



CERTIFICATE OF ANALYSIS

COA - 04103

Product Description: MULTI-SPOT[®] 96 Well, 4 Spot X-Talk CB Plate

Catalog #/Size: 3-2049-120002 / 1 Plate

Lot Number: 25DY6A

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
Preparer	Rolanda Spratley	Biologics/Chemistry		28 JAN 2014
Review/Approval	Michael Eager	Quality Control		11 FEB 2014