

# Certificate of Quality

## Luminex® 100/200 Performance Verification Kit LX200-CON-K25

REF Product Number

LOT Lot Number

Kit Expiration Date

### Kit Components

### Part Number

### Lot\*

xMAP Classification Control, Microspheres, 5 mL	L100-CON1	
MagPlex Classification Control, Microspheres, 5 mL	MCON1-05	
xMAP Reporter Control, Microspheres, 5 mL	L100-CON2	
xMAP Fluidics 1, Microspheres, 5 mL	FLUID1-05	
xMAP Fluidics 2, Microspheres, 5 mL	FLUID2-05	
8-Well Microtiter plate strips, Pack of 25	13-52047	N/A
LX100/200 Performance Verification Kit CD (contains Kit Lot Importable File, component CoQs, MSDS and Kit Instructions)	89-20192-00-001	

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 Performance Verification Kit, LX200-CON-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 verification with this kit lot.

# **Luminex**

**EN ISO 13485 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)