

# Certificate of Quality

## FLEXMAP 3D™ Performance Verification Kit F3D-PVER

Product Number	F3D-PVER-K25
Lot Number	A7282
Kit Expiration Date	2011-01-23

---

### Kit Components

Kit Components	Part Number	Lot*
FLEXMAP 3D Classification Verifier, Microspheres, 5 mL	F3DVER1-05	B11752
FLEXMAP 3D e Classification Verifier, Microspheres, 5 mL	F3DeVER1-05	B11988
FLEXMAP 3D Reporter Verifier, Microspheres, 5 mL	F3DVER2-05	B12297
xMAP Fluidics 1, Microspheres, 5 mL	FLUID1-05	B11582
xMAP Fluidics 2, Microspheres, 5 mL	FLUID2-05	B11583
8-Well Microtiter plate Strip wells, Pack of 28	13-52032	N/A
FLEXMAP 3D Performance Verification Kit CD (contains Kit Lot Importable File, component CoQs, and Kit Instructions)	89-20190-00-001	A7282

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 FLEXMAP 3D Performance Verification Kit, F3D-PVER, lot defined above conforms to all prescribed acceptance criteria and is fit for its intended use.



Travis Davies, Director of Quality Control  
08 June, 2009

\*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**®

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)