

# Certificate of Quality

## MagPlex™ Classification Calibrator Microspheres, MCAL1

Product Number	MCAL1-05
Lot Number	A6893
Date of Manufacture	07 JAN 2009
Expiration Date	07 JAN 2011
	<u>Target Values</u>
DD	5098
CL1	3710
CL2	3556

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 MagPlex Classification Calibrator Microspheres, MCAL1, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control  
23 January, 2009

# **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)