

Certificate of Quality

xMAP[®] Reporter Calibrator Microspheres, CAL2, 1000ml

Product Number	40-75662
Lot Number	A6004
Date of Manufacture	05 NOV 2007
Expiration Date	05 NOV 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
20 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