



### CERTIFICATE OF CONFORMANCE

Page 1 of 1

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

**Catalog Number:** L15XA-Series

**Lot Number:** 21K3RA

**Expiration Date:** 30 SEP 2022

**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
<b>Review/Approval</b>	Xiomara Talavera	Quality	<i>Talavera</i>	22 SEP 2020