

Certificate of Quality

xMAP® Reporter Control Microspheres, CON2

Product Number	L100-CON2
Lot Number	A6043
Date of Manufacture	28 NOV 2007
Expiration Date	28 NOV 2009
	Parameter A1
	<u>Target Value</u>
Reference 01	29.4
Reference 02	180
Reference 03	1894
Reference 04	16896

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 January, 2008

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