

# Certificate of Quality

## xMAP™ Classification Control Microspheres, CON1

Product Number L100-CON1  
Lot Number A4863  
Expiration Date 03 NOV 2007

	<u>Parameter A1 Target Value</u>	<u>Parameter B1 Target Value</u>	<u>Parameter C1 Target Value</u>
Reference 01	10129	721	217
Reference 02	10134	66.3	69.4
Reference 03	11095	24074	5776
Reference 04	8934	2384	17370
Reference 05	9832	225	5849

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  
19 December, 2005

# **Luminex**

An ISO 9001:2000 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)