



CERTIFICATE OF CONFORMANCE

Product Description: MULTI-ARRAY® 384-Well SECTOR Imager 6000 Streptavidin Plate
Catalog Number: L21SA-Series
Lot Number: Z0022153
Expiration Date: 30 NOV 2025
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		10 OCT 2023