



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: MULTI-ARRAY® 96-well High Bind Streptavidin Plate

Catalog Number: L15SB-Series

Lot Number: Z0021979

Expiration Date: 31 DEC 2024

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
Review/Approval	Xiomara Talavera	Quality		07 DEC 2022