



CERTIFICATE OF ANALYSIS

COA - 07627

Product Description: MULTI-ARRAY® 96-well High Bind Plate Pack
Catalog #/Size: L15XB-series
Lot Number: 22G10A
Expiration Date: 31 OCT 2018

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

BON (Electrical resistivity): < 200 microohms in all circuits

Leak test: No leaks

Free Tag 15,000 test

ECL (mean ± tolerance): 23 461 ± 4,692

Max Average Intraplate CV for any spot: 3%

Maximum Intraplate CV for any plate: 7%

Max Interplate CV for any spot: 7%

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

	Name	Function	Signature	Date
Preparer	Carissa Hubbard	Quality Control		13 DEC 2016
Review/Approval	Dawn Stollar	Quality Assurance		15 DEC 2016