



CERTIFICATE OF ANALYSIS

COA – 07986

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

Catalog #/Size: L15XA - Series

Lot Number: 21G45A

Expiration Date: 31 OCT 2018

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 \pm tolerance): 15 000 \pm 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>X Talavera</i>	10 MAR 2017
Review/Approval	Vanessa Castleberry	Quality Assurance	<i>Vanessa Castleberry</i>	13 MAR 2017