

Certificate of Quality

Luminex® 100/200 Performance Verification Kit, RUO LX2R-PVER-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)	89-20778-00-001	

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, RUO, LX2R-PVER-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