	Literature IC_{50} Value
SD 1008	5.6	Inhibition seen at 10 μM in as little as 30 minutes ¹⁴
Stattic	6.6	5.1 ¹⁵
Cryptotanshinone	6.1	4.6 ¹⁶

Table 2. IC_{50} values for known inhibitors of STAT3 activation.

Epigenetics

The DNA molecules in each cell are wrapped around histone proteins, and collectively referred to as chromatin. This serves to compact the DNA strand, and also controls gene expression. Changes to the chromatin structure are controlled by the addition or removal of epigenetic factors, such as methyl or acetyl groups, to histone tails. A relaxed chromatin structure allows for gene transcription, whereas the opposite silences that part of the genetic code. This is a normal process during cell differentiation. However, the loss of proper epigenetic control can lead to aberrant gene expression and the development of multiple diseases, including diabetes, cancer, or neurological disorders. This has led to much new research, and has made epigenetics one of the fastest growing drug discovery markets, with projections that the worldwide market will reach \$18.2 billion by 2015¹⁷.

Assessment of HDAC1 Inhibition using a Bioluminescent Histone Deacetylase I/II Assay

The HDAC-Glo™ I/II Assay from Promega Corporation (Madison, WI) is a single-reagent-addition, homogeneous, luminescent assay that measures the relative activity of histone deacetylase (HDAC) class I and II enzymes from cells, extracts or purified enzyme sources. The assay is broadly useful for class I and II enzymes, and uses an acetylated, live-cell-permeant, luminogenic peptide substrate that can be deacetylated by HDAC activities (Figure 16). Deacetylation of the peptide substrate is measured using a coupled enzymatic system in which a protease in the developer reagent cleaves the peptide from aminoluciferin, which is quantified in a reaction using Ultra-Glo™ Recombinant Luciferase. The HDAC-mediated luminescent signal is proportional to deacetylase activity.

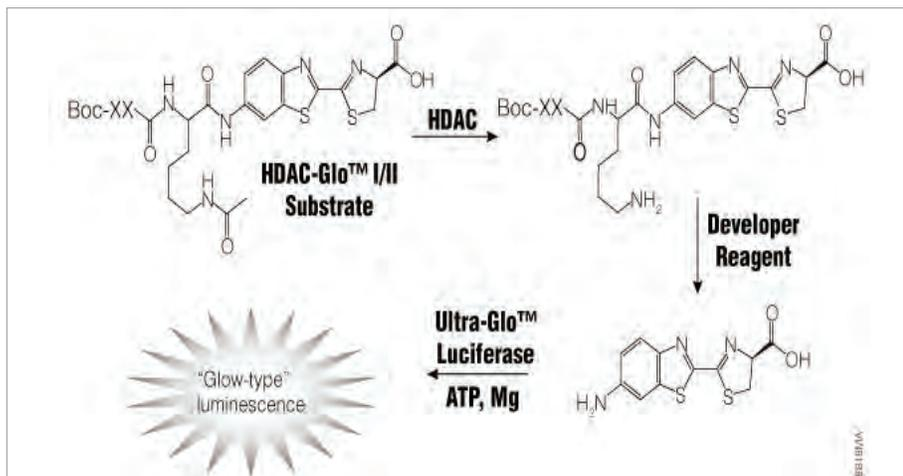


Figure 16. HDAC-Glo I/II Assay Principle. All three enzymatic modifications of the substrate occur in coupled, nearly simultaneous reactions upon addition of a single reagent.

Here we used the HDAC-Glo I/II Assay to screen for HDAC1 inhibitors in a low volume 384-well format. Initial experiments, using a pre-determined enzyme concentration of 10 ng/mL (data not shown), validated that the automated assay procedure delivers accurate and robust results (Figure 17). The IC_{50} value for Trichostatin A of 1.04 nM agrees with previously published data¹⁸, and the Z' -factor score of 0.87 confirms its suitability for screening.

