Development of Simple, Robust Bioassays for Biosimilar & Biobetter Development

Abhishek Saharia, Ph. D., Mong Saetern, Hanako Daino-Laizure, Sangeetha Gunthuri, Rajini Bompeli and Jane Lamerdin, Ph.D. DiscoveRx Corporation, Fremont, CA 94538-3142

Abstract One of the major bottlenecks in the development of biosimilars is the need for good bioassays to create potency, stability and neutralizing antibody (NAb) assays. Ideal bioassays need to reflect the clinical mechanism of action (MOA) of the biopha maceutical drug and should be simple, precise, reproducible and robust. To further enable NAb assays, there is an added need for serum/matrix tolerance and sensitivity. Here, we discuss the development and application of diverse PathHunter® cell-based assays that cover distinct cellular mechanisms either downstream of or at the level of receptor engagement by large molecules. Importantly many of these assays rely on expression of native receptors thereby reflecting the clinical MOA of the biopharmaceutical drug. These assays are highly specific, quantitative, scalable, robust and utilize a homogenous mix-and-read protocol, which facilitates rapid and reproduc ible detection of drug potency. The technology is also amenable to accurate and sensitive detection of neutralizing antibodies. even in high concentrations of human serum through a simple chemiluminescent output. The assays are developed in a convenient ready-to-use format that minimizes assay variability often occurring due to cell culture. The cell preparation, bioassay protocol and reagents have been optimized to provide superior bioassay performance with high reproducibility (<7% RSD). PathHunter Enzyme Fragment Complementation Active Enzyn Inactive Fragr Light coveRx's proprietary PathHunter Enzyme Fragment Complementation (EFC) technology consists of the β-galactosidas ponents, the enzyme donor peptide (ED) and enzyme acceptor (EA). When brought together in close proximity, ED cos a active enzyme catalyzes the substrate generating chemiluminescent light, providing a highly amplified ail and thus an assay of high sensitivity. e (β-gal) enzyme, split into two inactive mplements with EA forming active β-gal componer The active Assay Principle For Anti-TNFa Bioassay NF-κB Signaling - IκB Degradation TNF TNFF 22 Assay Principle. TNF α binds to and stimulates the TNF receptor, triggering a downstream signaling cascade that activates the NF κ B pathway. This leads to the degradation of the kB protein. Here, the PL-tagged kB protein is expressed in THP-1 cells. Stimulation of cells with TNF α results in dose-dependent degradation of the key proteins on the receptor can be detected via kB degradation in a very sensitive manner by the addition of the text of the substrate, resulting in chemiluminescent light that is proportional to the amount of PL-tagged kB present in the cells. Therefore, a higher activation of TNF will result in lower assay signal. Robust Bioassay Response with Anti-TNF drugs - Infliximab and Adalimumab TNFa Response in THP1 Cells Remicade & Humira Assay 2500000 150000 TNFα [M] Romicado Remicad Humira h 200000 10000

H JLC RLU 1000000 50000 50000 0+ 10-1 15 10-14 10-13 10-12 10-11 10-10 10-9 10-8 10-10-14 10-13 10-12 10-11 10-10 10-9 10 TNFα [M] Antibody [g/mL] +100pM TNFo HillSlope -0.7918 TNFα [M] Remicade [g/mL] Humira [g/mL] 6.001e-012 -1.099 1.085e-012 2.024 8.470e-010 EC₅₀ S/B 2.025 9.257e-010 9.7 S/B 5.6 3.2 4.8

Continuous culture cells for the Anti-TNF assay. Cells were cultured and plated overnight before the assay was conducted. These assays demonstrate a high reproducibility (<15% intra-plate CV) in response to TNF α (left), and when treated with the innovator molecules Infliximab (Remicade[®]) and Adalimumab (Humira[®]) (right).







Using EFC, assays have been created to measure neceptor heterodimerization, homodimerization or co-receptor meruitment. The two target neceptors are tagged with ProLink™ (PK) or Enzyme Acceptor (EA). Upon ligand-induced activation, the receptors naturally dimerize forcing the two β-gal components to complement and create an active enzyme. Active β-gal generates a chemiluminescent signal in the presence of substrate, signaling the formation of receptor dimers.

