

# Certificate of Quality

## xMAP<sup>®</sup> Reporter Control Microspheres, CON2

|                     |                     |
|---------------------|---------------------|
| Product Number      | L100-CON2           |
| Lot Number          | A5923               |
| Date of Manufacture | 25 SEP 2007         |
| Expiration Date     | 25 SEP 2009         |
|                     | Parameter A1        |
|                     | <u>Target Value</u> |
| Reference 01        | 29.4                |
| Reference 02        | 180                 |
| Reference 03        | 1894                |
| Reference 04        | 16896               |

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 Control Microspheres, CON2, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control  
08 October, 2007

# **Luminex**<sup>®</sup>

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](http://www.luminexcorp.com)