



CERTIFICATE OF ANALYSIS

COA - 04116

Product Description: MULTI-SPOT[®] 96, SI, 10 Spot HB Plate

Catalog #/Size: 3-2026-120003 / 1 Plate

Lot Number: 2CD3HA

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 15 000 test

ECL (mean \pm tolerance): 29 000 \pm 5,800 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
Preparer	Rolanda Spratley	Biologics/Chemistry		24 FEB 2014
Review/Approval	Michael Eager	Quality Control		24FEB2014