

Luminex



FLEXMAP 3D[®]

Multiplexed Genomic, Transcriptomic,
and Proteomic Biomarker Analysis



FLEXMAP3D®

Multiplexing Solutions for the High Throughput Laboratory

• Higher Multiplexing = Higher Throughput

Multiplex up to 500 analytes per well. Each microsphere is impregnated with different amounts of three internal dyes. Monitoring the relative intensity of the three signals allows the system to discriminate up to 500 different microsphere sets.

• Enhanced Pipetting = Fast Turnaround Time

The FLEXMAP 3D system has the capability to run 48,000 analytes in less than one hour. Dual sample fluidics paths and increased syringe injection speed facilitates a faster sample injection rate.

• 96 and 384 Well Compatibility

The FLEXMAP 3D accommodates 96 and 384-well formats for enhanced sample volume flexibility and increased throughput. Sample plate heating control also allows the user to run thermally sensitive applications such as nucleic acids.

• Automation/LIS Compatibility

The new xPONENT® 4.0 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.

• Simplified Maintenance

The FLEXMAP 3D system has multiple features allowing walk away start-up, shutdown and maintenance operations, and greatly enhances usability for High Throughput Screening (HTS) applications.

Luminex Instrumentation: Solutions you can rely on.

Since its incorporation in 1995, Luminex has become a global leader in multiplexing solutions. Luminex offers a full family of instruments designed to meet all your multiplexing needs. Whether it's the efficient MAGPIX system, the dynamic Luminex® 100/200™ system or the powerful FLEXMAP 3D® system, we're sure to have a solution that will help you optimize your research capabilities. Learn more about our suite of enabling multiplexing solutions. Only Luminex offers the piece of mind that comes from:

- Peer reviewed publications
- Patented xMAP technology
- Strong network of market leading partners
- Over 50 510k cleared assays
- Thousands of instruments placed worldwide

Ordering Information

Email: orders@luminexcorp.com

Americas: +1.512.336.3550 or +1.866.401.5450

Fax: +1.512.219.0544

EMEA and India: +31.162.408.333

Fax: +31.162.408.330

UNITED STATES
Luminex Corporation
Austin, Texas
Tel: +1.512.219.8020
Fax: +1.512.219.5195

CANADA
Luminex Molecular Diagnostics
Toronto, Ontario
Tel: +1.416.593.4323
Fax: +1.416.593.1066

EUROPE
Luminex B.V.
Oosterhout, The Netherlands
Tel: +31.162.408333
Fax: +31.162.408337

CHINA
Luminex Shanghai Trading Co.
Shanghai, China
Tel: +21.616.50809
Fax: +21.616.50811

JAPAN
Luminex Japan Corporation, Ltd.
Tokyo, Japan
Tel: +81.3.5545.7440
Fax: +81.3.5545.0451

AUSTRALIA
BSD Robotics, A Luminex Company
Brisbane, Queensland
Tel: +61.7.3273.0273
Fax: +61.7.3273.0274

www.luminexcorp.com

info@luminexcorp.com