





Certificate of Quality

xMAP® Classification Calibrator Microspheres, CAL1

| | | |
|---|---------------------|----------------------|
|  REF | Product Number | L100-CAL1 |
|  LOT | Lot Number | B15184 |
|  | Date of Manufacture | 2009-12-16 |
|  | Expiration Date | 2011-12-16 |
| | | <u>Target Values</u> |
| | DD | 10300 |
| | CL1 | 3730 |
| | CL2 | 3655 |

This certificate provides descriptive lot information and assurance that quality requirements have been met for the components, intermediates, and finished product. Luminex Corporation certifies that the xMAP Classification Calibrator Microspheres, CAL1, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control
15 January, 2010

Luminex

An ISO 13485:2003 Quality Certified Company

12212 Technology Blvd • Austin, Texas 78727-6115 USA

Phone: 1.512.381.4397 • Toll-free: 1.877.785.2323 • Fax: 512.219.5195 • www.luminexcorp.com