



# Solutions for Research and Analysis

Developing and Manufacturing Assay Products since 2003!

## BioAssay Systems

offers assay kits that are simple, convenient to use and superior in performance. With our assay kits, researchers need little-to-no time for assay optimization. We specialize in biochemical and cell-based assays for both routine laboratory tests and for high-throughput drug discovery applications with a focus on safe, nonradioactive assay formats such as absorbance, fluorescence and luminescence detection techniques. Our product portfolio consists of assay kits and compatible equipment for a broad range of research areas including:

- Blood/Urine Chemistry
- Energy Metabolism
- Enzyme Activity
- HTS Reagents
- Ion Assays
- Oxidative Stress
- Signal Transduction

For more information, please contact:

### BioAssay Systems

3191 Corporate Place  
Hayward, CA 94545  
USA

Toll free 1-877-782-3888

Tel: 1-510-782-9988

Fax: 1-510-782-1588

Email: [info@bioassaysys.com](mailto:info@bioassaysys.com)

Web: [www.bioassaysys.com](http://www.bioassaysys.com)

## Assay Kits

BioAssay Systems has developed over 160 different assay kits. Below is a list of all assay targets, for a full catalog please visit our website.

**Acetate**  
 Acetylcholine  
 Acetylcholinesterase  
 Aconitase  
 Adipolysis  
 ADP  
 ADP/ATP Ratio  
 Alanine (L-)  
 Alanine Transaminase  
 Albumin (BCG, BCP)  
 Alcohol dehydrogenase  
 Alkaline Phosphatase  
 Ammonia/Ammonium  
 AMPK Phosphorylation  
 Amylase ( $\alpha$ -)  
 Antioxidant  
 Arginase  
 Ascorbic Acid  
 Aspartate  
 Aspartate Transaminase  
 ATP  
 ATPase  
**Bile Acid**  
 Bilirubin  
**Calcium (Ca<sup>2+</sup>)**  
 Caspase-3  
 Catalase  
 Cell Viability MTT  
 Chloride (Cl<sup>-</sup>)  
 Cholesterol  
 Choline  
 Chromium  
 Citrate  
 Coenzyme A  
 Copper (Cu<sup>+</sup>/Cu<sup>2+</sup>)  
 Creatine  
 Creatine Kinase  
 Creatinine  
 Cytotoxicity LDH  
**DNA**  
**ERK Phosphorylation**

Ethanol  
**Fatty Acid**  
 Formaldehyde  
 Fructose  
 Fumarase  
 Fumarate  
**Galactose**  
 Glucose  
 Glucose Dehydrogenase  
 Glucose Oxidase  
 Glucose 6-Phosphate  
 G6P Dehydrogenase  
 Glucosidase ( $\alpha/\beta$ -)  
 Glutamate  
 Glutamate Dehydrogenase  
 Glutamine  
 Glutathione (GSH)  
 Glutathione Peroxidase  
 Glutathione Reductase  
 GSH/GSSG  
 Glycerol  
 Glycogen  
 Glyoxalase I  
 GTPase  
**HDL & LDL/VLDL**  
 Heme  
 Hemoglobin  
**Indican**  
 Indican (Obermeyer)  
 Invertase  
 Iron (Fe<sup>2+</sup>/Fe<sup>3+</sup>)  
 Isocitrate  
 Isocitrate Dehydrogenase  
**Ketone Body**  
 Kinase  
**Lactate (D/L-)**  
 Lactate Dehydrogenase  
 Lactose  
 Lipase  
 Luciferase Reporter  
 Luciferase Dual Reporter

**Magnesium (Mg<sup>2+</sup>)**  
 Malate  
 Malate Dehydrogenase  
 Mannosidase  
 Monoamine Oxidase  
**NAD/NADH**  
 NADP/NADPH  
 Neuraminidase  
 NF $\kappa$ B Phosphorylation  
 Nitric Oxide  
 Nitric Oxide Synthase  
**Oxalate**  
 Oxaloacetate  
**Peroxidase**  
 Peroxide  
 Phenylalanine  
 Phosphatase  
 Phosphate  
 Phospholipase D  
 Phospholipid  
 Protein  
 Pyruvate  
 Pyruvate Kinase  
**Sialic Acid**  
 Sorbitol  
 Sorbitol Dehydrogenase  
 Starch  
 Sucrase  
 Sucrose  
 Sulfate (SO<sub>4</sub><sup>2-</sup>)  
 Superoxide Dismutase  
**TBARS**  
 Total Urinary Protein  
 Triglyceride  
**Urea**  
 Urease  
 Uric Acid  
**Xanthine**  
 Xanthine Oxidase  
**Zinc (Zn<sup>2+</sup>)**

Convenient, Safe and Cost Effective!



## Technical Support

Our technical support team consists of real lab scientists! Your questions are answered directly by the same scientists who developed our assay kits.

## Services

BioAssay Systems' scientists have over 30 years of extensive experience in assay development.

## Analytical Services

If you need a sample to be characterized or tested, look no further than BioAssay Systems. No matter the project our experienced scientists can provide accurate and cost-effective solutions.

## Assay Customization

We can customize any of our existing assays or develop a new assay specifically for your purposes. Contact us for more information.

[www.bioassaysys.com](http://www.bioassaysys.com)