





Certificate of Quality

xMAP® Classification Calibrator Microspheres, CAL1

	Product Number	L100-CAL1
	Lot Number	B17522
	Date of Manufacture	2010-06-11
	Expiration Date	2012-06-11
		<u>Target Values</u>
	DD	10000
	CL1	3533
	CL2	3582

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
09 July, 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