



CERTIFICATE OF CONFORMANCE

Page 1 of 1

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

Catalog Number: L15XA-Series

Lot Number: 21M4PA

Expiration Date: 31 OCT 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 |
|-----------------|------------------|----------|-----------|-------------|
| Review/Approval | Xiomara Talavera | Quality | | 21 OCT 2022 |