





# Certificate of Quality

## xMAP® Classification Calibrator Microspheres, CAL1

 REF	Product Number	L100-CAL1
 LOT	Lot Number	B15918
	Date of Manufacture	2010-02-10
	Expiration Date	2012-02-10
		<u>Target Values</u>
	DD	10300
	CL1	3730
	CL2	3655

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  
26 February, 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](http://www.luminexcorp.com)