
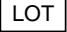
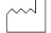



Certificate of Quality

MagPlex™ Classification Control Microspheres, MCON1

 REF	Product Number	MCON1-05
 LOT	Lot Number	B17084
	Date of Manufacture	2010-05-04
	Expiration Date	2012-05-04

	<u>DD (A1)</u> <u>Target Value</u>	<u>CL1 (B1)</u> <u>Target Value</u>	<u>CL2 (C1)</u> <u>Target Value</u>
Reference 01	5044	751	215
Reference 02	4956	70.7	69.1
Reference 03	5356	23786	5537
Reference 04	4749	2363	17745
Reference 05	4876	240	5840

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 Control Microspheres, MCON1, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control
24 May, 2010

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