

Certificate of Quality

xMAP[®] Reporter Calibrator Microspheres, CAL2

Product Number	L100-CAL2
Lot Number	A5178
Date of Manufacture	13 JUN 2006
Expiration Date	13 JUN 2008
	<u>Target Value</u>
RP1	3515

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 Calibrator Microspheres, CAL2, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control
21 June, 2006

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