



CERTIFICATE OF CONFORMANCE

COC - 10350

Product Description: MSD[®] Diluent 2

Catalog #/Size: R51BB-2/200 mL

Lot Number: M0020027

Expiration Date: 30 NOV 2011

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