



### CERTIFICATE OF CONFORMANCE

Page 1 of 1

**Product Description:** MULTI-ARRAY® 96-well Plate / Plate Pack

**Catalog Number:** L15XA-Series

**Lot Number:** 21M2RA

**Expiration Date:** 30 JUN 2024

**Storage Temperature:** Room Temperature

**Product Specifications:**

|                              |                                 |
|------------------------------|---------------------------------|
| BON (Electrical resistivity) | < 200 microohms in all circuits |
| Leak Test                    | No Leaks                        |

**Free Tag 15,000 Test Specifications:**

|                                            |                  |
|--------------------------------------------|------------------|
| ECL (mean ± tolerance)                     | 15000 ± 3000 cts |
| Maximum Average Intraplate CV for Any Spot | 4.0%             |
| Maximum Intraplate CV for Any Plate        | 4.0%             |
| Maximum Interplate CV for Any Spot         | 4.0%             |

This product was manufactured and tested according to MSD documents. This lot meets MSD's specifications. This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

|                        | Name             | Function | Signature | Date        |
|------------------------|------------------|----------|-----------|-------------|
| <b>Review/Approval</b> | Xiomara Talavera | Quality  |           | 27 JUN 2022 |