



## CERTIFICATE OF ANALYSIS

COA - 04411

**Product Description:** MULTI-SPOT<sup>®</sup> 96, SI, 10 Spot Std Plate

**Catalog #/Size:** 3-2026-120002 / 1 Plate

**Lot Number:** 2BD0SA

**Expiration Date:** 28 FEB 2015

**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  $\pm$  tolerance): 15 000  $\pm$  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
<b>Preparer</b>	Rolanda Spratley	Biologics/Chemistry		12 MAY 2014
<b>Review/Approval</b>	Michael Eager	Quality Control		14May2014