

High Throughput Screening

The FLEXMAP 3D™ system combines differentially dyed fluorescent microsphere sets with an innovative instrument design to take bead-based multiplexing technology to the next dimension. Dual sampling channels, increased operating pressure and 384-well plate compatibility deliver faster read times, while other features such as automated probe adjustment, straightforward routine maintenance operations and an intuitive automation and LIS-compatible software interface simplifies operation. The combined of these features make FLEXMAP 3D the platform of choice for flexible, simple high-throughput multiplex testing.



How is FLEXMAP 3D faster?

Improvements that significantly reduce assay runtime:

- Dual sample fluidics paths and increased syringe injection speed for faster sample injection rate
- Optimized sample flow for higher multiplex requirements
- Fast switching, low dead volume, multi-port valve to direct sample flow stream
- Detailed DSP algorithms to handle the high rate of real-time resultant data



Test (50 µl well volume)	Read Time*	Tests/hour
96 well (100 plex, 2500 beads per well)	18 min	32,000
384 well (100 plex, 2500 beads per well)	1 hour 15 min	30,700
96 well (500 plex, 2500 beads per well)	45 min	64,000
384 well (500 plex, 2500 beads per well)	2 hour 15 min	85,000

*Read times measured on 4 dual channel instruments. Actual results may vary.

384 Well Plate Compatibility

The FLEXMAP 3D is compatible with 96 and 384-well plate formats and offers sample plate heating control for thermally sensitive applications. These features combined with the foil piercing sampling probe provide reliable temperature control over a range of 35C to 60C.

Automation/LIS Compatibility

xPONENT® Software controls the FLEXMAP 3D system and offers interfacing options for laboratory information systems (LIS) as well as automation. An optional swivel base for the system further simplifies integration with third-party liquid handlers.

Simpler Maintenance

The FLEXMAP 3D system has multiple features which simplify start-up, shutdown and maintenance operations and greatly enhance usability for HTS applications.

- Automated probe height adjustment
- Automatic calibration for MicroPlex® and Magplex® microspheres, and enhanced PMT settings
- Calibration and verification kits with color coded reagents and CD with importable reagent target values
- System initialization routine combines system warm-up, washes, calibration and verification into one simple daily startup routine



UNITED STATES
Luminex Corporation
12212 Technology Blvd.
Austin, Texas 78727
United States
Tel: 512.219.8020
Fax: 512.219.5195

CANADA
Luminex Molecular Diagnostics
439 University Ave, Suite 900
Toronto, ON M5G 1Y8
Canada
Tel: 416.593.4323
Fax: 416.593.1066

EUROPE
Luminex B.V.
Krombraak 15
4906 CR Oosterhout
The Netherlands
Tel: +31.162.408333
Fax: +31.162.408337

www.luminexcorp.com
info@luminexcorp.com
investor@luminexcorp.com