

# Certificate of Quality

## xMAP® Classification Control Microspheres, CON1

Product Number      L100-CON1  
Lot Number            A5428  
Date of Manufacture   28 NOV 2006  
Expiration Date        28 NOV 2008

|              | <u>Parameter A1<br/>Target Value</u> | <u>Parameter B1<br/>Target Value</u> | <u>Parameter C1<br/>Target Value</u> |
|--------------|--------------------------------------|--------------------------------------|--------------------------------------|
| Reference 01 | 10017                                | 723                                  | 201                                  |
| Reference 02 | 10047                                | 66.1                                 | 67.6                                 |
| Reference 03 | 11098                                | 22834                                | 5412                                 |
| Reference 04 | 9071                                 | 2358                                 | 17622                                |
| Reference 05 | 9713                                 | 227                                  | 5753                                 |

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 Control Microspheres, CON1, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control  
07 December, 2006

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