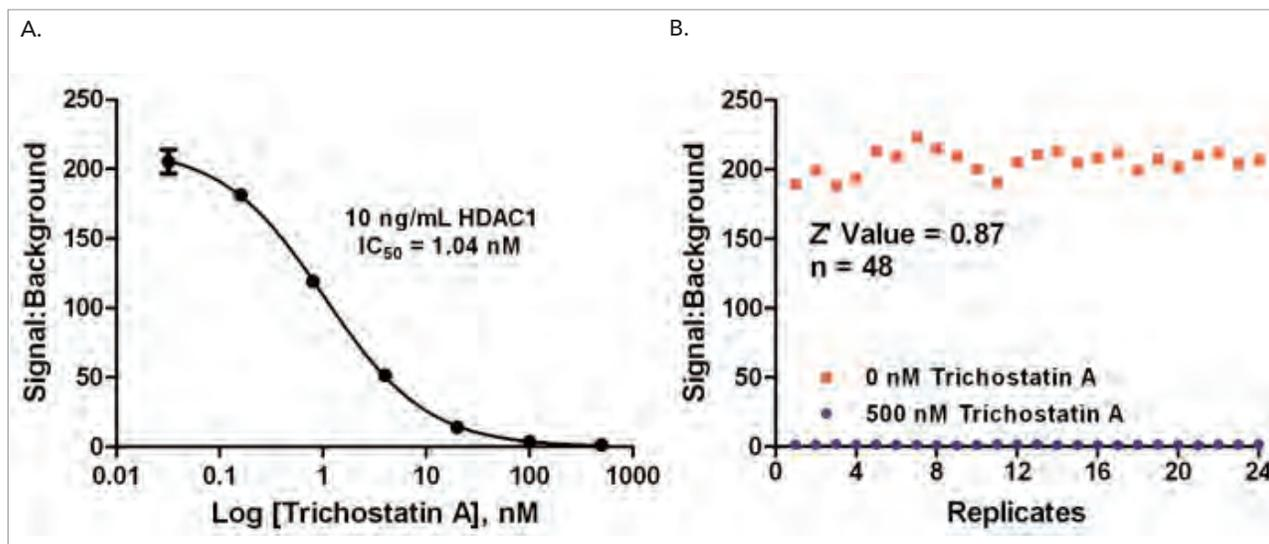


Figure 17. (A) Dose response curve showing inhibition of HDAC1 enzyme activity by Trichostatin A; (B) Z' -factor score from 24 replicate measurements each at [Trichostatin A] of 500 nM and vehicle.

A compound screen was then performed using the Screen-Well Epigenetics Library. Five additional compounds were also included which are shown to inhibit the activity of various epigenetic proteins. Compounds were diluted from the 10 mM, 100% DMSO stock to three intermediate concentrations and then screened at final concentrations of 20, 2, and 0.2 μ M. Percent inhibition (Figure 18) was calculated by comparing the luminescent values from each compound concentration tested to a no compound, uninhibited enzyme reaction control.

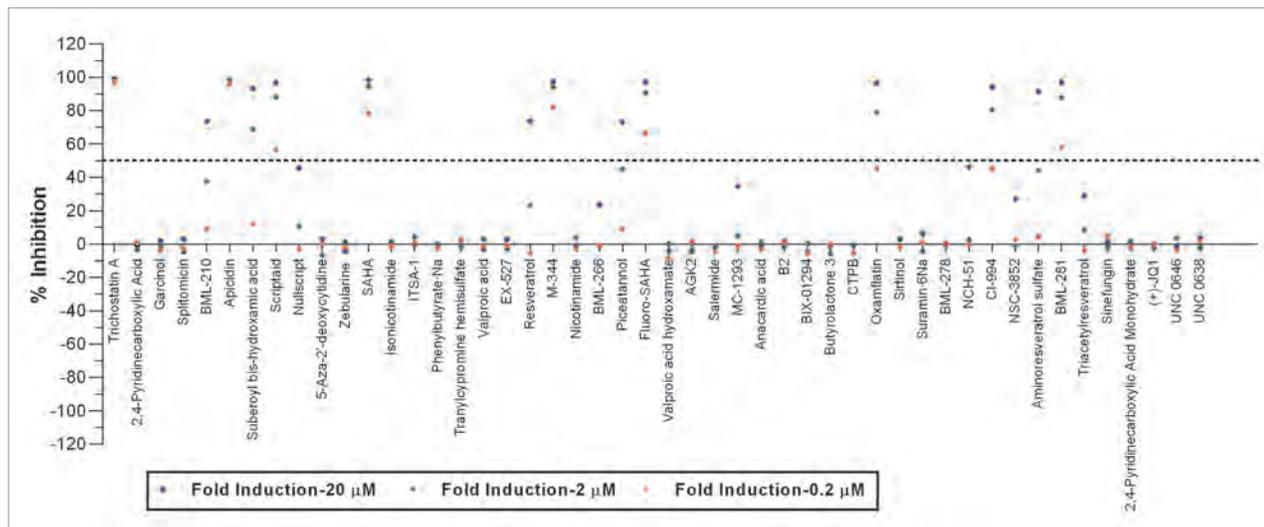


Figure 18. % Inhibition results from HDAC1 assay screen using Screen-Well Epigenetics Library.

