



CERTIFICATE OF ANALYSIS

COA - 06297

Product Description: Diluent 1

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



Lot Number: Y0220026

Expiration Date: 31 AUG 2016

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

pH: 7.7

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		09 DEC 2015
Review/Approval	Dawn Stollar	Quality Assurance		11 DEC 2015