

Certificate of Quality

xMAP[®] Reporter Calibrator Microspheres, CAL2

Product Number	L100-CAL2
Lot Number	A6063
Date of Manufacture	11 DEC 2007
Expiration Date	11 DEC 2009
	<u>Target Value</u>
RP1	3415

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
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