



CERTIFICATE OF CONFORMANCE

Page 1 of 1

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

Catalog #/Size: L15SB - Series

Lot Number: Z0021417

Expiration Date: 30 SEP 2020

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

Specifications:

Release method: Coating uniformity

Mean intraplate CV: $\leq 10.0\%$

Mean interplate CV: $\leq 12.0\%$

No plate with CV $> 14.0\%$

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	13 NOV 2018