



CERTIFICATE OF ANALYSIS

COA – 07156

Product Description: MULTI-ARRAY® 96-well Plate Pack

Catalog #/Size: L15XA-1 / 1 Plate; L15XA-3 / 10 Plates; L15XA-6 / 100 Plates; L15XA-7 / 500 Plates

Lot Number: 21F4TA

Expiration Date: 31 DEC 2017

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

BON (Electrical resistivity): < 200 microohms in all circuits

Leak test: No leaks

Free Tag 15 000 test

ECL (mean ± tolerance): 15 000 ± 3,000 cts

Max Average Intraplate CV for any spot: 4%

Maximum Intraplate CV for any plate: 7%

Max Interplate CV for any spot: 7%

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Xiomara Talavera	Quality Control	<i>Talavera</i>	25 JUL 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	25 JUL 2016