



### CERTIFICATE OF CONFORMANCE

Page 1 of 1

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

**Catalog Number:** L15XB-Series

**Lot Number:** 22L10A

**Expiration Date:** 30 NOV 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 ± tolerance)                     | 23461 ± 4692 cts |
| Maximum Average Intraplate CV for Any Spot | 3.0%             |
| 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  | <i>Karen Hamilla</i> | 13 JAN 2022 |