

Certificate of Quality

xMAP[®] Classification Calibrator Microspheres, CAL1, 1000ml

Product Number	40-75660
Lot Number	A5667
Date of Manufacture	26 APR 2007
Expiration Date	26 APR 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 May, 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