



## CERTIFICATE OF ANALYSIS

COA - 04457

**Product Description:** Diluent 1

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



**Lot Number:** Y0220023

**Expiration Date:** 30 APR 2015

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

pH: 7.5

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

	Name	Function	Signature	Date
Preparer	Rolanda Spratley	Biologics/Chemistry		24 SEP 2014
Review/Approval	Anastasia Belilovets	Quality Control		25 SEP 2014