

Eric Johnson - Abstract

Title: Application of TR-FRET Systems in Kinome Profiling, Mechanistic Studies, and as an Alternative to Conventional Fluorescence Polarization

Abstract: Rationale and strategy for kinome profiling including inhibitor and kinase selection, HTRF vs ³³P-based platform comparison, and assay development will be discussed. Examples of mechanistic studies on kinase inhibitors utilizing the HTRF platform will be given. The development of TR-FRET binding assays to replace/compliment conventional fluorescence polarization methods and the increased utility of the TR-FRET method will be described.