

Certificate of Quality

xMAP® Classification Control Microspheres, CON1

Product Number L100-CON1
Lot Number A5922
Date of Manufacture 25 SEP 2007
Expiration Date 25 SEP 2009

	<u>Parameter A1 Target Value</u>	<u>Parameter B1 Target Value</u>	<u>Parameter C1 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
08 October, 2007

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