

# Certificate of Quality

## xMAP® Classification Calibrator Microspheres, CAL1

Product Number	L100-CAL1
Lot Number	A5718
Date of Manufacture	23 MAY 2007
Expiration Date	23 MAY 2009
	<u>Target Values</u>
DD	10050
CL1	3570
CL2	3645

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 Calibrator Microspheres, CAL1, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control  
08 June, 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](http://www.luminexcorp.com)