



CERTIFICATE OF ANALYSIS

COA - 11269

Product Description: Diluent 1

Catalog #/Size: R50CK-2 / 200 mL; R50CK-4 / 50 mL

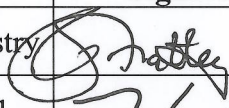

Lot Number: Y0220021

Expiration Date: 31 JUL 2014

Statement: This product was manufactured and tested according to approved MSD documents.

pH: 7.4

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		31 JUL 2013
Review/Approval	Michael Eager	Quality Control		07 JUL 2013