



CERTIFICATE OF ANALYSIS

COA - 03742

Product Description: QUICKPLEX™ 96 Plate,1 Spot, STD, Plate Assay

Catalog #/Size: 3-2079-120002-B / 1 Plate

Lot Number: 2MD02A

Expiration Date: 31 JAN 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 ± tolerance): 15 000 ± 3,000 cts

Max Average Intraplate CV for any spot: 4%

Maximum Intraplate CV for any plate: 7%

Max Interplate CV for any spot: 7%

The above product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		08 Jul 2013
Review/Approval	Pu Liu	LSR Qualified Kit Development		08 Jul 2013