



## CERTIFICATE OF ANALYSIS

COA - 11208

**Product Description:** Diluent 100

**Catalog #/Size:** R50AA-3 / 1000 mL; R50AA-2 / 200 mL; R50AA-4 / 50 mL

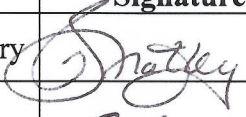
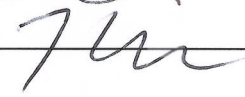
**Lot Number:** Y0270187

**Expiration Date:** 30 JUN 2014

**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		19 APR 2013
Review/Approval	Michael Eager	Quality Control		24 MAY 2013