Bioassay for Avastin and Other Anti-VEGF-A Drugs



Avastin Assay Provides Highly Reproducible Data. VEGF-A is known to cause homodimerization of VEGFR2 (KDR), as the first step in the activation cascade of these receptors. Anti-VEGF-A antibodies such as Avastin prevent this dimer formation, leading to inhibition of VEGF-A dependent signaling. Here we have tested the VEG-FR2 homodimer assay with VEGF-A and Avastin, demonstrating a robust response and a high level of reproducibility with multiple runs. This assay provides a much simpler method for testing Avastin potency and NAb assay development than the available HUVEC cell proliferation assay.

Bioassay for Anti-IL-17A and Anti-IL-17F Antibodies



Highly Reproducible IL17RA/IL17RC Assay to Develop Anti-IL-17 Drugs. IL-17A or IL-17F will cause heterodimerization of IL-17RA and IL-17RC receptors, as the first step in their activation cascade. Anti-IL-17 antibodies will prevent the formation of this heterodimer, leading to inhibition of IL-17 receptor signaling. Here, we have developed and tested the IL-17A and IL-17F in this assay, demonstrating a robust response and a highly reproducible response. This assay provides a much simpler method for testing IL-17 potency and NAb assay development than the available assays.

imple, Robust Assay for Insulin Biosimilar Development





Vial#3

4110

42297

-7.090

1.008

10.0

Bioassay development for Insulin Biosimilars and Biobetters – (Left) Insulin binds and activates the Insulin receptor, leading to conformation change and trans-phosphorylation. This results in the recruitment of an SH2-domain protein to the receptor, as part of the natural signaling cascade. By expressing a tagged Insulin receptor and an EA-tagged SH2-domain protein, we can monitor target-proximal receptor activation with a simple chemiluminescent read (Right) Two different sample of Insulin provide an appropriate response in the Insulin bioassay. sing a PK-





Frozen Ready-to-Assay Cells Provide Highly Reproducible Data. Anti-TNFα bioassay from frozen ready-to-use cells, shows excellent sensitivity and reproducibility in response to TNFα (left), and when treated with innovator molecules Remicade and Humira (right). As noted above, frozen cells were plated directly onto the assay plate post thaw, incubated overnight to allow them to attach and assayed the next day with TNFα and the inhibitor molecules.

Robust Bioassay For Anti-RANKL Antibody - Denosumab



	RANKL	Denosumal
HillSlope	-2.628	3.565
EC 50	3.329e-010	2.816e-007
S/B	2.5	3.3

Assay Principle & Robust Response to Denosumab (Anti-RANKL) – RANKL activates TNFR superfamily member, RANK, which leads to the activation on the NFxB response. By expressing PL-tagged IkB in U2OS cells that express RANK on their surface, we can monitor receptor activation, in a similar manner to TNFR activa-tion. 24 hours after plating the engineered cells, Denosumab was mixed with 10MR RANK ligand (RANKL) and added to the cells for 1hr. Following the addition of PathHunter detection reagent we observe a robust response to both agonist and inhibitor drug.

PathHunter bioassay kits use a simple homogenous protocol with rapid results. (A) Ready-to-assay cells from DiscoveRx are plated on a 96-well plate or 384-well plate and incubated for about 24 hours at 37°C. The agonist/test molecule is added to the plate and incubated for 1-3 hours. The detection reagents are added sequentially in two additions steps and the chemiluminescent signal can be detected on any plate-reading luminometer. (B) Cells manufactured in bioassay kits are meant for single use in ready-to-assay vials. The fozen cells are plated directly onto plates to run the assay and this format has several advantages as outlined above.

Summary & Conclusions

Potency and NAb assays available for > 30 biosimilar targets.

• To minimize assay variability, the bioassays are prepared in ready-to-assay kits with all required assay components, including frozen, single use cell vials

- Cells in ready-to-assay format minimize assay variability stemming from the use of continuous culture cells

- Each kit contains enough material for 1000 datapoints (10 x 96-well plates)

- Customized lots of up to 1000 vials are prepared for each project



tel | 1.866.448.4864 tel | +44.121.260.6142 e | info@discoverx.com

© 2014 DiscoveRx Corporation. All Rights Reserved. 20239 090314