



CERTIFICATE OF CONFORMANCE

Page 1 of 1

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

Catalog Number: L15XB-Series

Lot Number: 22L0BA

Expiration Date: 30 APR 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% |
| Maximum Intraplate CV for Any Plate | 4% |
| Maximum Interplate CV for Any Spot | 4% |

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 | | 11 MAY 2021 |