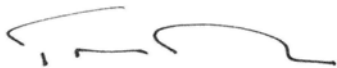


# Certificate of Quality

xMAP® Reporter Calibrator Microspheres, CAL2, 1000ml

Product Number	40-75662
Lot Number	A5978
Date of Manufacture	24 OCT 2007
Expiration Date	24 OCT 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  
06 November, 2007

## **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](http://www.luminexcorp.com)