

# Solutions for Research & Discovery

Convenient, Safe and Cost-effective

www.bioassaysys.com



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, non-radioactive assay formats such as absorbance, fluorescence and luminescence detection techniques. Our product portfolio consists of assay reagents 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

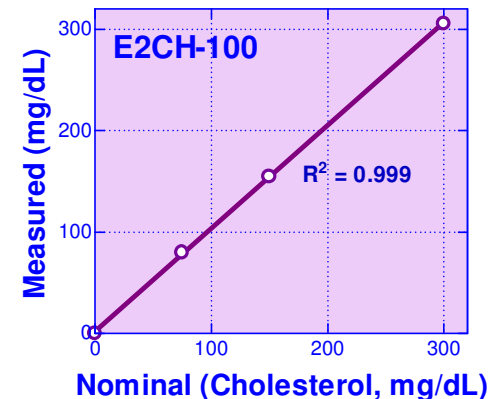
For more information, please contact:

**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

## Handheld Readers

Portable and light-weight, our handheld readers provide exceptional sensitivity, rapid-reading, and reliability at a much lower cost than other portable readers on the market. These convenient readers can be used for applications in many fields such as biology, chemistry, medicine, food safety and environmental monitoring. Built-in memory saves data, which can be retrieved for data analysis by a USB interface. Reads out direct concentration or signal intensity.

| Product Name     | Catalog # | Features   |
|------------------|-----------|--|
| Fluorometers     | FL360450  | Compact handheld fluorometer (exc.360 and em.450nm). Compatible with BioAssay Systems' DNA, ammonia, ammonium, formaldehyde and protein/peptide reagents.  |
|                  | FL530590  | Compact handheld fluorometer (exc.530 and em.590nm). Compatible with BioAssay Systems' 35 assays that measures fluorescence intensity at 530/590nm such as cholesterol, HDL/LDL.                                   |
|                  | FL480530  | Compact handheld fluorometer with exc. 480 and em. 530nm.  |
| Colorimeter      | CL570     | Compact handheld colorimeter with 570nm filter. Compatible with BioAssay Systems' 51 assays that measure 570nm absorbance such as cholesterol, LDL/HDL. Colorimeter with other wavelengths available upon request. |
| Mini glass tubes | MGLTB100  | The transparent high-quality mini glass tubes are designed for use on the handheld readers. Reaction volume: 100 to 200 $\mu$ L.   |



### Cholesterol Assay Performance

Linear Detection Range:

0 to 300 mg/dL

Detection Limit:

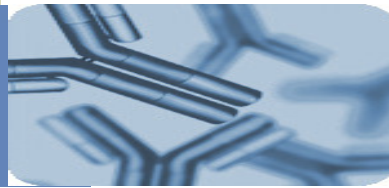
1 mg/dL

Typical Precision (CV%):

<1% at 300 mg/dL;

<1% at 75 mg/dL.

Reads out direct cholesterol concentration or signal intensity.



# Solutions for Research & Discovery

*Convenient, Safe and Cost-effective*

www.bioassaysys.com



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, non-radioactive assay formats such as absorbance, fluorescence and luminescence detection techniques. Our product portfolio consists of assay reagents 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

## Plate Readers

### *Chameleon V: The Most Versatile Multilabel Microplate Reader*

Chameleon V is the only commercially available multilabel microplate reader that offers Liquid Scintillation Counting (LSC) and represents the latest technology in multilabel detection instruments. Having a compact design which includes up to 6 different measurement modes to provide a powerful and flexible platform for performing assays using three different detectors on a unique and patented optical design, Chameleon V achieves optimal performance for filter-based assays, such as absorbance and fluorescence, as well as ultra-high sensitivity single-photon detection for luminescence and liquid scintillation counting.

Chameleon V uses programmed filter slides for both fluorescence and absorbance that allow fast and easy filter exchange by the user. Up to 20 different excitation and 20 different emission filters can be stored for easy access to multiple assays. As an option, Chameleon V can accommodate up to 2 automated injectors within its compact design. Chameleon V offers the greatest flexibility to researchers by providing access to any number of different assays. Choose from several configurations and options to suit your specific requirements:

**Chameleon V/6:** FI, TRF, FP, LSC, LUM, OD

**Chameleon V/5:** FI, TRF, FP, LUM, OD

**Chameleon V/3:** LSC, LUM, OD

**Chameleon V/2:** LSC, LUM

**Chameleon V/1:** LUM

#### Applications

- Receptor binding assays
- Cell proliferation assays (DNA quantification)
- Cytotoxicity assays
- Enzyme assays
- Reporter gene assays
- ATP quantification
- HPLC fraction counting
- RIA, LIA, FIA
- Cell surface assays
- Dot blots
- Coated-well binding
- Organic extraction
- DNA hybridization
- Swipe tests & analysis
- Protein-protein interactions
- More.....

#### Chameleon V Key Features

- Small footprint and lightweight (60 lbs.)
- Uses external PC and Software
- Compatible with virtually any brand of microplate
- 3 Detectors: FL/ABS/ LSC-LUM for max performance in each measurement mode
- Programmable filter slides: 4 filters per slide up to 5 different slides for both excitation and emission
- Can export or directly into Excel
- Upgradeable: from one or more, up to six modes
- Chameleon Options: up to 2 dispensers (automatic injectors), temperature control, cell harvester, robotic arm, and Workstation

#### Affordable, Six-in-One Versatility!

- Liquid Scintillation Counting (LSC)
  - of  $^3\text{H}$ ,  $^{14}\text{C}$ ,  $^{32}\text{P}$ ,  $^{33}\text{P}$ ,  $^{35}\text{S}$ , etc.
- Luminescence (LUM)
- Absorbance (OD or ABS)
- Fluorescence Intensity (FI)
- Time-Resolved Fluorescence (TRF)
- Fluorescence Polarization (FP)

**And its footprint is only slightly larger than a laptop computer!**

