

BioAssay Systems

QuantiQuik™ Test Strips









Rapid, Convenient, Food and Beverage Analysis



- Cost Effective
- Semi-quantitative
- Results in under 15 minutes
- Minimal equipment required



**Ideal for brewing, wine making, and at home food analysis.
Reliable results without expensive lab equipment.**

<u>Quick Test Strip</u>	<u>Detection Range</u>	<u>Samples</u>
 L-Lactic Acid	0 - 360 mg/L or 0 - 4.0 mM	Wine, beer, milk, etc
 D-Lactic Acid	0 - 180 mg/L or 0 - 2.0 mM	Wine, beer, milk, etc
 Malic Acid	0 - 500 mg/L or 0 - 3.7 mM	Wine, beer, milk, etc
 Glucose	0 - 720 mg/L or 0 - 4.0 mM	Wine, beer, juice, etc
 Pyruvate	0 - 131 mg/L or 0 - 1.5 mM	Wine, beer, juice, etc
 Glutamate	0 - 300 mg/L or 0 - 2.0 mM	Soups, sauces, milk, etc
 Histamine	0 - 200 mg/L or 0 - 1.8 mM	Seafood, wine, milk, etc
 Urea	0 - 1500 mg/L or 0 - 24 mM	Milk, serum, urine, etc

More Quick Test Strips available soon!

Custom development available upon request.



BioAssay Systems
3191 Corporate Place
Hayward, CA 94545
U. S. A.

Tel: 510-782-9988
Toll Free: 1-877-782-3888
Email: info@bioassaysys.com
Web: www.bioassaysys.com

Also available at
 