Compounds exhibiting greater than 50% inhibition of enzyme activity at any of the three concentrations tested, including known HDAC inhibitors, and the SIRT activators Resveratrol, Piceatannol, and Aminoeresveratrol Sulfate, were carried forward for dose response testing. Two different tests were completed. 12-point titrations were created for each test using a 1:4 dilution scheme starting at a 1X concentration of 100 μ M. The first determined the specific inhibitory characteristics of each lead compound (Figure 19A). IC_{50} values were calculated from each inhibition curve. The second test determined the inhibitory potential of each compound on the components used in the detection step (Figure 19B). A nonacetylated HDAC-Glo I/II control substrate was added to the detection reagent. The control substrate does not require deacetylation to produce luminescence, and therefore is not affected by HDAC inhibitors.

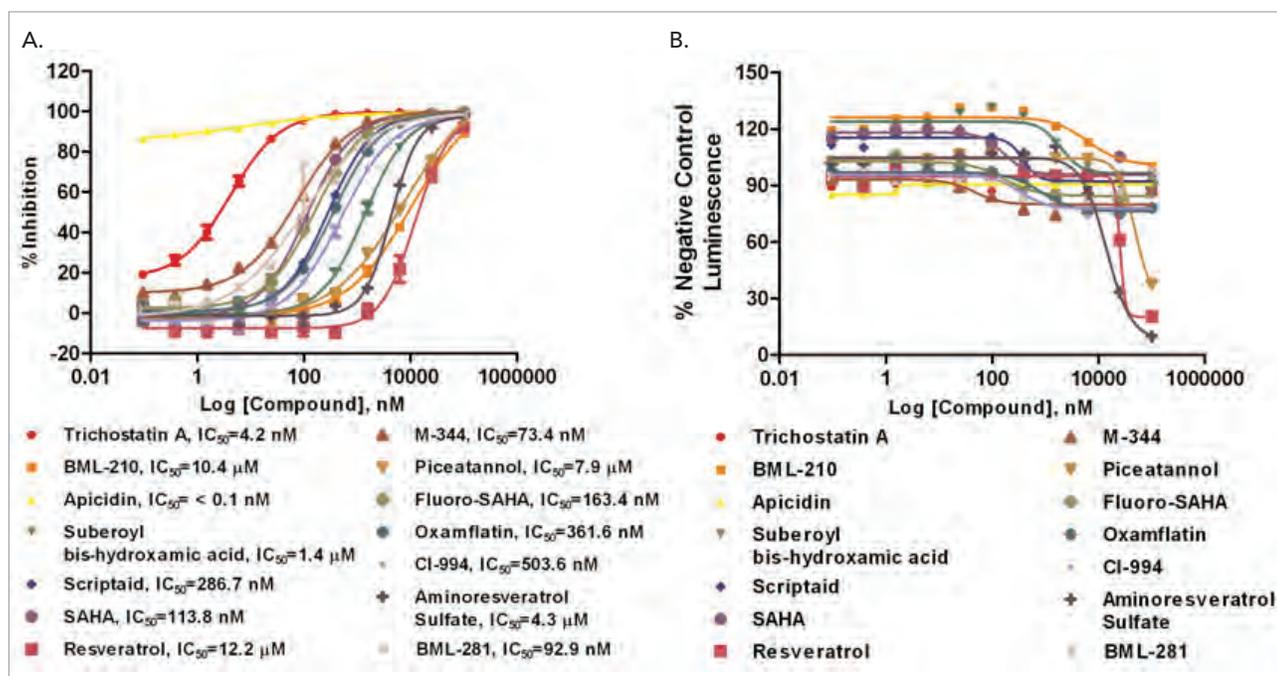


Figure 19. Dose response curves generated for "hit" compounds using (A) standard detection reagent components, or (B) detection reagent plus control substrate.

The dose response tests confirmed that each known HDAC inhibitor demonstrated true HDAC1 activity inhibition, as no significant negative change in luminescent signal was seen when each compound was tested with the control substrate. This was in contrast to the three SIRT activator compounds, which showed evidence of a negative effect on the detection reagent components at high compound concentration. Therefore these compounds should not be included with the others as actual inhibitors of HDAC1 activity.

Conclusions

While the high throughput screening market continues to feel great pressure, leading to a number of changes in method and ways of thinking, the tools available for use in HTS also continue to adapt and evolve. New assay chemistries, both biochemical and cell-based, give more precise answers using increasingly *in-vivo*-like models. BioTek's liquid handling and detection instrumentation also meets the new demands of the industry. The Precision™, EL406™ and MultiFlo™ can be used to automate numerous assay procedures, including intricate tasks such as medium exchanges and buffer washes for cell-based assays. The appropriate Synergy™ Multi-Mode Microplate Reader can then be incorporated to detect the emission signal from the specific assay. Together, the combination of assay chemistry and instrumentation create an ideal solution to meet the needs of today's high throughput screening laboratory.

