



## CERTIFICATE OF ANALYSIS

COA - 04237

**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:** 21D2FA

**Expiration Date:** 30 JUN 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  $\pm$  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.**

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		18 MAR 2014
<b>Review/Approval</b>	Michael Eager	Quality Control		18 MAR 2014