



## CERTIFICATE OF ANALYSIS

COA – 06335

**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:** 21F0GA

**Expiration Date:** 28 FEB 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  $\pm$  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 |  | 16 DEC 2015 |
| <b>Review/Approval</b> | Dawn Stollar     | Quality Assurance   |  | 17 DEC 2015 |