





Certificate of Quality

xMAP® Classification Control Microspheres, CON1

 REF	Product Number	L100-CON1
 LOT	Lot Number	B16125
	Date of Manufacture	2010-03-05
	Expiration Date	2012-03-05

	<u>Parameter A1 Target Value</u>	<u>Parameter B1 Target Value</u>	<u>Parameter C1 Target Value</u>
Reference 01	9917	694	207
Reference 02	9928	68.0	65.7
Reference 03	10717	23112	5580
Reference 04	9542	2426	17632
Reference 05	9700	217	5728

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
30 March, 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