



CERTIFICATE OF CONFORMANCE

COC - 10420

Product Description: MSD® Diluent 3

Catalog #/Size: R51BA-4/5 mL; R51BA-5/25 mL; R51BA-6/125 mL

Lot Number: M0030049

Expiration Date: 30 APR 2014

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

Storage conditions: $\leq -10^{\circ}\text{C}$

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

	Name	Function	Signature	Date
Preparer	Meredith Yeager	Inventory Technician		03/14/2012
Review/Approval	Staci Leffel	Manufacturing Manager		03/13/2012
Review/Approval	Richard Mann	Quality Assurance		03/14/2012