



CERTIFICATE OF CONFORMANCE

COC - 10295

Product Description: MSD® Diluent 3

Catalog #/Size: R51BA-5/25mL

Lot Number: M0030036

Expiration Date: 31 DEC 2012

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		10/18/2010
Review/Approval	Staci Leffel	Manufacturing Manager		10/18/2010
Review/Approval	Richard Mann	Quality Assurance		10/18/2010