



CERTIFICATE OF CONFORMANCE

Page 1 of 1

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

Catalog Number: L15XA-Series

Lot Number: 21J3ZA

Expiration Date: 30 SEP 2021

Storage Temperature: Room Temperature

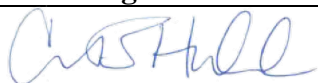
Product Specifications:

BON (Electrical resistivity)	< 200 microohms in all circuits
Leak Test	No Leaks

Free Tag 15,000 Test Specifications:

ECL (mean \pm tolerance)	15 000 \pm 3 000 cts
Maximum Average Intraplate CV for Any Spot	4%
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	Carissa Hubbard	Quality		03 OCT 2019