



### CERTIFICATE OF ANALYSIS

COA - 03648

**Product Description:** MULTI-ARRAY<sup>®</sup> 96-well Plate Pack  
**Catalog #/Size:** L15XA-1 / 1 Plate; L15XA-3 / 10 Plates; L15XA-6 / 100 Plates; L15XA-7 / 500 Plates  
**Lot Number:** 21D0XA  
**Expiration Date:** 31 MAY 2015  
**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	Rolanda Spratley	Biologics/Chemistry		07 JUN 2013
Review/Approval	Michael Eager	Quality Control		03 JUL 2013