

# Certificate of Quality

## **Luminex**<sup>®</sup> 100/200 Calibration Kit LX200-CAL-K25

**REF** Product Number

**LOT** Lot Number

 Kit Expiration Date

### **Kit Components**

xMAP Classification Calibrator, Microspheres, 5 mL

MagPlex Classification Calibrator, Microspheres, 5 mL

xMAP Reporter Calibrator, Microspheres, 5 mL

8-Well Microtiter plate strips, Pack of 25

LX100/200 Calibrator Kit CD (contains Kit Lot Importable File,  
component CoQs, MSDS and Kit Instructions)

### **Part Number**

L100-CAL1

MCAL1-05

L100-CAL2

13-52047

89-20191-00-001

### **Lot\***

N/A

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-K25, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.

\*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>

EN ISO 13485 **C**ertified 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)