



CERTIFICATE OF CONFORMANCE

Page 1 of 1

Product Description: MULTI-SPOT® 96, SI, 7 Spot Std, Plate Assy, Lipless Top

Catalog Number: N/A

Lot Number: 29I05A

Expiration Date: 29 FEB 2020

Storage Temperature: Room Temperature

Product Specifications:

BON (Electrical resistivity)	< 200 microohms in all circuits
Leak Test	No Leaks

Free Tag 15,000 Test Specifications:

ECL (mean ± tolerance)	15 000 ± 3,000 cts
Maximum Average Intraplate CV for Any Spot	5%
Maximum Intraplate CV for Any Plate	7%
Maximum Interplate CV for Any Spot	8%

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	Dawn Stollar	Quality		06 MAY 2019