

Certificate of Quality

xMAP[®] Reporter Calibrator Microspheres, CAL2

| | |
|---------------------|---------------------|
| Product Number | L100-CAL2 |
| Lot Number | A6692 |
| Date of Manufacture | 14 OCT 2008 |
| Expiration Date | 14 OCT 2010 |
| | <u>Target Value</u> |
| RP1 | 3415 |

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 Reporter Calibrator Microspheres, CAL2, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control
07 November, 2008

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