



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

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

**Catalog Number:** L15XA-Series

**Lot Number:** 21L10A

**Expiration Date:** 31 MAR 2023

**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 $\pm$ tolerance)                 | 15000 $\pm$ 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        |
|-----------------|---------------|----------|-----------|-------------|
| Review/Approval | Karen Hamilla | Quality  |           | 16 MAR 2021 |