


# Certificate of Quality

## Luminex 100/200 Calibration Kit LX200-CAL

REF	Product Number	LX200-CAL-K25
LOT	Lot Number	B13899
	Kit Expiration Date	2011-08-11

### Kit Components

Kit Components	Part Number	Lot*
xMAP Classification Calibrator, Microspheres, 5 mL	L100-CAL1	B13503
MagPlex Classification Calibrator, Microspheres, 5 mL	MCAL1-05	B13269
xMAP Reporter Calibrator, Microspheres, 5 mL	L100-CAL2	B13665
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	B13899

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  
21 October, 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)