



**CERTIFICATE OF CONFORMANCE**

**Product Description:** MULTI-ARRAY® 96-well High Bind Streptavidin Plate  
**Catalog Number:** L15SB-Series  
**Lot Number:** Z0022256  
**Expiration Date:** 30 APR 2026  
**Storage Temperature:** 2-8°C

**Uniformity Results:**

Metric	Specification
Average Intraplate CV	≤ 10.0%
Number of plates with Intraplate CV > 14%	0 Plates
Interplate CV	≤ 12%

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		03 JAN 2024