



CERTIFICATE OF CONFORMANCE

COC - 10382

Product Description: MSD® Diluent 101

Catalog #/Size: R51AD-7/700 mL

Lot Number: M1010002

Expiration Date: 30 JUN 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		10/27/2011
Review/Approval	Staci Leffel	Manufacturing Manager		10/27/2011
Review/Approval	Richard Mann	Quality Assurance		10/27/2011