



CERTIFICATE OF ANALYSIS

COA - 11225

Product Description: Diluent 100

Catalog #/Size: R50AA-2 / 200 mL; R50AA-4 / 50 mL

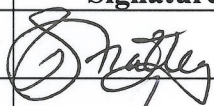

Lot Number: Y0270188

Expiration Date: 30 SEP 2015

Statement: This product was manufactured and tested according to approved MSD documents. This lot meets MSD's specifications.

pH: 7.4 \pm 0.1

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		17 MAY 2013
Review/Approval	Michael Eager	Quality Control		24 MAY 2013