

# Ultra high throughput GPCR screening utilizing CCD imaging systems with HTRF<sup>®</sup>

Scott Feighner

*Merck, Linden, NJ USA*

## Abstract

The Merck Rahway Facility for Automation and Screening Technologies (RY-FAST) is a group with three main functions: a.) compound management for lead optimization programs, b.) provide assay support for medicinal chemistry programs and c.) function as a liason between the uHTS group and therapeutic groups by miniaturizing assays to 1536 and 3456 well formats. This presentation will outline the steps to define a robust protocol for a Gs-coupled GPCR from 384 well format through 1536 well and up to 3456 well microtiter plates using a Viewlux multimode CCD-based plate reader.

## Resume

Dr Feighner received his Ph.D from the Medical College of Virginia/Virginia Commonwealth University in Microbiology in 1979 and completed a Post-doc. at Michigan State University in 1981. He joined Merck Research Laboratories in Rahway immediately following his Post-doc. He have been involved in prokaryotic and human drug discovery projects throughout his career at MRL. He became interested in GPCRs as drug targets in early 1990s and have been involved in de-orphanizing several GPCRs including the ghrelin and motilin receptors. In 2005 he was recruited to help form a new group to centralize lead optimization, compound management and focused screening for all Medicinal Chemistry projects in Rahway.