References

1. PRWeb. http://www.prweb.com/releases/high_throughput_screening/drug_discovery/prweb8512562.html (accessed Dec 21, 2012). Global Industry Analysts, Inc., press release.
2. Xiaohong, X.; Lei, Y.; Luo, J.; Wang, J.; Zhang, S.; Yang, X.; Sun, M.; Nuwaysir, E.; Fan, G.; Zhao, J.; Lei, L.; Zhong, Z. Prevention of β -amyloid induced toxicity in human iPS cell-derived neurons by inhibition of Cyclin-dependent kinases and associated cell cycle events. *Stem Cell Res* [In Press, Accepted Manuscript]. <http://dx.doi.org/10.1016/j.scr.2012.11.005>. Published Online: Dec 7, 2012. (accessed Dec 21, 2012).
3. Robinson-Rechavi, M.; Garcia, H.; Laudet, V. The Nuclear Receptor Family. *J Cell Sci.* **2003**, *116*(4), 585-586.
4. Via, M. *Nuclear Receptors: The Pipeline Outlook*; Market report, Insight Pharma Reports: Needham, MA, 2010.
5. Zhang, J.; Chung, T.; Oldenburg, K. A Simple Statistical Parameter for Use in Evaluation and Validation of High Throughput Screening Assays. *J Biomol Screen.* **1999**, *4*(2), 67-73.
6. Kroeze, W.; Sheffler, D.; Roth, B. G-protein-coupled receptors at a glance. *J Cell Sci.* **2003**, *116*(24), 4867-4869.
7. Suwa, M.; Sugihara, M.; Ono, Y. Functional and Structural Overview of G-Protein-Coupled Receptors Comprehensively Obtained from Genome Sequences. *Pharmaceuticals.* **2011**, *4*(4), 652-664.
8. Rubenstein, K. *GPCRs: Dawn of a New Era?*; Market report, Insight Pharma Reports: Needham, MA, 2008.
9. *G-Protein Coupled Receptors (GPCRs) Market - Global Industry Size, Market Share, Trends, Analysis and Forecast, 2012 - 2018*; Market report, Transparency Market Research: Albany, NY, 2012.
10. Cheng, Y.; Prusoff, W. Relationship between the inhibition constant (K_i) and the concentration of inhibitor which causes 50 percent inhibition (I_{50}) of an enzymatic reaction. *Biochem Pharmacol.* **1973**, *22*(23), 3099-3108.
11. Jorgensen, R.; Martini, L.; Schwartz, T.; Elling, C. Characterization of Glucagon-Like Peptide-1 Receptor β -Arrestin 2 Interaction: A High-Affinity Receptor Phenotype. *Mol Endocrinol.* **2005**, *19*(3), 812-823.
12. Knudsen, L.; Pridal, L. Glucagon-like-peptide-1-(9-36) amide is a major metabolite of glucagon-like-peptide-1-(7-36) amide after *in vivo* administration to dogs, and it acts as an antagonist on the pancreatic receptor. *Eur J Pharmacol.* **1996**, *318*(2-3), 429-435.

13. PR Newswire. <http://www.prnewswire.com/news-releases/kinase-inhibitors-market-to-reach-116-billion-in-us-403-billion-globally-by-2016-161937715.html> (accessed Dec 19, 2012). Global Information Inc., press release.
14. Duan, Z.; Bradner, J.; Greenberg, E.; Mazitschek, R.; Foster, R.; Mahoney, J.; Seiden, M. 8-Benzyl-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.
15. Schust, J.; Sperl, B.; Hollis, A.; Mayer, T.; Berg, T. Stattic: A Small-Molecule Inhibitor of STAT3 Activation and Dimerization. *Chem Biol.* **2006**, *13*(11), 1235-42.
16. Shin, D.; Kim, H.; Shin, K.; Yoon, Y.; Kim, S.; Han, D.; Kwon, B. Cryptotanshinone Inhibits Constitutive Signal Transducer and Activator of Transcription 3 Function through Blocking the Dimerization in DU145 Prostate Cancer Cells. *Cancer Res.* **2009**, *69*(1), 193-202.
17. PR Newswire. http://www.prweb.com/releases/epigenetics/drug_discovery/prweb3553614.htm (accessed Dec 21, 2012). Global Industry Analysts Inc., press release.
18. Viquushin, D.; Ali, S.; Pace, P.; Mirsaide, N.; Ito, K.; Adcock, I.; Coombes, R. Trichostatin A is a histone deacetylase inhibitor with potent antitumor activity against breast cancer in vivo. *Clin Cancer Res.* **2001**, *7*(4), 971-976.