

Jennifer Brogdon<sup>1</sup>, Susanne Szabo<sup>1</sup>, Sylvie Sulocha<sup>2</sup> and Fred King<sup>3</sup>

<sup>1</sup> Novartis

<sup>2</sup> CisBio

<sup>3</sup> GNF

Abstract:

The aberrant production of T cell cytokines can lead to autoimmunity or asthma. Identification of small molecule inhibitors for cytokine production can be challenging due to the lack of appropriate cell lines which mimic T cell cytokine regulation. We developed an HTRF-based method to measure cytokine production from primary human T cells and adapted the assay to automation for ultra high throughput screening in 1536 well format. Assay development, optimization and screening of primary cells will be discussed.

Resume:

Jennifer Brogdon is currently a Research Investigator at the Novartis Institutes for Biomedical Research working on early drug discovery efforts in the Department of Developmental and Molecular Pathways. She obtained her PhD from the Immunology Department at Duke University and expanded her cellular and molecular immunology background with a postdoctoral position in the Immunobiology Department at Yale University. In her three year tenure at Novartis, her research has focused on the development and implementation of two ultra high throughput screens using primary human cells, hit selection and SAR studies, followed by target id and validation with lead compounds.