



CERTIFICATE OF CONFORMANCE

COC - 10312

Product Description: MSD® Diluent 3
Catalog #/Size: R51BA-4/5 mL; R51BA-5/25 mL
Lot Number: M0030038
Expiration Date: 31 MAR 2013
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	Manufacturing Associate		12/09/2010
Review/Approval	Staci Leffel	Manufacturing Manager		12/09/2010
Review/Approval	Richard Mann	Quality Assurance		12/09/2010