

Certificate of Quality

xMAP® Classification Calibrator Microspheres, CAL1

Product Number	L100-CAL1
Lot Number	A5990
Date of Manufacture	30 OCT 2007
Expiration Date	30 OCT 2009
	<u>Target Values</u>
DD	10000
CL1	3430
CL2	3505

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
12 November, 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