



Certificate of Quality

xMAP[®] Reporter Control Microspheres, CON2

REF	Product Number	L100-CON2
LOT	Lot Number	B16126
	Date of Manufacture	2010-03-19
	Expiration Date	2012-03-19
		Parameter A1
		<u>Target Value</u>
	Reference 01	30.0
	Reference 02	192
	Reference 03	1927
	Reference 04	16119

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 Reporter Control Microspheres, CON2, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control
04 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