

## **Denise Williams--Medarex**

### **Title:**

The use of HTRF<sup>®</sup> in an automated workflow for Human MAb screening at Medarex

Please include the following names as second authors on the talk  
Bill Koukis, Claudia Melara, and Richard Theolis

### **Abstract**

Screening for monoclonal antibodies at Medarex traditionally involved the use of ELISA which included several time consuming steps not ideal for automation. Due to our throughput requirements and the need to standardize for increased efficiency, we have developed an automated high throughput screening system at Medarex which utilizes an HTRF<sup>®</sup> assay that works flawlessly integrated with our custom robotics and data management system. Our simple and robust HTRF<sup>®</sup> assay uses commercially available antibodies which are custom labeled for us by Cisbio. The decision to include the HTRF<sup>®</sup> in our automated process has enabled us to screen a much higher number of plates and has shown a steady increase in our antibody output over the last several years.

Medarex is using the HTRF assay in other areas of research and development as well. As a service to our research department, we use the HTRF assay to quantitate the antibody concentration in our Hybridoma supernatants once hits are identified and have also transferred our knowledge to our cell line development group where our lead candidates are transfected into CHO cell lines.

### **Resume**

Denise Williams has been working with Medarex now for about 9 years. She spent the first 4 years as a scientist in the hybridoma development department and then joined the automation effort that was started in 2002. She is currently the assistant director of Research Automation Services at Medarex which is a service based department that works closely with IT and scientific computing to support research in their efforts to make lead antibodies for commercialization.