



## CERTIFICATE OF CONFORMANCE

Page 1 of 1

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


**Catalog Number:** L15XA - Series

**Lot Number:** 21N3ZA

**Expiration Date:** 31 AUG 2025

**Storage Temperature:** Room Temperature

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.

	<b>Name</b>	<b>Function</b>	<b>Signature</b>	<b>Date</b>
<b>Review/Approval</b>	Xiomara Talavera	Quality		12 SEP 2023