



## CERTIFICATE OF ANALYSIS

COA - 03677

**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:** 21C2ZA

**Expiration Date:** 31 DEC 2014

**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


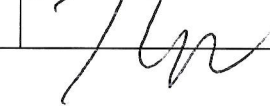
ECL (mean ± tolerance): 15 000 ± 3,000 cts

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		18 Jun 2013
Review/Approval	Michael Eager	Quality Control		03 JUL 2013