

Certificate of Quality

Bio-Plex™ Reporter Calibrator Microspheres, CAL2

| | |
|--------------------------|-------------|
| Product Number | 40-75013 |
| Lot Number | A6279 |
| Expiration Date | 07 APR 2010 |
| Concentration (beads/mL) | 4.415E+05 |
| Volume (mL) | 10.0 |
| High RP1 Target Value | 15538 |
| Low RP1 Target Value | 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 Bio-Plex 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
28 April, 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