



## CERTIFICATE OF ANALYSIS

COA - 04104

**Product Description:** MULTI-SPOT<sup>®</sup> 96 Well, 7 Spot X-Talk CB Plate

**Catalog #/Size:** 3-2059-120002 / 1 Plate

**Lot Number:** 29DY6A

**Expiration Date:** 30 JUN 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 300 000 test



ECL:  $\geq$  175 000 cts in every well

Max Average Intraplate CV for any spot: 10%

Maximum Intraplate CV for any plate: 15%

Max Interplate CV for any spot: 12%

**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		28 JAN 2014
<b>Review/Approval</b>	Michael Eager	Quality Control		10 FEB 2014