



## CERTIFICATE OF ANALYSIS

COA - 05849

**Product Description:** Diluent 1

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



**Lot Number:** Y0220025

**Expiration Date:** 31 MAR 2016

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

pH: 7.6

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		26 AUG 2015
Review/Approval	Monika Jain	Quality Assurance		26 AUG 2015