


Certificate of Quality

xMAP™ Reporter Control Microspheres, CON2

Product Number	L100-CON2
Lot Number	A4729
Expiration Date	26 JUL 2007
	Parameter A1 <u>Target Value</u>
Reference 01	28.4
Reference 02	187
Reference 03	1936
Reference 04	15988

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
09 August, 2005

Luminex®

An ISO 9001:2000 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