



CERTIFICATE OF ANALYSIS

COA – 07613

Product Description: MULTI-ARRAY® MA6000 384-Well Avidin Plate

Catalog #/Size: L21AA-1

Lot Number: Z0010610

Expiration Date: 28 FEB 2017

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

Specifications:

Release method: Coating uniformity

Mean Intraplate CV: $\leq 10.0\%$

Mean Interplate CV: $\leq 12.0\%$

No more than one plate with CV $> 12.0\%$

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
Preparer	Xiomara Talavera	Quality Control	<i>X Talavera</i>	09 DEC 2016
Review/Approval	Dawn Stollar	Quality Assurance	<i>Dawn Stollar</i>	13 DEC 2016