





# Certificate of Quality

## MagPlex™ Classification Calibrator Microspheres, MCAL1

	Product Number	MCAL1-05
	Lot Number	B15534
	Date of Manufacture	2010-01-14
	Expiration Date	2012-01-14
		<u>Target Values</u>
	DD	5049
	CL1	3730
	CL2	3655

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  
11 February, 2010

**Luminex**<sup>®</sup>

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)