

# Certificate of Quality

## Luminex 100/200 Calibration Kit LX200-CAL

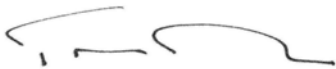
Product Number	LX200-CAL-K25
Lot Number	A7263
Kit Expiration Date	02 APR 2011

---

### Kit Components

Kit Components	Part Number	Lot*
xMAP Classification Calibrator, Microspheres, 5 mL	L100-CAL1	A7167
MagPlex Classification Calibrator, Microspheres, 5 mL	MCAL1-05	A7168
xMAP Reporter Calibrator, Microspheres, 5 mL	L100-CAL2	A7127
8-Well Microtiter plate strips, Pack of 28	13-52032	N/A
LX100/200 Calibrator Kit CD (contains Kit Lot Importable File, component CoQs, MSDS and Kit Instructions)	89-20191-00-001	A7263

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 Luminex 100/200 Calibration Kit, LX200-CAL, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control  
29 May, 2009

\*Reagent target values may vary lot to lot. Only the reagent lots listed should be used when performing system calibration with this kit lot.

# **Luminex**<sup>®</sup>

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)