



CERTIFICATE OF CONFORMANCE

Product Description: MULTI-ARRAY® 96-well High Bind Streptavidin Plate
Catalog #: L15SB - Series
Lot Number: Z0021541
Date of Manufacture: 10 JUN 2019
Expiration Date: 30 NOV 2021

Statement: This product was manufactured and tested according to approved MSD Standard Operating Procedures. This lot meets MSD Specifications

Metric	Specification
Mean Intraplate CV	≤ 10 %
Interplate CV	≤ 12 %
Number of plates with intraplate CV ≥14	0 Plates

The above product is intended for research use only. Not for use in diagnostic procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality	<i>Karen Hamilla</i>	25 JUN 2020