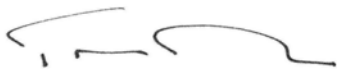


Certificate of Quality

Bio-Plex™ Classification Calibrator Microspheres, CAL1

Product Number	40-75012
Lot Number	A5857
Expiration Date	10 SEP 2009
Concentration (beads/mL)	3.871E+05
Volume (mL)	10.0
DD Target Value	5829
CL1 Target Value	3570
CL2 Target Value	3645

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 Bio-Plex Classification Calibrator Microspheres, CAL1, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control
20 September, 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