

Certificate of Quality

xMAP™ Classification Calibrator Microspheres, CAL1, 1000ml

Product Number	40-75660
Lot Number	A5069
Expiration Date	06 APR 2008
	<u>Target Values</u>
DD	10100
CL1	3515
CL2	3651

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
18 April, 2006

Luminex

An ISO 9001:2000 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