Comprehensive Cell-based Assays for Characterization & Development of Agonistic Antibodies as Therapeutics for Cancer & Autoimmune Diseases

Venkatesh Chari, Luhan Yang, Jane E. Lamerdin, Jennifer Lin-Jones, & Gaurav Agrawal Eurofins DiscoverX[®] | Fremont, CA USA

Abstract

While checkpoint blockade antibodies are established cancer therapeutics, they have emerged as promising therapeutics for suppressing inflammation in autoimmune diseases. Their ability to regulate immune function targeting both co-stimulatory & co-inhibitory receptors renders them as emerging therapeutics in autoimmune diseases & for restoring tolerance in graft-versus-host disease (GVHD) complications. Existing assays frequently fail to accurately assess the agonistic activity of antibodies as they cannot replicate the necessary physiologically relevant conditions.

We present case studies demonstrating the pivotal role of cell-based assay approaches developed by Eurofins DiscoverX in driving the accurate characterization & clinical development of agonistic antibody therapeutics targeting checkpoint receptors. Additionally, we demonstrate the indispensable role of Fcy Receptor (FcyR) mediated antibody clustering as part of the assay design in capturing the agonist potential of antibodies.

Active Areas of Clinical Development for Agonistic Antibodies



Functional Cell-based Assays for Checkpoint Receptor Targets Based on EFC Technology



Figure 1. Eurofins DiscoverX's Enzyme Fragment Complementation (EFC) Assay Design for the CD40 Checkpoint Receptor. The EFC technology is based on a split β -galactosidase (β -gal) enzyme system. Individually, the enzyme donor (ED) peptide & enzyme acceptor (EA) fragment are inactive. When brought in proximity, the ED complements with EA forming an active β -gal enzyme that hydrolyzes substrate resulting in a chemiluminescent signal. The PathHunter NIK Signaling Cell Line was engineered to express ED-tagged NIK protein. Addition of CD40L (ligand) leads to activation of the NF-KB pathway resulting in stabilization & accumulation of the NIK protein with increasing concentrations of CD40L. Addition of substrate, EA, & buffer, lyses the cells & forces complementation of the ED & EA enzyme fragments. This results in the formation of a functional enzyme that hydrolyzes substrate to generate a chemiluminescent signal.

FcyR Mediated Clustering Enhances Agonistic Effect of Antibodies Targeting Co-stimulatory Checkpoint Receptors Assay Design with FcyR Clustering Cells Significantly Augments the Agonistic Effect of Anti Co-stimulatory Checkpoint Antibodies FcyR Cells Antigen Agonistic Antibody Candidates Receptors Add Detection Incubate Reagent Read Plate Plate Target Cells ++++ FcyR Cells Target Cells

Figure 2. Modified Assay Design Including FcyR Clustering Cells. Using FcyR clustering cells (FcyR1a, 2a, 2b, or 3c) in the assay design aids in clustering the antibodies & presenting to the target cells resulting in enhanced signaling compared to standard assay design without FcyR mediated clustering.



Figure 3. Augmented Agonism of Target Receptors. A. & B. Enhanced agonism observed with mouse FcyR2b (mFcyR2b) clustering cells in the PathHunter CD137 & OX40 Signaling Assays, compared to the anti-CD137 or anti-OX40 alone, respectively. C. & D. Agonistic activities of anti-CD40 & anti-TWEAKR antibodies were assessed in the PathHunter NIK Signaling Assay, where the target receptors, CD40 & TWEAKR, are endogenously expressed. Adding mFcyR2b clustering cells to the assay design significantly enhances the efficacy & potency of these antibodies compared to the antibody alone.



Introducing FcyR Clustering in the Assay Design Helps Identify a Strong Agonistic Effect in Antibodies Previously Established as Antagonists



Figure 4. Checkpoint Receptors Signaling using Modified Assay Design for Capturing Enhanced Agonistic Activity of Antibodies. A. The PathHunter PD-1 Signaling Assay measures SHP1 recruitment to PD-1 ITIM motifs. When PD-L1 Ligand Cells are co-cultured with PD-1 Signaling Cells, PD-L1 activates the PD-1 receptor & SHP1 or SHP2. SH2-EA enzyme fragment fusion protein is recruited to PD-1 tagged with the complementary ED tag. Anti-PD-1 antibody disrupts the interaction between PD-1 & PD-L1, inhibiting PD-1 signaling, resulting in a loss of chemiluminescent signal. B. Results from the PathHunter PD-1 (SHP1) Signaling Assay. Anti-PD-1 antibody (NAT105) was used to block PD-1 activation mediated by PathHunter PD-L1 Ligand Cell Line in co-culture. C. Assay design representing PD-1 Signaling driven by SHP1 or SHP2 recruitment in agonist mode when including FcyR clustering cells, in the absence of PD-L1 cells, results in enhanced signaling. D. Addition of FcyR2a cells in the assay design significantly enhances the agonistic activity of Pembrolizumab (anti-PD-1 antibody) thereby resulting in suppression of T-cell activation. E. & F. Results from the PathHunter BTLA & CTLA4 Signaling Assays, which measure SHP recruitment as an activation mechanism. mFcyR2b cells enhance the agonistic effect of anti-BTLA & anti-CTLA4 antibodies as observed by the increased efficacy & potency in the presence of clustering cells compared to the antibody alone.



Primary & Immortalized Immune Cells with Endogenous Fcy Receptor Expression also Augments Agonistic Activity of Pembrolizumab

HB14, QA18A47) in the absence of FcyR clustering cells. B. Enhanced agonism for all three clones of anti-CD40 antibody was observed with clustering mediated by PathHunter mFcyR2b cells, exhibiting a rank order potency with HB14 being most potent clone & the 5C3 clone being the least potent clone.

Summary

- FcyR mediated clustering significantly augments agonistic activity of anti-checkpoint antibodies offering a new avenue to design therapeutics against auto-immune diseases
- Recording direct activation of checkpoint receptors through SHP recruitment helps identify a previously unregistered agonistic activity of co-inhibitory checkpoint antibodies
- Availability of a comprehensive portfolio of cell-based assays for co-inhibitory & co-stimulatory checkpoint receptors along with all major FcyR clustering cell lines enables screening & characterization towards clinical development of agonistic antibodies