



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

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

**Catalog Number:** L15XB-Series

**Lot Number:** 22L02A

**Expiration Date:** 31 JAN 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)                 | 23461 $\pm$ 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 | Xiomara Talavera | Quality  |           | 18 FEB 2021 |