

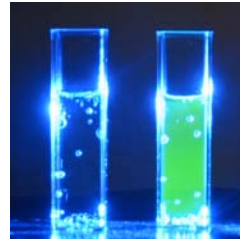
What We Bring To Your Project:



ARBOR
ASSAYS

OUR ABILITIES

- Activity, Immunoassay, and Chemical detection formats
- Low, Medium and High Throughput systems
- Methods for molecular weights 30 to 750,000



HISTORY

- Leaders in developing assays for In Vitro Diagnostics and Life Science assay systems
- FDA In Vitro Diagnostics experience: Abbott Diagnostics, Kallestad, Sanofi, Roche, etc.
- Life Science experience: Founders of Assay Designs - > 150 different assays developed
- Specialized Assay Testing Services
- Small molecule chemical conjugation methods
- Inventor of Acridinium ester chemiluminescent technology used in multiple IVD assay systems.
- Developer of Beckman Access™ immunoassay system
- Extensive knowledge of scale up processes, GMP manufacturing and documentation systems

CURRENT

- Multiple contract assay development and rare reagent research projects
- Developed immunoassays, colorimetric and fluorimetric detection kits being sold worldwide
- Kit to measure palladium contamination levels in API's using technology exclusively licensed
- Universal Histone Demethylase assay patented
- Novel detection and immunoassay kits

ASSAY DEVELOPMENT EXPERTISE

RARE REAGENT GENERATION

The scientists at Arbor Assays are specialized in the design of peptides and organic derivatives to allow for the generation of high quality polyclonal and monoclonal antibodies as well as suitable enzyme and luminescent detection conjugates for assay use. The employees of Arbor Assays have extensive experience in developing some of the best kits currently in use for the measurement of such molecules as hydrogen peroxide, cyclic AMP, oxytocin, phosphokinases, nitric oxide, and IVD molecules.

COMPETITIVE IMMUNOASSAYS

For the development of a competitive assay we will carry out the development of a suitable chemistry route to conjugate the antigen to a carrier protein for antibody development, and/or attach the antigen to a suitable readout system. We are experts in the development of super-fast and/or super-sensitive and multiplexed assays for small molecules. We can guide you on the selection of antibody sources, such as monoclonal or polyclonal and which species to consider.

SANDWICH IMMUNOASSAYS

We will select suitable solid phase antibodies based upon performance criteria developed from your assay specifications. The readout can be from traditional HRP or AP enzymatic colorimetric detection to fluorescent, chemiluminescent or bioluminescent. We also have experience in developing microparticle-based assays from the ground up. The scientists at Arbor Assays have extensive experience with each of these detection methods.

ACTIVITY ASSAYS

Activity assays are substantially different from the basic immunological assays. These assays utilize either a chemical change to determine the presence or amount of a particular molecule, or they can monitor the enzyme activity of a specific protein utilizing some novel read out. We will research various options available and design, synthesize and validate new activity methods. With the expertise in solid phases, homogenous and heterogeneous formats and detection technologies we have the ability to design an assay to meet your needs.

ASSAY TESTING SERVICES

We offer full service testing for samples utilizing our developed assays and other commercially available assays on a wide range of sample types. Full confidentiality is guaranteed with speedy turnaround.

“A customer is the most important visitor on our premises. He is not dependent on us. We are dependent on him. He is not an interruption in our work. He is the purpose of it. He is no outsider in our business. He is part of it. We are not doing him a favor by serving him. He is doing us a favor by giving us an opportunity to do so.”— Mahatma Gandhi

**1514 Eisenhower Place
Ann Arbor, Michigan 48108 USA
Tel: 734-677-1774
Fax: 734-661-1884**



For details concerning these services please contact us:
734-677-1774 • info@ArborAssays.com • www.ArborAssays.com