



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MULTI-ARRAY® 96-well Plate / Plate Pack
Catalog Number: L15XA-Series
Lot Number: 21M4SA
Expiration Date: 31 OCT 2024
Storage Temperature: Room Temperature

Product Specifications:

Table with 2 columns: Specification Name, Value. Rows: BON (Electrical resistivity) < 200 microohms in all circuits; Leak Test No Leaks

Free Tag 15,000 Test Specifications:

Table with 2 columns: Specification Name, Value. Rows: 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.

Table with 5 columns: Review/Approval, Name, Function, Signature, Date. Row: Review/Approval, Xiomara Talavera, Quality, [Signature], 26 OCT 2022