



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
 Web: 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	Ethanol	Magnesium (Mg ²⁺)
Acetylcholine	Fatty Acid	Malate
Acetylcholinesterase	Formaldehyde	Malate Dehydrogenase
Aconitase	Fructose	Mannosidase
Adipolysis	Fumarase	Monoamine Oxidase
ADP	Fumarate	NAD/NADH
ADP/ATP Ratio	Galactose	NADP/NADPH
Alanine (L-)	Glucose	Neuraminidase
Alanine Transaminase	Glucose Dehydrogenase	NFκB Phosphorylation
Albumin (BCG, BCP)	Glucose Oxidase	Nitric Oxide
Alcohol dehydrogenase	Glucose 6-Phosphate	Nitric Oxide Synthase
Alkaline Phosphatase	G6P Dehydrogenase	Oxalate
Ammonial/Ammonium	Glucosidase (α/β-)	Oxaloacetate
AMPK Phosphorylation	Glutamate	Peroxidase
Amylase (α-)	Glutamate Dehydrogenase	Peroxide
Antioxidant	Glutamine	Phenylalanine
Arginase	Glutathione (GSH)	Phosphatase
Ascorbic Acid	Glutathione Peroxidase	Phosphate
Aspartate	Glutathione Reductase	Phospholipase D
Aspartate Transaminase	GSH/GSSG	Phospholipid
ATP	Glycerol	Protein
ATPase	Glycogen	Pyruvate
Bile Acid	Glyoxalase I	Pyruvate Kinase
Bilirubin	GTPase	Sialic Acid
Calcium (Ca ²⁺)	HDL & LDL/VLDL	Sorbitol
Caspase-3	Heme	Sorbitol Dehydrogenase
Catalase	Hemoglobin	Starch
Cell Viability MTT	Indican	Sucrose
Chloride (Cl-)	Indican (Obermeyer)	Sucrose
Cholesterol	Invertase	Sulfate (SO ₄ ²⁻)
Choline	Iron (Fe ²⁺ /Fe ³⁺)	Superoxide Dismutase
Chromium	Isocitrate	TBARS
Citrate	Isocitrate Dehydrogenase	Total Urinary Protein
Coenzyme A	Ketone Body	Triglyceride
Copper (Cu ²⁺ /Cu ²⁺)	Kinase	Urea
Creatine	Lactate (D/L-)	Urease
Creatine Kinase	Lactate Dehydrogenase	Uric Acid
Creatinine	Lactose	Xanthine
Cytotoxicity LDH	Lipase	Xanthine Oxidase
DNA	Luciferase Reporter	Zinc (Zn ²⁺)
ERK Phosphorylation	Luciferase Dual Reporter	

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