



CERTIFICATE OF CONFORMANCE

COC - 10352

Product Description: MSD[®] Diluent 3
Catalog #/Size: R51BA-4/5 mL; R51BA-5/25 mL; R51BA-6/125 mL
Lot Number: M0030046
Expiration Date: 30 NOV 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		08/24/2011
Review/Approval	Staci Leffel	Manufacturing Manager		08/19/2011
Review/Approval	Richard Mann	Quality Assurance		08/23/2011