Measuring Receptor Dimerization To Create Functional Cell-Based Assays for ~85% of the Interleukin Receptor Family

Scott Gridley, Ph. D.*, Abhishek Saharia, Ph. D., Hyna Dotimas, Sangeetha Gunthuri, Hanako Daino-Laizure Ph. D., Albert Doan, Phil Achacoso, and Jane Lamerdin Ph. D.

DiscoveRx Corporation, Fremont, CA 94538-3142

Abstract

Ligand-induced receptor dimerization is an early functional step in receptor activation, representing the most proximal, functional readout for receptor activation. It is well understood that the family of Interleukin receptors will dimerize with the other members of its family leading to a complicated signaling cascade that is critically involved in a variety of auto-immune, inflammatory and oncogenic diseases. Surprisingly, existing cellular assays have been unable to faithfully monitor these interactions in a proximal manner to the receptor. Here we present a novel application of the Enzyme Fragment Complementation system to monitor receptor-receptor interactions at the surface of intact live cells, applicable to diverse receptor types such as Interleukin receptors, BMP receptors, receptor tyrosine kinases and cytokine receptors, with a specific focus on the interleukin family of receptors. The high specificity, simplicity of the assay protocol, large signal to noise ratio, serum tolerance and reproducibility of these assays has enabled their use in cell-based screening, functional characterization, QC lot release assays and neutralizing antibody studies. Examples discussed here include assays for the IL-1, IL-2, IL-6, IL-10 receptor families, with assays developed or in development for ~85% of the Interleukins and their receptors.



Structura une style in interesting in interesting is because the starting of the starting of the starting starting is a starting starting in the starting starting is a starting starting in the starting starting is a starting starting in the starting starting is a starting starting is a starting starting is a starting starting starting is a starting starting starting is a starting starting is a starting starting starting starting starting is a starting sta



PathHunter dimerization assay technology has been utilized to construct assays to interrogate an early event in the interleukin pathway signaling cascade. In brief, one receptor in the dimer pair is tagged with the PK fragment of the β -pal reporter, while the second subunit is fused to the EA component of β -pal. When the ligand engages the high affinity receptor, it causes it to functionally dimerize with its heterodimer partner, bringing the fragments of the β -pal reporter into close proximity which forces them to reconstitute enzymatic function, which can be detected by the addition of enzyme substrate.



Representative examples of assays for interleukin receptors from 6 different families of interleukins / receptors. Each plot shows a dose response for the relevant ligand(s) in a given assa from the indicated family. Data plotted are mean RLU and standard deviation from at least triplicate wells for each dose. These assays are characterized by robust assay windows and

imerization of inflammation targets IL-23R and IL-12R is highly specif



Dimerization of IL-23R and IL-12R is highly specific. A: IL23R and IL12R share a common subunit (IL12RB1). Each receptor binds a heterodimeric ligand, which share the p40 subunit. Both receptors have been linked to Crohn's disease by GWAS studies. B. Dimerization of IL-12R (IL12RB1/IL12RB2) is stimulated by IL-12, but not by IL-23 or p40 monomer. C. Dimerization of IL-23R (IL23R /IL12RB1) is stimulated by IL-23 or p40 monomer.



DISCOVER, www.discoverx.com