



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MULTI-SPOT® 96-well 7 Spot Plate Pack

Catalog #/Size: 3-2025-020002-A

Lot Number: 29H10A

Expiration Date: 30 JUN 2019

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): 15 000 ± 3,000 cts

Max Average Intraplate CV for any spot: 5%

Maximum Intraplate CV for any plate: 7%

Max Interplate CV for any spot: 8%

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>	15 JAN 2018